



nanjion

SOT

56th
Annual Meeting
and ToxExpo™

Baltimore,
Maryland

Axiogenesis & Nanion: Exhibitor-Hosted Session

Monday, March 13,
10:30 AM – 11.30 AM, Room 337

Title:

Mechanistic Insights into
Contractility and Electrophysiology
Complemented by Optical
Stimulation of iPS Cardiomyocytes

nanjion

axio GENESIS



Stem Cell Networking Reception

Monday, March 13,

5:00 PM – 8:00 PM

Pratt Ale House, 206 W. Pratt St.