

Early identification of hERG liability in drug discovery programs by automated patch clamp

Timm Danker and Clemens Möller

Journal Name:	Frontiers in Pharmacology
ISSN:	1663-9812
Article type:	Methods Article
Received on:	27 Feb 2014
Accepted on:	14 Aug 2014
Provisional PDF published on:	14 Aug 2014
www.frontiersin.org:	www.frontiersin.org
Citation:	Danker T and Möller C(2014) Early identification of hERG liability in drug discovery programs by automated patch clamp. <i>Front. Pharmacol.</i> 5:203. doi:10.3389/fphar.2014.00203
Copyright statement:	© 2014 Danker and Möller. This is an open-access article distributed under the terms of the Creative Commons Attribution License (CC BY) . The use, distribution or reproduction in other forums is permitted, provided the original author(s) or licensor are credited and that the original publication in this journal is cited, in accordance with accepted academic practice. No use, distribution or reproduction is permitted which does not comply with these terms.

This Provisional PDF corresponds to the article as it appeared upon acceptance, after rigorous peer-review. Fully formatted PDF and full text (HTML) versions will be made available soon.

1
2
3
4
5
6
7
8
9
10
11
12
13
14
15
16
17
18
19
20
21
22
23
24
25
26
27
28
29
30
31

Early identification of hERG liability in drug discovery programs by automated patch clamp

Timm Danker(1), Clemens Möller(2)

Abstract

Blockade of the cardiac ion channel coded by hERG can lead to cardiac arrhythmia, which has become a major concern in drug discovery and development. Automated electrophysiological patch clamp allows assessment of hERG channel effects early in drug development to aid medicinal chemistry programs and has become routine in pharmaceutical companies. However, a number of potential sources of errors in setting up hERG channel assays by automated patch clamp can lead to misinterpretation of data or false effects being reported. This article describes protocols for automated electrophysiology screening of compound effects on the hERG channel current. Protocol details and the translation of criteria known from manual patch clamp experiments to automated patch clamp experiments to achieve good quality data are emphasized. Typical pitfalls and artifacts that may lead to misinterpretation of data are discussed. While this article focuses on hERG channel recordings using the QPatch (Sophion A/S, Copenhagen, Denmark) technology, many of the assay and protocol details given in this article can be transferred for setting up different ion channel assays by automated patch clamp and are similar on other planar patch clamp platforms.

Key Words

hERG, Torsades de Pointes, cardiac arrhythmia, electrophysiology, automated patch clamp, safety pharmacology, ion channel, ADMET

Author Affiliations

(1) NMI-TT GmbH, 72770 Reutlingen, Germany

(2) Corresponding author. Albstadt-Sigmaringen University of Applied Sciences, Anton-Günther-Str. 51, 72488 Sigmaringen, Germany. Phone: +49.(0)7571.732 8247. Email: moeller.clemens@gmail.com.

32

33 **1. Introduction**

34 A number of drug discovery and development programs have been hampered by issues with drug
35 induced cardiac arrhythmia. This is particularly well-known for histamine receptor antagonists, e.g. the
36 potent hERG channel blocker Terfenadine (Teldane[®], Seldane[®]). Terfenadine is an H1 receptor
37 antagonist that was launched in 1982 and was later withdrawn from the market because it potentially
38 caused a life-threatening ventricular tachyarrhythmia, torsades de pointes. The potential to affect
39 cardiac ion channel currents and thereby potentially induce cardiac arrhythmia can occur for compounds
40 from many different chemical classes and in very different therapeutic areas.

41 **1.1 hERG channel blockade can cause cardiac arrhythmia**

42 Nearly 20 years ago it was found that mutations in the human ether-à-gogo-related gene (hERG) could
43 cause long QT syndrome. This inherited disorder can be observed in the electrocardiogram as prolonged
44 QT interval and is correlated to torsades de pointes ventricular tachyarrhythmia (Curran *et al.* 1995;
45 Sanguinetti and Tristani-Firouzi 2006; Sanguinetti *et al.* 1995). The protein coded by hERG was later
46 identified as a potassium-selective cardiac ion channel, Kv11.1, which plays an important role during
47 cardiac repolarization. Consequently, malfunction of this ion channel may cause a delay in cardiac
48 repolarization. The importance of the hERG ion channel for drug discovery programs stems from the
49 observation that not only mutations, but also drug induced blockade of the channel, may cause
50 repolarization abnormalities. Many drugs from different chemical classes and therapeutic areas have
51 been shown to block the hERG-coded ion channel and may in turn potentially trigger torsades (Bruin *et al.*
52 2005). Other reasons for cardiac arrhythmias have been identified (Roden *et al.* 1996; Gintant and
53 Hoffman 1984; Gintant *et al.* 2011; Fenichel *et al.* 2004), but hERG channel blockade has become the
54 most frequent single cause for drug withdrawals (Fenichel *et al.* 2004), and many drug discovery
55 programs have been delayed (imposing significant costs on the pharmaceutical company) or stopped due
56 to hERG channel liabilities of potential drug candidates.

57 Examples of compounds that show undesired hERG channel activity include Terfenadine, Astemizole
58 (Zhou *et al.* 1999), brompheniramine (Park *et al.* 2008) and many other drugs as listed on e.g.
59 <http://crediblemeds.org/everyone/composite-list-all-qt-drugs/>, a website maintained by the independent
60 non-profit organization AZCERT, Arizona, and currently sponsored by the Science Foundation Arizona.

61 Reduction of hERG channel liabilities is continuously and extensively discussed in the literature (Tye *et al.*
62 2011; Moorthy *et al.* 2014; Bahl *et al.* 2012; Hudkins *et al.* 2012; Becknell *et al.* 2012; Levoine *et al.* 2011;
63 Davenport *et al.* 2010; Anderson *et al.* 2010; Ito 2009). While compounds from very different chemical
64 classes may interact with the hERG channel due to its relatively large hydrophobic pore, the property of
65 hERG channel liability has been observed especially often for histamine receptor antagonists. The reason
66 for this is that the pharmacophores of the hERG channel and the histamine receptor show remarkable
67 similarities (Davenport *et al.* 2010).

68 **1.2 Automated patch clamp electrophysiology is an emerging technique for identifying cardiac ion** 69 **channel liabilities**

70 To identify potential hERG liabilities early in drug discovery programs and thus avoid problems with hERG
71 channel interactions for late-stage compounds, it has become common practice in drug discovery
72 programs to start testing compounds relatively early during the drug discovery process on potential
73 hERG channel blockade. Finally, before administration to humans, the ICH S7B guideline requests
74 compounds to be tested on potential repolarization issues under GLP, typically employing
75 electrophysiological patch clamp assays which are considered the “gold-standard” for ion channel
76 investigations.

77 A number of assays have been developed to gain a picture of compound effects on the cardiac action
78 potential and in particular on repolarization effects. These assays include hERG ion channel assays
79 employing (but not being limited to) fluorescence, binding, atomic absorption or electrophysiological
80 techniques measuring interaction with, or the function of, the hERG ion channel, with the different
81 throughput inherent to these techniques. In addition, assays have been developed that assess the effects
82 of compounds on other cardiac ion channels. These data may add important information and draw a
83 more complete picture on cardiac effects than possible with hERG channel data alone, since effects on
84 other cardiac ion channels may potentiate or reduce hERG channel effects.

85 Also, assays that could potentially deliver more physiologically relevant data than assays relying on only
86 one ion channel type are used. In the past years, techniques based on stem-cell derived cardiomyocytes
87 particularly gained significant interest (Hamdam *et al.* 2013). Such assays integrate effects on several ion
88 channels beyond the hERG channel. Therefore effects on several channels and interactions of such
89 effects might be detected in assays using these preparations.

90 Despite these efforts in developing, validating and employing more complex assays (which might be
91 considered especially important for valuable late-stage compounds), the hERG ion channel remains a
92 major potential trouble maker in drug discovery programs. It is therefore not surprising that interest in
93 this ion channel was one of the major drivers in developing automated electrophysiology patch clamp
94 instrumentation (Dunlop *et al.* 2008). Automated patch clamp assays have been successfully used in a
95 number of drug discovery programs, both for identifying compound effects on ion channel targets as
96 well as for identifying undesired off-target effects. Technical advances in the automated patch clamp
97 technique and in cell preparations have facilitated employing automated patch clamp assays with
98 increasing success rates (Finkel *et al.* 2006; Zeng *et al.* 2008), high parallelization (and, consequently,
99 increased throughput), at (or near) physiological temperature (Polonchuk 2012), at increased ligand
100 application rates, and also using more challenging cell preparations such as primary cells on these
101 instruments (Jones *et al.* 2009; Stoelzle *et al.* 2011a; Stoelzle *et al.* 2011b; Haythornthwaite *et al.* 2012;
102 Farre *et al.* 2009; Golden *et al.* 2011; Milligan *et al.* 2009; Milligan and Möller 2013; Gillie *et al.* 2013;
103 Becker *et al.* 2013). The development of higher throughput automated patch clamp instruments has
104 allowed moving electrophysiological identification of hERG ion channel assays earlier into the drug
105 discovery process and has made electrophysiology measurements compatible with medicinal chemistry
106 iteration cycles. This is especially important for those drug discovery programs in which hERG channel
107 interactions could be expected from earlier experiences with certain target (such as histamine receptor
108 programs, as noted above) or compound classes.

109 In this manuscript we discuss protocols particularly suited for measuring hERG ion channel effects during
110 the Hit-to-Lead and Lead Optimization phases in drug discovery programs. While during high-throughput
111 screening a small percentage of false negatives or false positives could potentially be tolerated, this is
112 generally less acceptable during later stages of compound development, as false or misinterpreted data
113 might guide medicinal chemistry programs in the wrong direction. Therefore, for this article, we focus on
114 using medium throughput automated patch clamp instrumentation, such as the QPatch™ (Sophion
115 Biosciences A/S, Copenhagen) or the Patchliner (Nanion Technologies GmbH, Munich, Germany)
116 automated electrophysiology platforms, to deliver high data quality. These instruments allow 2 to 8
117 (Patchliner) or 8 to 48 (QPatch) parallel high quality measurements of individual cells in the gigaseal
118 configuration. Importantly, many of the notes and caveats discussed in this manuscript stem from
119 general (and partially basic) electrophysiological considerations. Therefore, this basic electrophysiology,
120 as well as cell preparation details, applies to most of the other automated patch clamp platforms (e.g.,

121 SyncroPatch 96 and SyncroPatch 384 PE (Nanion Technologies GmbH, Munich); PatchXpress, IonWorks
122 Quattro, Barracuda and Barracuda Plus (MDS, Sunnyvale); IonFlux HT (Fluxion Bioscience Inc, San
123 Francisco), Qube (Sophion Biosciences A/S, Copenhagen)), and required validation steps are very similar
124 between the different instruments. Other protocol details for measuring more challenging cell types, in
125 particular for cardiac safety evaluations using stem-cell derived cardiomyocytes using the Patchliner,
126 may be found in other articles (Stoelzle *et al.* 2011a; Milligan and Möller 2013). Protocol optimization for
127 other ion channels can be guided along some of the notes discussed in this article.

128

129

130 **2. Materials**

131 **Cell line**

132 A commercially available cell line ("CHO hERG DUO", BSys, Switzerland) was used (Note 1)

133 **Cell culture and cell harvesting reagents**

134 *Cell culture medium:*

135 450 ml HAM's F-12 + Glutamax (Invitrogen, 31765)

136 + 50 ml FBS Gold (PAA, A15-151)

137 + 1 ml G418 = 100 µg/ml (PAA, P11-012)

138 + 1 ml Hygromycin 100 µg/ml (PAA, P02-015)

139

140 *Serum free medium:*

141 25 ml CHO-S-SFM I (Invitrogen, 12052)

142 + 25 mM HEPES

143 + 0.04 mg/ml Soy bean trypsin inhibitor (Sigma, T65222),

144 + 100 unit/ml Penicillin/Streptomycin (P/S) Invitrogen15140)

145

146 *Thawing medium:*

147 450 ml HAM's F-12 + Glutamax (Invitrogen, 31765)

148 + 50 ml FBS Gold (PAA, A15-151)

149

150 *Harvesting agent:*

151 Accutase (PAA (L11-007))

152

153 **Recording buffers**

154 Extracellular recording buffer (EC), in mM: 145 NaCl, 4 KCl, 1MgCl₂, 2 CaCl₂, 10 HEPES, 10 Glucose;
155 adjust pH to 7.4 with NaOH.

156 Intracellular recording buffer (IC), in mM: 120 KCl, 10 HEPES, 5 CaCl₂, 1.7 MgCl₂, 4 K₂ATP, 10 EGTA;
157 adjust pH to 7.2 with KOH, osmolarity to 292 mOsm with Saccharose.

158

159 **3. Methods**

160 **3.1 Cell culture protocols**

161 **Culturing of cells from frozen vials**

162 1. Prepare 35 ml cell culture thawing medium.

163 1. Thaw the vial quickly in a 37 °C water bath and add to thawing medium.

164 2. Transfer the thawing medium with the cells to a T175 flask containing pre-heated culture
165 medium.

166 3. After 3-4 h, exchange thawing medium with cell culture medium.

167 4. Sub-culture after 2 days.

168 **Sub-culturing**

169 1. Remove culture medium.

170 2. Wash with 10 ml PBS.

171 3. Remove PBS and add 2 ml Accutase.

172 4. Incubate at room temperature (approx. 4 minutes).

- 173 5. Make sure that the cells have a round shape before tapping.
- 174 6. Gently tap on the side of the flask and add 5-7 ml culture medium and resuspend the cells by
175 working the cell suspension up and down 5-10 times.
- 176 7. Determine the cell density and viability by counting the cells in a Hemocytometer using Trypan
177 Blue.
- 178 8. Add the number of cells to the mother flask and the experiment flasks according to the
179 subculturing plan below.
- 180 9. Grow the cells at 37 °C, 5% CO₂ until roughly 80% of the available surface is covered with cells,
181 corresponding to a confluency of the cells of 80%.

182

183 **Sub-culturing plan for making mother flasks and experiment flasks**

- 184 1. Add 1.85×10^6 cells per T175 flask for sub-culturing/experiments after 48 hours.
- 185 2. Add 0.8×10^6 cells per T175 flask for sub-culturing/experiments after 72 hours.
- 186 3. Add 0.3×10^6 cells per T175 flask for sub-culturing/experiments after 96 hours.

187

188 **Cell harvesting for automated electrophysiology patch clamp experiments (for T175 flask)**

189 The cell harvesting for the automated patch clamp experiment should be carried out right before the
190 start of the experiment. Typically, good results can be obtained until up to 4 hours after preparation.

191 The aim of the cell harvesting procedure is to bring the adherently growing cells into suspension while
192 fully maintaining their viability and physiological properties. After preparation, the cell suspension
193 should consist of isolated, single cells. The abundant presence of cell clusters as well as fragments from
194 dead cells would inevitably lead to unsatisfying success rates.

- 195 1. Remove culture medium and wash with 7 ml PBS.
- 196 2. Add 2 ml Accutase per T175 culture flask.
- 197 3. Incubate the culture flask at room temperature for ~4 min (ensure that the cells have reached a
198 round shape and begin to detach before proceeding to the next step) (see note 2).
- 199 4. Gently tap on the side of the culture flask a few times. This should at this time be sufficient to
200 detach almost all of the cells from the flask bottom.

- 201 5. Add 8 ml serum free medium and resuspend the cells by slowly working the cell suspension up
202 and down 5-10 times. This step is crucial to allow a precise counting in the next step. Resuspend
203 just as vigorous as needed to obtain single cells and to avoid cell clusters. Don't do more than
204 needed since this might damage some of the cells. Avoid air bubbles.
- 205 6. Determine the cell density and viability by diluting an aliquot 1:2 in Trypan Blue and count the
206 cells in a Hemocytometer. Also determine the number of cell clusters vs. number of single cells.
- 207 7. Adjust cell density to 3 million cells per ml. You should gain at least 12 ml of cell suspension
208 which are then placed in the cell container of the QPatch ("QStirrer").

209

210 **3.2 Preparation of test compound sample dilutions**

211 For testing with the automated patch clamp systems, dilutions of the test samples in the final test
212 concentration have to be prepared and placed in a 96-well plate. This so called "compound plate"
213 should be prepared immediately before the experiment. The plate can be either a plastic disposable or a
214 reusable tray, accommodating glass inserts. The latter is recommended, since certain types of test
215 compounds tend to adhere to the walls of plastic ware, which may reduce the concentration actually
216 being tested (see also note 7).

217 **3.3 Automated electrophysiological recordings**

218 After loading the compound plates and the harvested cells into the instrument, a previously defined
219 assay protocol is started to carry out the measurements. The assay protocol contains cell type specific
220 settings for establishing the patch clamp recording configuration, as well as "voltage protocols" to elicit
221 the ion channel currents and "application protocols" that determine the order and timing of drug
222 applications.

223 While the settings to achieve the recording configuration for most common hERG expressing cell lines
224 are typically provided by the instrument supplier, and are therefore not discussed here, we will provide
225 detailed information about setting up the voltage and application protocols in the following sections.

226 **3.3.1 Voltage stimulation protocol for hERG**

227 As a voltage gated channel, the hERG channel can be opened and closed by varying the membrane
228 potential. This is accomplished by applying appropriate voltage stimulation protocols, which has to be
229 set up in the appropriate software section of the instrument.

230 Typical for voltage gated potassium channels, the hERG channels opens and partially inactivates at
231 positive voltages. However, in contrast to most other potassium voltage gated channels, the hERG
232 channel shows a very prominent tail current at repolarization which can be even larger than the current
233 recorded during depolarization. The protocol shown in figure 1A shows a widely used stimulation
234 protocol.

235 From a holding potential of -80 mV, the cell is briefly clamped at -50 mV to test leak current at this
236 potential with closed hERG channels (“baseline step”). This is followed by a depolarization to +40 mV,
237 where the hERG channels are alternating between the open and the inactivated state. After returning to
238 -50 mV, the inactivated subpopulation of hERG channels rapidly recovers from inactivation and switches
239 into the open state, from where the channels slowly close. This results in a large tail current to be
240 observed in the second -50 mV phase.

241 A cell expressing hERG channels will respond to this stimulation protocol with a typical current signature,
242 which can be unequivocally identified as a hERG current, as shown in figure 1B.

243 To quantify the hERG current amplitude, values for “baseline” and “peak tail current” are determined by
244 averaging the recorded currents during the baseline step and at the beginning of the second – 50 mV
245 period (these time intervals are marked by the red dotted lines (“cursors”) in figure 1B). The hERG
246 channels are in different states during the two -50 mV phases. In the phase preceding the baseline
247 measurement, the cell is clamped at holding potential and the hERG channels are fully closed. The
248 potential of -50 mV is not sufficient to open closed hERG channels, so the entire channel population of
249 the cell remains closed. Any current recorded during this period is therefore considered not originating
250 from HERG channels (“baseline current” or “leak current”). In contrast, after the depolarization to
251 +40 mV, some of the hERG channels are in an open state, but most of them are in a non-conducting,
252 inactivated state. After repolarizing back to -50 mV, recovery from inactivation is a very fast process.
253 Therefore, the hERG current reaches its highest peak at the beginning of the repolarization to -50 mV
254 (“peak tail current”). The “baseline” level measured before, representing the isolated leak current, will
255 be subtracted from the peak tail current for correction.

256

257 **3.3.2 Application Protocols**

258 For evaluation of drug effects, after establishment of the recording configuration the voltage stimulation
259 protocols are performed in equidistant time intervals of 15 seconds while different liquid solutions are
260 applied to the cell. For analysis, the current amplitudes of each recorded stimulation protocol can be
261 plotted versus time. Figure 2 shows an example of a standard application protocol. After a stabilization
262 period, fresh saline is applied to the cell and a control recording period of at least 3 minutes should be
263 recorded. Current amplitude at beginning and end of this control period should show no significant
264 change. The current amplitude at the end of the control period is then used as a reference point for the
265 measured currents after drug application. (It is good practice to apply each drug concentration multiple
266 times, e.g. twice, as shown in figure 2; see notes 3, 6 and 7.)

267

268

269 **3.3.3 Data filtering and dose response analysis**

270 A typical dose response curve for hERG channel testing will be calculated from a minimum of 3 and up to
271 8 different drug concentrations, where each drug concentration will be tested on at least 3 different
272 cells. While it is common practice to test more than one drug concentration on a single cell (as shown in
273 figure 2), the number of concentrations applied to a cell is limited by the life time of the
274 electrophysiological recording, which is often not more than 30-40 minutes. After this period, relevant
275 quality criteria of the recording (see notes 4, 5) tend to degrade. On the other hand, each drug
276 application should be given enough time to allow the drug effect to reach steady state (see note 6).
277 Therefore, the number of test concentrations that are applied to a cell is limited. The combination of the
278 number of drug concentrations to be tested and a sufficiently long drug exposure time must not exceed
279 the average life time for a stable, high quality recording in whole cell configuration.

280 The test concentrations can be distributed over several cells and the data recombined during analysis.
281 The built in data analysis of the QPatch software is capable to automatically group all data from cells that
282 where treated with the same compound and to calculate the dose response curves accordingly (see
283 figure 3). However, a thorough quality control which filters out all low quality recordings should precede
284 this step.

285 A recording from a cell that qualifies for inclusion into further data analysis, e.g. for dose response curve
286 (DRC) fitting (figure 3) should meet at least the following criteria (see also note 9):

- 287 1) The membrane resistance (R_m) should not fall below a given threshold (e.g. 500 M Ω) at any
288 time (see note 4)
- 289 2) The series resistance (R_s) should not exceed a given threshold (e.g. 10 M Ω) at any time (see
290 note 5)
- 291 3) The initial I_{hERG} peak tail current should not be smaller than a given value (e.g. 250 pA)
- 292 4) The leak current should not be greater than 20% of the initial hERG current at any time
- 293 5) The change in current during the control phase ("run down" or "run up"), should not exceed a
294 given value (e.g. +/- 10%), see note 8.

295

296 To demonstrate the effect of quality control using the key parameters R_m and R_s , we analysed the
297 percentage of remaining hERG current after treatment with 100 nM Terfenadine in a test data set of 22
298 recordings (see table 1). Quality control parameters for R_m and R_s where defined and used to divide the
299 test set into an "accepted" and a "rejected" subgroup, and the mean and standard deviation (SD) of
300 relative remaining currents for each subgroup was calculated. Table 1 shows the effects of applying
301 quality criteria for R_m , R_s , or both. The subsets with good R_m and R_s values show reduced standard
302 deviation and more reliable data. Therefore, especially during routine screens with small sample sizes,
303 excluding data with poor R_s or R_m is highly recommended.

304
305
306
307
308
309
310
311

Table 1: Influence of quality control on the relative remaining hERG current after block with 100 nM Terfenadine. A test set of 22 cells was treated with 100 nM Terfenadine and the relative remaining hERG current was measured. Different quality criteria were applied to divide the test set into two subgroups, named “accepted” and “rejected”. For each criterium, the results of the subgroups are given as mean +/- SD (number of cells).

Quality criterium	Relative remaining current after application of 100 nM Terfenadine	
	Accepted	Rejected
Rm > 150 MΩ	60% ± 13% (n=18)	80% ± 78% (n=4)
Rs < 15 MΩ	61% ± 15% (n=16)	70% ± 60% (n=6)
Rm > 150 MΩ and Rs < 15 MΩ	57% ± 11% (n=14)	74% ± 52% (n=8)

312
313
314

315 On the QPatch and Patchliner assay softwares, these quality criteria can be conveniently checked by
316 setting up a customized plot which shows I_hERG, Rm, Rs on a combined panel. Alternatively, these can
317 be set up as automated quality controls, based on user defined so called “data filters”. Such data filters
318 enable the user to deal with voluminous amounts of data typical for large projects.

319 Of course, rejecting a large number of cells comes at a significant cost. In large screens, a statistical
320 analysis of test datasets may help to find the exact thresholds for quality criteria in order to balance data
321 quality against economics.

322

323

324 4. Notes

325 Note 1: Cell line

326 A number of excellent validated hERG channel expressing cell lines are commercially available (e.g., from
327 ChanTest, Ohio, Cleveland; Millipore Merck KGaA, Darmstadt, Germany; CCS Cell Culture Service,
328 Hamburg, Germany; bSys, Basel, Switzerland; Anaxon, Berne, Switzerland), or can be constructed using
329 well-established molecular biology techniques. Also, cell lines conveniently provided as frozen cells
330 (Donovan *et al.* 2011) are available (e.g., from CytoCentrics, Rostock, Germany and CCS Cell Culture
331 Service, Hamburg, Germany). Especially with the possibility to patch cells directly from frozen stocks,
332 these greatly ease the requirements for cell culture works making cell culture a much more manageable
333 process. When deciding on a cell line, the channel expression rate in suspension should be considered, as
334 this may be significantly different from the expression rate in adherent cells. When channel expression is
335 low after cell harvesting, FCS in the medium as well as incubating and growing the cells at lower
336 temperature can help to recover the current. With some cell lines we have made good experience with

337 incubating the cells at 8 °C for 1 – 3 hours after harvesting before the patch clamp experiment, which has
338 increased the hERG current amplitudes. Vendors of automated electrophysiology instruments make
339 recommendations on cell lines suitable for their instruments, and cell line vendors provide validation
340 data of their lines used on planar patch clamp systems. It has been discussed whether cell lines based on
341 CHO cells or HEK cells provide more relevant data (Wiśniowska and Polak 2009). Independent of which
342 cell type is used, however, it is vital to validate the channel response well, including data of reference
343 compounds, and to make sure that no non-hERG background currents confound the data.

344 **Note 2: Cell detachment**

345 Quality of the cells will critically depend on incubation time in the cell detachment agent before
346 harvesting. When the incubation is too short, cells will not detach from the flask easily and detachment
347 will require a great extent of mechanical interaction, like knocking on the side of the flask. As a result,
348 cells will be damaged when they detach from the flask, resulting in poor success rates in the patch clamp
349 experiments. The amount of cell damage can also easily assessed by a viability staining of the cell
350 preparation with trypan blue and should be less than 5%. When cells are incubated for too long, the seal
351 rate in experiments is often observed to be excellent; however seal stability often becomes poor.

352 **Note 3: Use of double drug applications**

353 In the example given in figure 2, each drug concentration is applied twice (“double drug application”).
354 From the example given, it is easy to see that extrapolation of the time course of compound effect
355 following the first drug application (dashed lines) shows a significant difference to the actual time course
356 after the second drug application of the same concentration, demonstrating that double drug
357 applications can give more accurate results compared to single drug applications. This is particularly
358 important for hydrophobic (“sticky”) compounds, where even more (sometimes more than 4) compound
359 applications may be required before steady state of compound effect is reached, see note 7 and figures 2
360 and 6.

361 **Note 4: Problems introduced by insufficient membrane resistance**

362 The total membrane resistance (R_m) of an intact, healthy cell with closed ion channels is in the order of
363 several G Ω . When the cell is placed on a patch clamp chip and recording configuration is established, a
364 fraction of the cells membrane is sucked into a small hole in the surface of the chip. While the cell
365 membrane in contact with the rim of the hole ideally forms an electrically tight seal with the material of
366 the chip, the membrane in the center of the hole finally becomes ruptured, which allows contact of the
367 amplifier to the interior of the cell through the hole. In this recording configuration (the so called “whole
368 cell” configuration), a high value of the measured membrane resistance indicates that the rest of the cell
369 membrane as well as the cells contact to the chip are intact. This is the prerequisite of high quality
370 recordings. Under conditions where the hERG channels are closed, for example at holding potential or
371 during the baseline step, almost no “leak” currents should be visible in the recording, indicating a high
372 R_m .

373 In figure 4, two example recordings from the same cell, with different R_m , are compared. The blue trace
374 is recorded shortly after obtaining the whole cell configuration. It represents a cell with a good
375 membrane resistance in the $G\Omega$ range. At holding potential (begin and end of the trace) and during the
376 baseline step the recorded “leak” current is close to zero. In contrast, the red trace shows a hERG
377 recording from the same cell after R_m had dropped to $\sim 100\text{ M}\Omega$. The uncorrected peak tail current has
378 dropped from 1600 pA to 1300 pA. At holding potential and during the baseline step, the “leak” currents
379 deviate significantly from zero. The peak tail current can be corrected with the baseline level (as
380 recommended), the corrected peak tail current will be less affected. However, with increasing
381 degradation of R_m the results will finally become highly inaccurate.

382 **Note 5: Problems introduced by large series resistance**

383 The series resistance (R_s), also commonly referred to as access resistance, is the electrical resistance
384 between the amplifier input and the cell membrane of the recorded cell. In automated patch clamp, it is
385 largely determined by the opening diameter of the hole in the patch clamp chip that makes contact to
386 the cell. However, R_s may further be increased when the membrane that spans the hole is only partially
387 ruptured, or for example when cell organelles are being drawn into the hole. With R_s being very large,
388 the amplifier is not able to fully control the electrical potential of the cell membrane. As a result, the
389 actual cell membrane potential may deviate significantly from the potentials defined by the voltage
390 stimulation protocol. The ion channels will therefore not be accurately stimulated, and thus their current
391 responses will be altered.

392 Figure 5 shows an example of the typical artifacts introduced by large R_s when recording hERG currents.

393 As a rule of thumb, when doing dose response recordings, R_s should be as small as possible (preferably
394 $< 10\text{ M}\Omega$) and stable over time.

395

396 **Note 6: Incomplete steady state of drug effects**

397 The time needed for a compound effect to reach steady state varies from compound to compound.

398 Failure to achieve steady state before applying the next compound concentration will lead to
399 underestimation of the drug effect and therefore to artificially left-shifted dose response curves.

400 Whether a sufficient steady state has been reached can be judged by examining the I/t plot of the
401 recording.

402 Terfenadine is an example of a compound for which the results of IC_{50} measurements critically depend
403 on experimental conditions. For manual patch clamp measurements, 23-fold differences have been
404 found in literature values (Kirsch *et al.* 2004). Due to its hydrophobic or “sticky” nature, the compound is
405 particularly difficult to measure accurately with automated patch clamp devices (see note 7). Among
406 several known hERG inhibitors tested (Guo and Guthrie 2005), Terfenadine exhibited one of the largest

407 discrepancies (10 nM vs. 77 nM) in hERG IC₅₀ values when comparing standard patch clamp data to
408 automated data. The compound also reaches steady state of drug block very slowly, so that prolonged
409 drug exposure times have to be considered for accurate dose response assessments.

410 In the “bad practice” example shown in figure 6A, the blocking effect of Terfenadine is evaluated with
411 the same protocol as in figure 2. With the first applied drug concentration of 30 nM, almost no block is
412 reported by the assay. Analysis of the I/t plot clearly reveals that the drug effect develops very slowly, so
413 that steady state is not reached before the next drug concentration is applied. Increasing the drug
414 application time increases current blockade, and repeated drug applications improve the actual drug
415 concentration at the target. A “good practice” example is shown in figure 6B, where the number of
416 repeated drug applications and the drug exposure time is doubled, an acceptable steady state is
417 reached, and a significant drug induced block of 30% is found for the same concentration.

418 Therefore, the assay scheme in the first example leads to an underestimation of the compound effect. A
419 prolongation of the time periods between the drug applications leads to more accurate IC₅₀
420 determinations for such slowly acting compounds.

421 To demonstrate the effect of incomplete steady state, we analyzed the effect of a prolonged drug
422 exposure to Terfenadine.

423 The recording shown in figure 6A is representative for a test data set of 7 recordings featuring a drug
424 exposure time of 150 s per compound. The average IC₅₀ for Terfenadine in this data set is 180 nM (n=7).
425 In contrast, when we doubled the drug exposure time, as shown in figure 6B, we obtained an average
426 IC₅₀ of 72 nM under otherwise similar conditions (see figure 3).

427
428
429

430 **Note 7: Recommendations for sticky compounds**

431 Some compound classes are difficult to dissolve in aqueous solutions and also known for their notorious
432 “sticky” behavior, typically a result of their hydrophobic nature. It has been commonly observed that
433 with these compounds, the actual applied concentration is in fact reduced by adherence of compound
434 molecules to the walls of the containers where the compound dilutions are stored in. The result of such
435 effects would be an artificially right-shifted dose response curve, and therefore may lead to failure of
436 detection of problematic hERG effects.

437 To reduce such unwanted effects, the following precautions are recommended:

- 438 • Use glassware as a standard for compound plates, vials, and even pipette tips.
- 439 • Set up the dilutions in rather large vials (at least 1 ml).
- 440 • Prepare a DMSO stock solution at a 1000 fold higher concentration than the maximum tested
441 concentration.

442 • To prepare the various test concentrations, dilute the stock solution first with DMSO, then dilute
443 1:1000 with EC, thus keeping the final DMSO concentration at a constant level of 0.1%. An
444 increased DMSO concentration of up to 1% may be tolerated in your assay.

445 • If feasible prepare compound dilutions manually and closely look for precipitation.

446 • Prepare the compound plate immediately before running the experiment.

447 It may be required to analyze solubility and final concentrations of compounds in buffer solutions
448 using e.g. LC/MS (when available).

449

450 **Note 8: Run down**

451 From time to time, a downward trend of the hERG current amplitudes during the control phase, .i.e.
452 independent of any drug application, may be observed. This phenomenon is commonly referred to as
453 “run down”. Including cells exhibiting run down into dose response analysis should be strictly avoided.
454 Otherwise, it will result in artificially left shifted dose response curves, and therefore may lead to false
455 positive identification of hERG blockade.

456 The main focus to fix run down problems should be the cell culture. Cell density in the culture flasks
457 should not exceed 80% confluency, and CO₂ concentration be kept sufficiently constant. Using an
458 incubator that is frequently opened (i.e. because it accommodates cell cultures from different projects)
459 may lead to CO₂ fluctuations that might negatively affect cell quality.

460 It is also a good advice to check pH and osmolarity of IC and EC recording buffers before use, especially if
461 frozen stocks are employed.

462

463 **Note 9: Hill coefficient**

464 From dose-response (or IC₅₀)-fits with a sufficient number of concentrations, the Hill coefficient can also
465 be calculated with standard assay software packages. For many channel-compound interactions, the Hill
466 coefficient would be expected to be close to 1. While other Hill coefficients are possible, these should be
467 critically reviewed. Typical problems in experiments that can lead to Hill coefficients incorrectly deviating
468 from one, and, consequently would produce incorrect IC₅₀ values, are:

469 A Hill coefficient greater than one (steep IC₅₀ curve) is often caused by underestimation of compound
470 effects in particular at small concentrations, e.g., due to no steady state in compound effects (see note
471 6), absorption of compound to e.g. tube materials (see note 7), and other reasons.

472 A Hill coefficient smaller than one often stems from underestimation of compound effects at high
473 concentrations (e.g., due to poor voltage clamp conditions at large currents or large series resistance,

474 see note 5), or overestimation of compound effects at low concentrations, e.g. due to run down effects
475 (see note 8), that might be falsely interpreted as slow compound effects.

476

477 **5. Outlook**

478 This article focuses on protocols to identify cardiac safety issues caused by hERG ion channel blockade
479 early in drug discovery programs by high quality medium throughput automated electrophysiology
480 screening using cells expressing the hERG ion channel. The physiological significance of these recordings
481 can be improved, at reduced throughput, by performing recordings at or near physiological temperature
482 (Polonchuk 2012). In addition to screening on the particularly important hERG ion channel, assessing the
483 effects of compounds on other cardiac ion channels may be required. Recently, automated patch clamp
484 screening of compound effects on the action potential in stem cell derived cardiomyocytes has become
485 possible (Stoelzle *et al.* 2011a; Milligan and Möller 2013). These assays provide information to further
486 assess the cardiac safety of compounds. However, currently more validation of stem cell based assays is
487 required before these can be routinely used for cardiac safety investigations.

488 Advances in the parallelization of patch clamp electrophysiology robots will further increase the
489 throughput of patch clamp screening. A robot allowing parallel measurements of up to 768 wells with
490 gigaseal and corresponding low leak currents is now available (Syncropatch PE, Nanion Technologies,
491 Munich), and the throughput of patch clamp electrophysiology screening can be expected to increase
492 even further.

493

494

495

496 **Figure captions**

497

498 **Figure 1:** A, Stimulation protocol for hERG channel measurements. B, Example of hERG recordings
499 showing the effect of Cisapride. Four representative traces from the same cell are superimposed to
500 visualize the effect of increasing concentrations of the known hERG blocker Cisapride.

501

502 **Figure 2:** Current vs time (I/t) plot of a 3 point dose response assessment of Cisapride on hERG currents.
503 Downward arrows at the top of the figure mark time points where saline or test compound is applied,
504 and horizontal bars indicate the duration of exposure. The last 2 data points before the next application
505 (marked by colored points) are averaged to calculate the effect of the preceding drug application. Drug
506 applications are performed twice to ensure saturation of drug effect, where only the second drug
507 application is used for data analysis.

508

509 **Figure 3:** Dose response fit through the data of 9 cells (each identified by a unique color) with 2
510 concentrations of Terfenadine applied to each cell. For each cell, hERG tail current values representative
511 for each compound concentration were determined as shown in figure 3, and normalized to the
512 unblocked hERG current value which was recorded at the end of the saline period before the first
513 compound application. The combined data of all 9 cells was then fitted to the hill equation $y=c^h / IC_{50}^h+c^h$
514 to estimate the half maximal inhibition concentration (IC_{50} value) and hill coefficient h for this compound
515 ($IC_{50} = 72$ nM, $h = 0.8$).

516

517

518

519

520 **Figure 4:** Effect of small membrane resistance (R_m). Two superimposed hERG recordings from a cell with
521 instable membrane resistance are shown. The blue trace represents an early recording from this cell,
522 where R_m has a high value in the $G\Omega$ range. The red trace was recorded from the same cell after R_m
523 dropped to ~ 100 M Ω . The baseline (or "leak") currents recorded during the brief -50 mV "baseline step"
524 are indicated by horizontal dotted lines. The inset in the upper right corner provides a zoomed view on
525 how the baseline current is determined during the baseline step. Due to the drop in R_m , baseline current
526 and peak tail current are shifted to a comparable extent. The difference between baseline and peak tail
527 current, indicated by the double headed vertical arrows, is more robust to changes in R_m than
528 uncorrected peak tail current alone.

529

530 **Figure 5:** Effect of series resistance increase on hERG measurements. In this drastic example, a change in
531 series resistance (R_s), combined with a stable, but critically low membrane resistance ($R_m=120$ M Ω),
532 leads to a significant distortion in the readout.

533

534 **Figure 6:** Good and bad practice examples for steady state of drug induced block by Terfenadine. A, Bad
535 practice example of an I/t plot showing insufficient steady state of block, which would lead to
536 underestimation of the drug effect. B, Good practice example where the drug exposure time as well as
537 the number of drug application repeats has been doubled, leading to an acceptable steady state.

538

539

540

541

542

543

544 References

- 545 Anderson JT, Campbell M, Wang J, Brunden KR, Harrington JJ, and Stricker-Krongrad A et al. 2010.
546 Investigation of 4-piperidinols as novel H3 antagonists. *Bioorg. Med. Chem. Lett.* 20: 6246–6249.
- 547 Bahl A, Barton P, Bowers K, Brough S, Evans R, and Luckhurst CA et al. 2012. The discovery of CCR3/H1
548 dual antagonists with reduced hERG risk. *Bioorg. Med. Chem. Lett.* 22: 6688–6693.
- 549 Becker N, Stoelzle S, Göpel S, Guinot D, Mumm P, and Haarmann C et al. 2013. Minimized cell usage for
550 stem cell-derived and primary cells on an automated patch clamp system. *J Pharmacol Toxicol*
551 *Methods* 68: 82–87.
- 552 Becknell NC, Dandu Rr, Lyons JA, Aimone LD, Raddatz R, and Hudkins RL. 2012. Synthesis and evaluation
553 of 4-alkoxy-[1'-cyclobutyl-spiro(3,4-dihydrobenzopyran-2,4'-piperidine)] analogues as histamine-3
554 receptor antagonists. *Bioorg. Med. Chem. Lett.* 22: 186–189.
- 555 Bruin ML, Pettersson M, Meyboom R, Hoes AW, and Leufkens HG. 2005. Anti-HERG activity and the risk
556 of drug-induced arrhythmias and sudden death. *Eur. Heart J.* 26: 590–597.
- 557 Curran ME, Splawski I, Timothy KW, Vincent GM, Green ED, and Keating MT. 1995. A molecular basis for
558 cardiac arrhythmia: HERG mutations cause long QT syndrome. *Cell* 80: 795–803.
- 559 Davenport AJ, Möller C, Heifetz A, Mazanetz MP, Law RJ, and Ebneith A et al. 2010. Using
560 electrophysiology and in silico three-dimensional modeling to reduce human Ether-a-go-go related
561 gene K(+) channel inhibition in a histamine H3 receptor antagonist program. *Assay. Drug Dev. Technol.*
562 8: 781–789.
- 563 Donovan BT, Bakshi T, Galbraith SE, Nixon CJ, Payne LA, and Martens SF. 2011. Utility of frozen cell lines
564 in medium throughput electrophysiology screening of hERG and NaV1.5 blockade. *J Pharmacol*
565 *Toxicol Methods* 64: 269–276.
- 566 Dunlop J, Bowlby M, Peri R, Vasilyev D, and Arias R. 2008. High-throughput electrophysiology: an
567 emerging paradigm for ion-channel screening and physiology. *Nat. Rev. Drug Discov.* 7: 358–368.
- 568 Farre C, Haythornthwaite A, Haarmann C, Stoelzle S, Kreir M, and George M et al. 2009. Port-a-patch and
569 patchliner: high fidelity electrophysiology for secondary screening and safety pharmacology.
570 *Comb. Chem. High Throughput. Screen.* 12: 24–37.
- 571 Fenichel RR, Malik M, Antzelevitch C, Sanguinetti M, Roden DM, and Priori SG et al. 2004. Drug-induced
572 torsades de pointes and implications for drug development. *J. Cardiovasc. Electrophysiol.* 15: 475–
573 495.
- 574 Finkel A, Wittel A, Yang N, Handran S, Hughes J, and Costantin J. 2006. Population patch clamp improves
575 data consistency and success rates in the measurement of ionic currents. *J. Biomol. Screen.* 11: 488–
576 496.
- 577 Gillie DJ, Novick SJ, Donovan BT, Payne LA, and Townsend C. 2013. Development of a high-throughput
578 electrophysiological assay for the human ether-à-go-go related potassium channel hERG. *J Pharmacol*
579 *Toxicol Methods* 67: 33–44.
- 580 Gintant GA, Gallacher DJ, and Pugsley MK. 2011. The 'overly-sensitive' heart: sodium channel block and
581 QRS interval prolongation. *Br. J. Pharmacol.* 164: 254–259.

582 Gintant GA and Hoffman BF. 1984. Use-dependent block of cardiac sodium channels by quaternary
583 derivatives of lidocaine. *Pflugers Arch.* 400: 121–129.

584 Golden AP, Li N, Chen Q, Lee T, Nevill T, and Cao X et al. 2011. IonFlux: a microfluidic patch clamp system
585 evaluated with human Ether-à-go-go related gene channel physiology and pharmacology. *Assay Drug*
586 *Dev Technol* 9: 608–619.

587 Guo L and Guthrie H. 2005. Automated electrophysiology in the preclinical evaluation of drugs for
588 potential QT prolongation. *J Pharmacol Toxicol Methods* 52: 123–135.

589 Hamdam J, Sethu S, Smith T, Alfirevic A, Alhaidari M, and Atkinson J et al. 2013. Safety pharmacology--
590 current and emerging concepts. *Toxicol. Appl. Pharmacol.* 273: 229–241.

591 Haythornthwaite A, Stoelzle S, Hasler A, Kiss A, Mosbacher J, and George M et al. 2012. Characterizing
592 human ion channels in induced pluripotent stem cell-derived neurons. *J Biomol Screen* 17: 1264–
593 1272.

594 Hudkins RL, Zulli AL, Dandu Rr, Tao M, Josef KA, and Aimone LD et al. 2012. 4-phenoxy piperidine
595 pyridazin-3-one histamine H(3) receptor inverse agonists demonstrating potent and robust wake
596 promoting activity. *Bioorg. Med. Chem. Lett.* 22: 1504–1509.

597 Ito C. 2009. Histamine H3-receptor inverse agonists as novel antipsychotics. *Cent Nerv Syst Agents Med*
598 *Chem* 9: 132–136.

599 Jones KA, Garbati N, Zhang H, and Large CH. 2009. Automated patch clamping using the QPatch.
600 *Methods Mol. Biol.* 565: 209–223.

601 Kirsch GE, Trepakova ES, Brimecombe JC, Sidach SS, Erickson HD, and Kochan MC et al. 2004. Variability
602 in the measurement of hERG potassium channel inhibition: effects of temperature and stimulus
603 pattern. *J Pharmacol Toxicol Methods* 50: 93–101.

604 Levoine N, Labeeuw O, Calmels T, Poupardin-Olivier O, Berrebi-Bertrand I, and Lecomte J et al. 2011.
605 Novel and highly potent histamine H3 receptor ligands. Part 1: withdrawing of hERG activity. *Bioorg.*
606 *Med. Chem. Lett.* 21: 5378–5383.

607 Milligan CJ, Li J, Sukumar P, Majeed Y, Dallas ML, and English A et al. 2009. Robotic multiwell planar
608 patch-clamp for native and primary mammalian cells. *Nat Protoc* 4: 244–255.

609 Milligan CJ and Möller C. 2013. Automated planar patch-clamp. *Methods Mol. Biol.* 998: 171–187.

610 Moorthy NS, Ramos MJ, and Fernandes PA. 2014. Predictive QSAR models development and validation
611 for human ether-a-go-go related gene (hERG) blockers using newer tools. *J Enzyme Inhib Med Chem*
612 29: 317–324.

613 Park S, Kim K, and Kim E. 2008. Blockade of HERG K⁺ channel by an antihistamine drug brompheniramine
614 requires the channel binding within the S6 residue Y652 and F656. *J Appl Toxicol* 28: 104–111.

615 Polonchuk L. 2012. Toward a New Gold Standard for Early Safety: Automated Temperature-Controlled
616 hERG Test on the PatchLiner. *Front Pharmacol* 3: 3.

617 Roden DM, Lazzara R, Rosen M, Schwartz PJ, Towbin J, and Vincent GM. 1996. Multiple mechanisms in
618 the long-QT syndrome. Current knowledge, gaps, and future directions. The SADS Foundation Task
619 Force on LQTS. *Circulation* 94: 1996–2012.

620 Sanguinetti MC, Jiang C, Curran ME, and Keating MT. 1995. A mechanistic link between an inherited and
621 an acquired cardiac arrhythmia: HERG encodes the IKr potassium channel. *Cell* 81: 299–307.

622 Sanguinetti MC and Tristani-Firouzi M. 2006. hERG potassium channels and cardiac arrhythmia. *Nature*
623 440: 463–469.

624 Stoelzle S, Haythornthwaite A, Kettenhofen R, Kolossov E, Bohlen H, and George M et al. 2011a.
625 Automated patch clamp on mESC-derived cardiomyocytes for cardiotoxicity prediction. *J Biomol*
626 *Screen* 16: 910–916.

627 Stoelzle S, Obergrussberger A, Brüggemann A, Haarmann C, George M, and Kettenhofen R et al. 2011b.
628 State-of-the-Art Automated Patch Clamp Devices: Heat Activation, Action Potentials, and High
629 Throughput in Ion Channel Screening. *Front Pharmacol* 2: 76.

630 Tye H, Mueller SG, Prestle J, Scheuerer S, Schindler M, and Nosse B et al. 2011. Novel 6,7,8,9-tetrahydro-
631 5H-1,4,7,10a-tetraaza-cyclohepta[f]indene analogues as potent and selective 5-HT(2C) agonists for
632 the treatment of metabolic disorders. *Bioorg.Med.Chem.Lett.* 21: 34–37.

633 Wiśniowska B and Polak S. 2009. hERG in vitro interchange factors--development and verification.
634 *Toxicol. Mech. Methods* 19: 278–284.

635 Zeng H, Penniman JR, Kinose F, Kim D, Trepakova ES, and Malik MG et al. 2008. Improved throughput of
636 PatchXpress hERG assay using intracellular potassium fluoride. *Assay Drug Dev Technol* 6: 235–241.

637 Zhou Z, Vorperian VR, Gong Q, Zhang S, and January CT. 1999. Block of HERG potassium channels by the
638 antihistamine astemizole and its metabolites desmethylastemizole and norastemizole. *J. Cardiovasc.*
639 *Electrophysiol.* 10: 836–843.

640

641

642

Figure 1.TIF

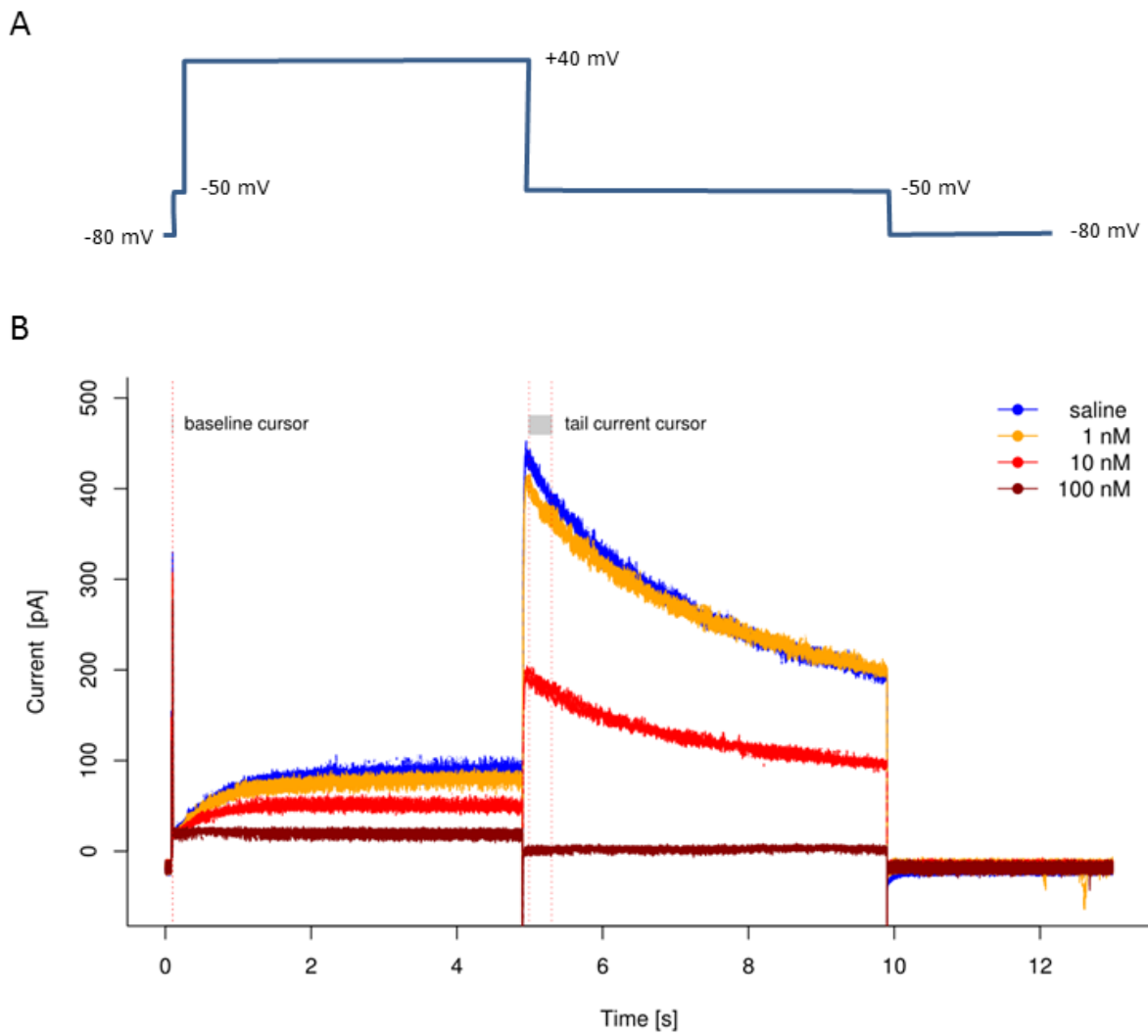


Figure 2.TIF

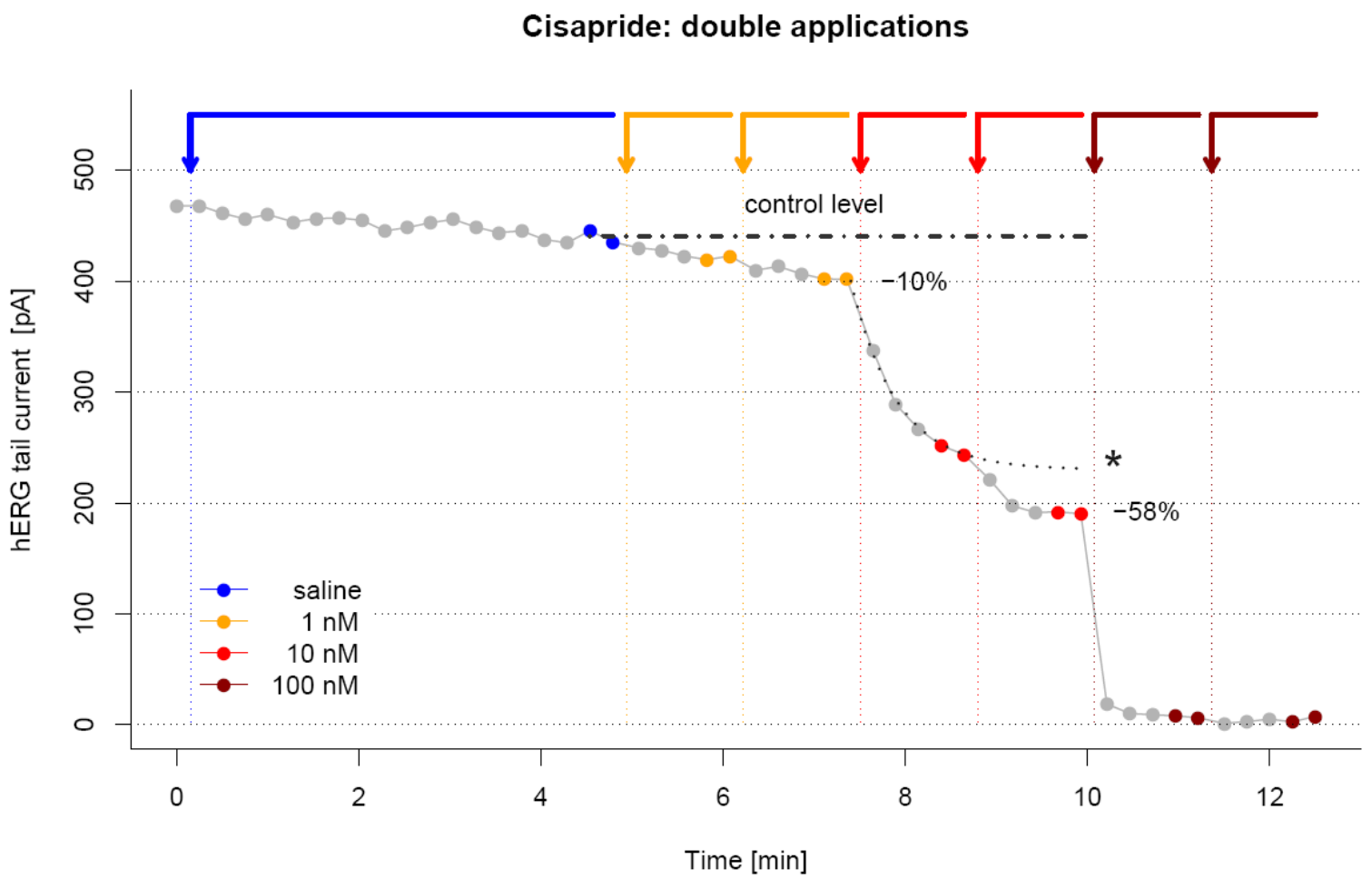


Figure 3.TIF

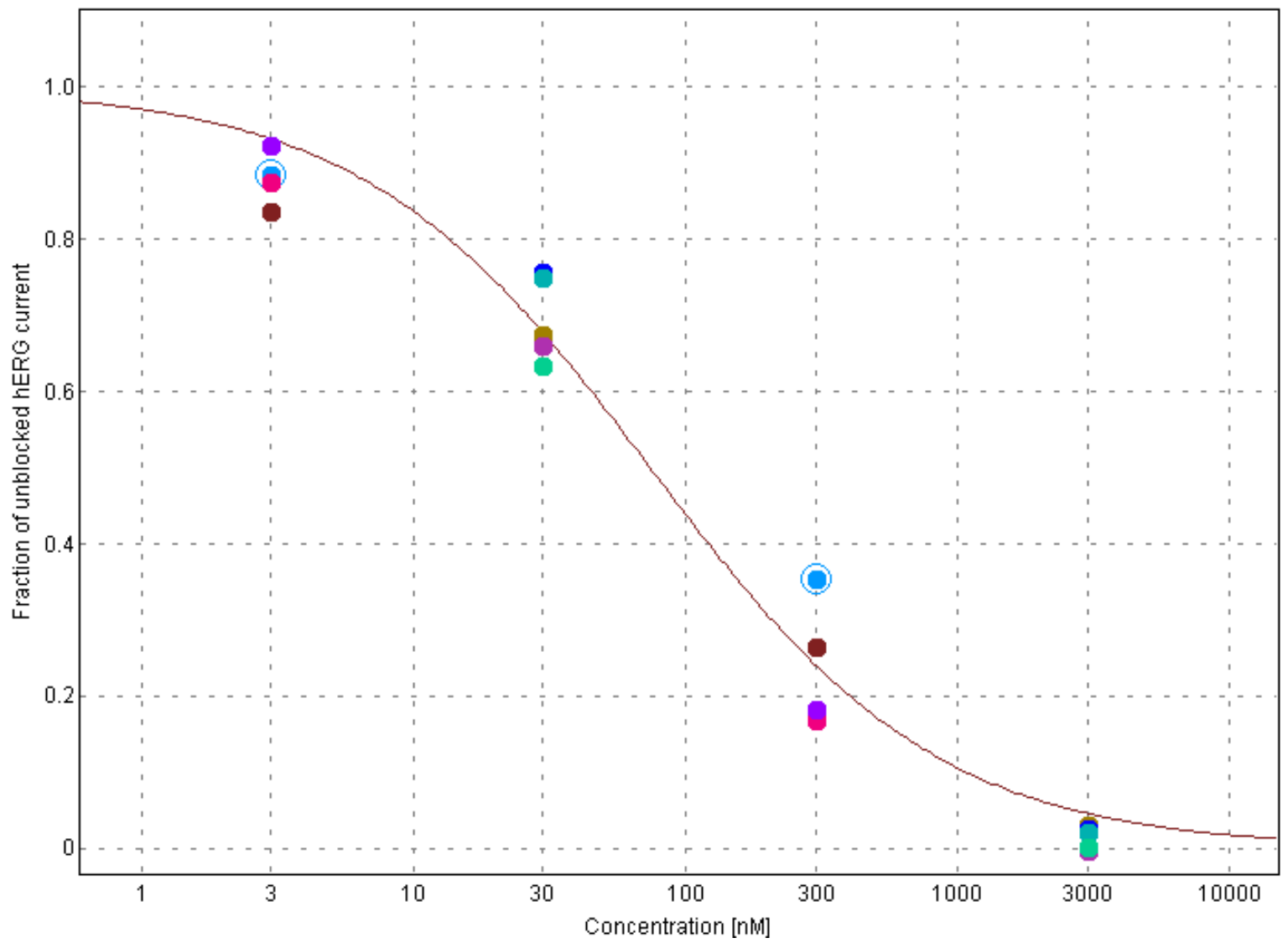


Figure 4.TIF

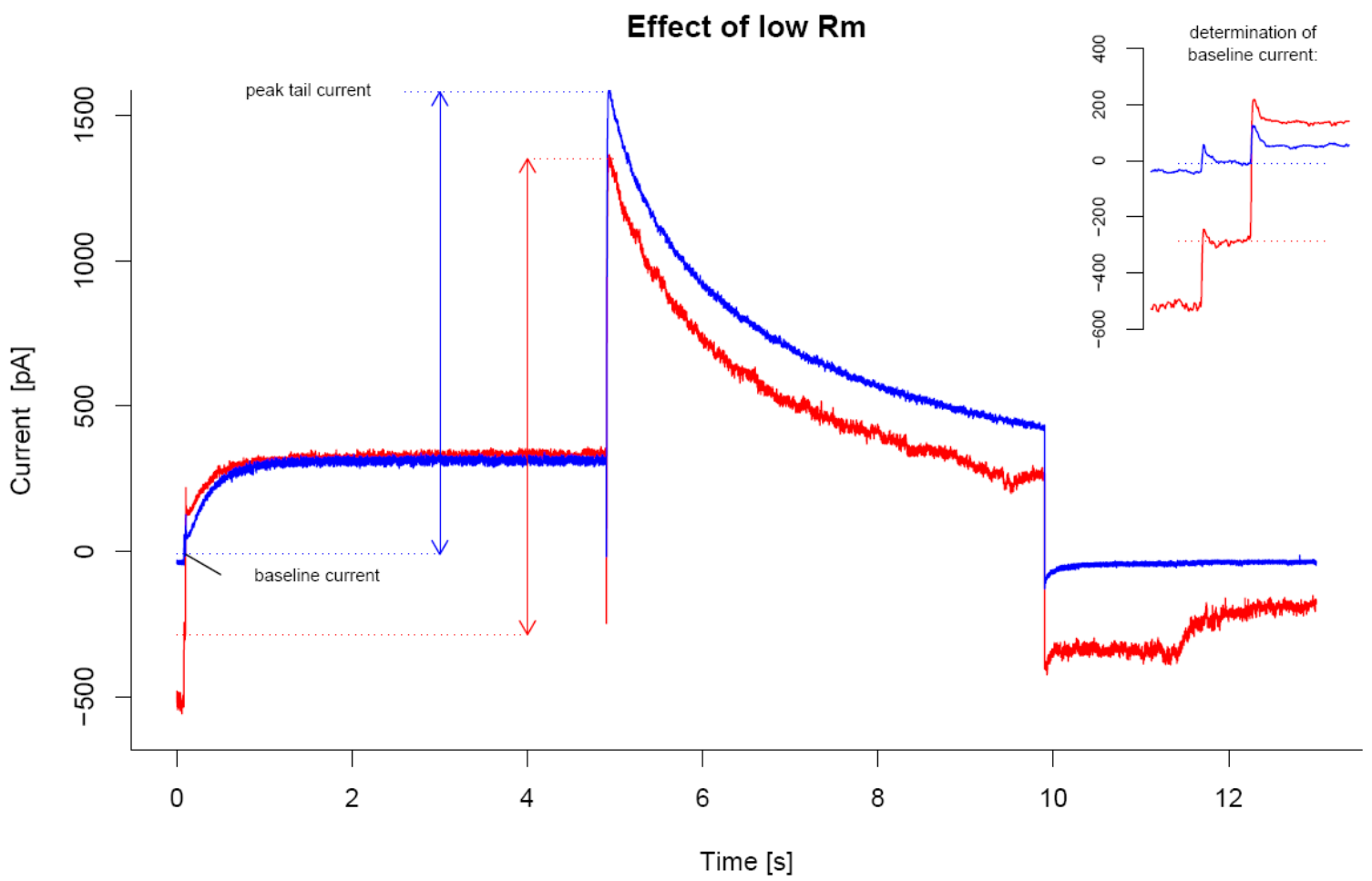


Figure 5.TIF

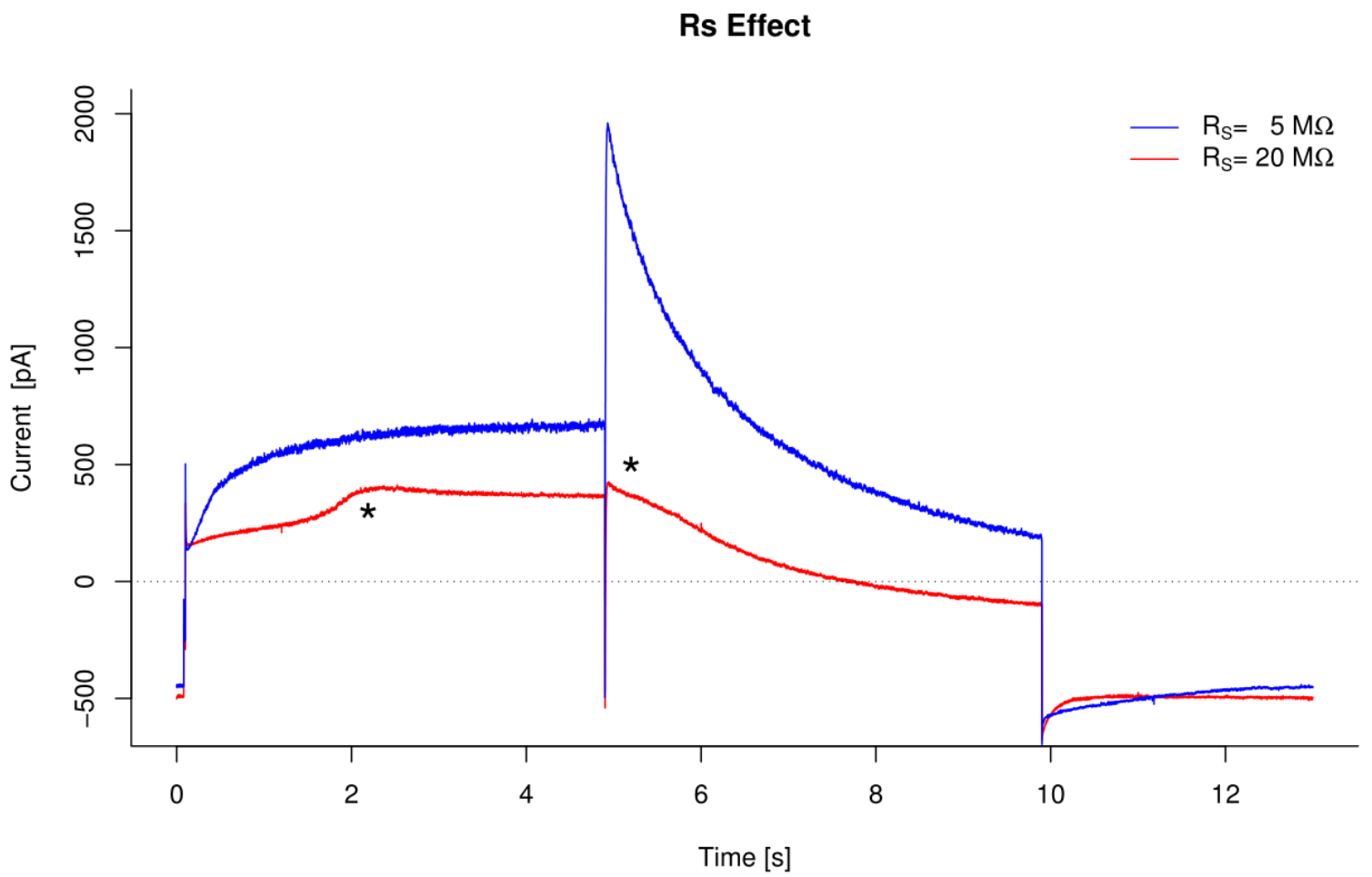
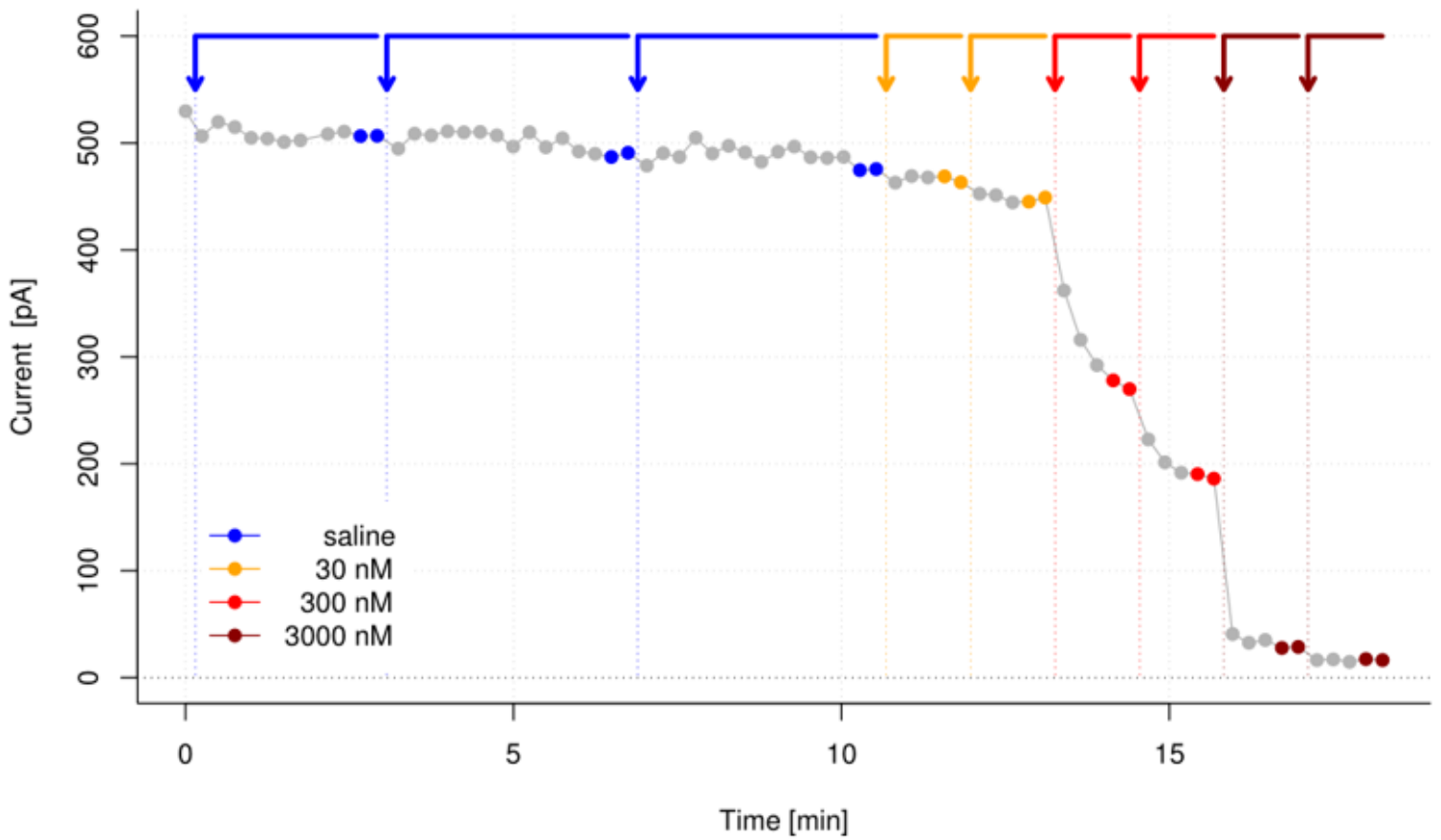


Figure 6.TIF

A Terfenadine: incomplete steady state



B Terfenadine: acceptable steady state

