

Electrophysiological recordings of CNT1 (SLC28A1) activity on Nanion's SURFE²R N1

The electrophysiology team at Nanion Technologies GmbH, Munich.

Summary

The concentrative nucleoside transporter 1 (CNT1) is a sodium-dependent uptake transporter encoded by the SLC28A1 gene¹. CNT1 functions as a co-transporter, coupling the uphill nucleoside transport into the cells to the electrochemical gradient of sodium². The stoichiometry of transport is proposed to be 1:1³, but a stoichiometry of 2 Na⁺: 1 nucleoside has also been suggested⁴. CNT1 is an electrogenic transporter, generating a net charge flow. It plays a major role in the uptake of pyrimidines, including uridine and cytidine, from the extracellular milieu into the cytoplasm¹ in nucleoside salvage pathways which is the first step of nucleoside biosynthesis². The transporter is expressed in epithelial tissues including liver, kidney and small intestine where it is localized to the apical membrane². CNTs are important targets for many antiviral and anticancer agents^{5,6}, and CNT1 has been proposed to play a role in tumor biology via a mechanism beyond nucleoside transport⁷. In fact, tumors expressing high levels of CNT1 can indicate a higher risk of relapse for breast cancer patients⁸, suggesting that nucleoside salvage may interfere with chemosensitivity⁸. On the other hand, high expression of the CNT1 protein could promote drug-induced cytotoxicity if patients were treated with suitable hCNT substrates¹. In any case, hCNT1 is an important mediator in the transport of anticancer and antiviral nucleoside drugs^{1,3,5,6} by mechanisms that require further study.

Here we present CNT1 activity measurements on the SURFE²R N1 instrument using purified plasma membrane of CHO cells expressing CNT1.

Download more Application Notes from www.nanion.de

Results

To activate CNT1 on the SURFE²R N1, a sensor with attached CNT1-containing membrane fragments was inserted into the device and perfused with a buffer containing NaCl and uridine. When the nucleoside is present, sodium movement across the membrane can be observed until an electrochemical equilibrium is reached. To generate sodium gradients, necessary as a driving force, the sensor was flushed with KCl before and after activation of CNT1 (Figure 1). Uridine elicits a larger response of CNT1 compared with cytidine but CNT1 has a higher apparent affinity for cytidine vs uridine (Figure 2).

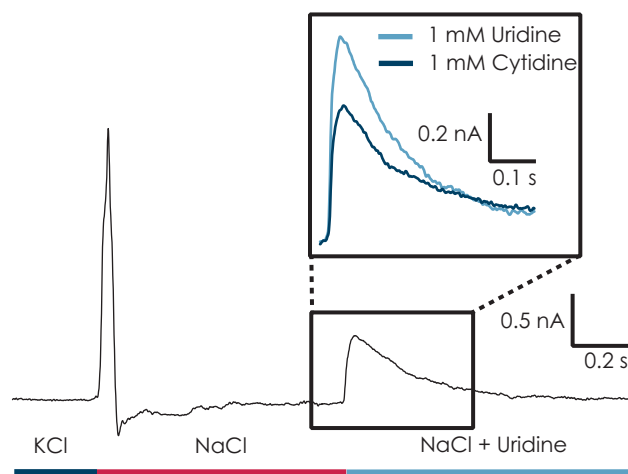


Figure 1: Typical CNT1 current response on the SURFE²R N1. Uridine was used as a substrate for CNT1. After establishing a sodium gradient (first peak), uridine was applied to the sensor (second peak). Uridine or cytidine can be used as the substrate, uridine eliciting a larger response than cytidine (inset).

Application Note

The apparent affinity of CNT1 to the substrates uridine and cytidine was investigated. Apparent K_m values of $78.67 \pm 21.16 \mu\text{M}$ ($n = 6$) for uridine, and $8.24 \pm 1.38 \mu\text{M}$ ($n = 5$) for cytidine were determined (Figure 2) in good agreement with range found in the literature^{3,4,9,10}.

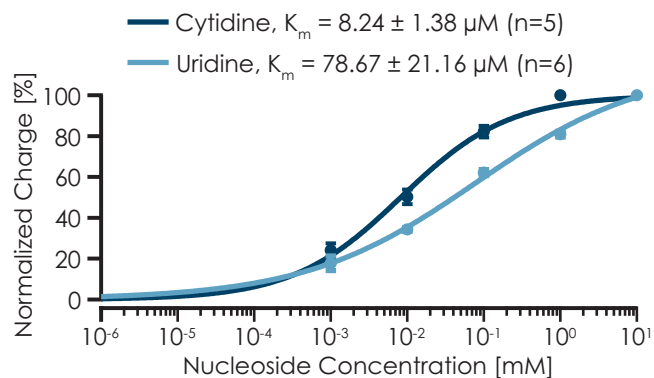


Figure 2: Increasing concentrations of 2 different substrates were added cumulatively on the same population of membrane fragments and the average concentration response curves for cytidine ($n = 5$) and uridine ($n = 6$) were constructed. When fitted with a Hill equation, apparent K_m values of $8.24 \pm 1.38 \mu\text{M}$ ($n = 5$) and $78.67 \pm 21.16 \mu\text{M}$ ($n = 6$) for cytidine and uridine, respectively, were obtained.

The effect of different ion gradients on CNT1 activity was also investigated. Figure 3 shows charge values for CNT1 under different ionic conditions: Na^+ , K^+ ,

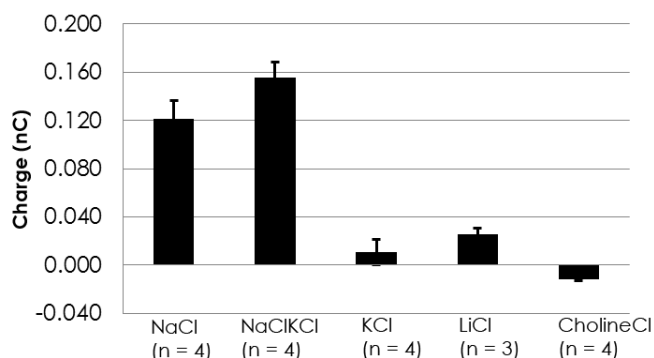


Figure 3: CNT1 co-transporters Na^+ , Li^+ acts as a weak substrate, K^+ and choline cannot be co-transported by CNT1. A Na^+ gradient (NaCl/KCl) increases the signal amplitude.

Li^+ and choline were used as the co-transported ion in symmetrical conditions. Significant ionic transfer by CNT1 only occurs in the presence of Na^+ , demonstrating strict coupling of nucleotide and sodium transport and a high Na^+ specificity of the transporter. Application of a sodium gradient (NaCl/KCl) increases the current amplitude, illustrating the resulting increase of driving force.

In conclusion, the SURFE²R N1 can be used to reliably measure CNT1 activity. This has important implications for drug discovery targeting CNT1 because nucleoside analogues are used as anticancer and antiviral therapies.

References

1. Pastor-Anglada, M., et al., 2008. *Xenobiotica*. 38(7-8):972-994
2. Gray, J.H., et al., 2004. *Pflügers Arch*. 447:728-734
3. Smith, K.M., et al., 2004. *J. Physiol*. 558.3:807-823
4. Larráyoz, I.M., et al., 2004. *JBC*. 279(10): 8999-9007
5. Pastor-Anglada, M., et al., 1998. *TIPS*. 19:424-430
6. Pastor-Anglada, M., et al., 2005. *Vir. Res*. 107:151-164
7. Pérez-Torras, S., et al., 2013. *Cell Death and Disease* 4, e648; doi:10.1038/cd-dis.2013.173
8. Gloeckner-Hofmann, K., 2006. *Oncology*. 70:238-244
9. Graham, K.A., et al., 2000. *Nucleosides Nucleotides Nucleic Acids*. 19(1-2):415-34
10. Griffith, D.A., & Jarvis, S.M. 1996. *Biochimica et Biophysica Acta* 1286:153-181

Methods

Plasma membrane preparation

According to the Nanion's standard procedure ("Quickguide Membrane Preparation from CHO cells"). Total protein concentration was between 5 - 10 $\mu\text{g}/\mu\text{l}$.

Buffers

CNT1 experiments were performed by the exchange of a sodium and nucleoside free ("resting") buffer for a sodium containing ("control") buffer and afterwards a nucleoside and sodium containing ("activating") buffer. Resting buffer contained: 140 mM KCl, 5 mM MgCl_2 , 30 mM HEPES, pH 7.4 with NMG. Activating buffer contained: 140 mM NaCl, 5 mM MgCl_2 , 30 mM HEPES, pH 7.4 with NMG, x mM uridine/cytidine.

SURFE²R sensor preparation

According to the Nanion standard procedure "SURFE²R Sensor Preparation". Sensors are prepared in resting buffer, membrane is diluted 1:10 with resting buffer.

SURFE²R N1 measurement workflow

CNT1 can be activated by providing uridine and cytidine as a nucleoside. A sodium gradient must be established in advance of nucleoside addition. Therefore, any 3-buffer Nanion standard protocol is suitable.