

## Electrophysiological recordings of CNT1 (SLC28A1) activity on Nanion's SURFE<sup>2</sup>R N1

The electrophysiology team at Nanion Technologies GmbH, Munich.

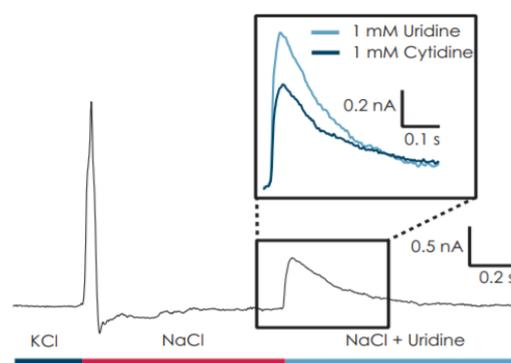
### Summary

The concentrative nucleoside transporter 1 (CNT1) is a sodium-dependent uptake transporter encoded by the SLC28A1 gene<sup>1</sup>. CNT1 functions as a co-transporter, coupling the uphill nucleoside transport into the cells to the electrochemical gradient of sodium<sup>2</sup>. The stoichiometry of transport is proposed to be 1:1<sup>3</sup>, but a stoichiometry of 2 Na<sup>+</sup>: 1 nucleoside has also been suggested<sup>4</sup>. 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 cytoplasm<sup>1</sup> in nucleoside salvage pathways which is the first step of nucleoside biosynthesis<sup>2</sup>. The transporter is expressed in epithelial tissues including liver, kidney and small intestine where it is localized to the apical membrane<sup>2</sup>. CNTs are important targets for many antiviral and anticancer agents<sup>5,6</sup>, and CNT1 has been proposed to play a role in tumor biology via a mechanism beyond nucleoside transport<sup>7</sup>. In fact, tumors expressing high levels of CNT1 can indicate a higher risk of relapse for breast cancer patients<sup>8</sup>, suggesting that nucleoside salvage may interfere with chemosensitivity<sup>8</sup>. On the other hand, high expression of the CNT1 protein could promote drug-induced cytotoxicity if patients were treated with suitable hCNT substrates<sup>1</sup>. In any case, hCNT1 is an important mediator in the transport of anticancer and antiviral nucleoside drugs<sup>1,3,5,6</sup> by mechanisms that require further study.

Here we present CNT1 activity measurements on the SURFE<sup>2</sup>R N1 instrument using purified plasma membrane of CHO cells expressing CNT1.

### Results

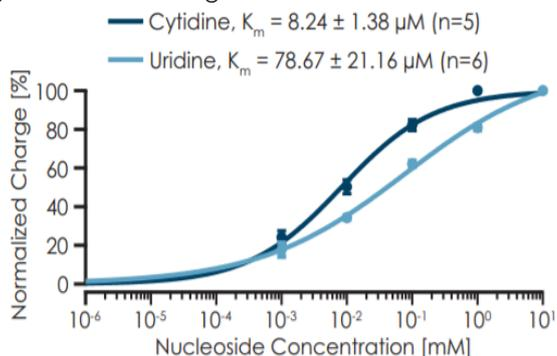
To activate CNT1 on the SURFE<sup>2</sup>R N1, a sensor with attached CNT1-containing membrane fragments was inserted into the device and perfused with a buffer containing NaCl and uridine. When the nucleoside is present, sodium movement across the membrane can be observed until an electrochemical equilibrium 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 SURFE<sup>2</sup>R 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).

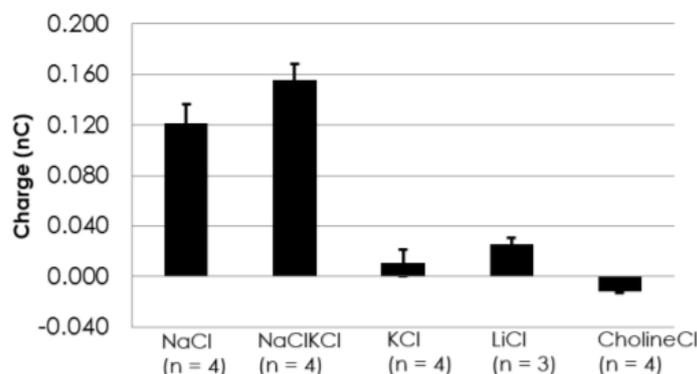
# Application Note

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<sup>3,4,9,10</sup>.



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



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

In conclusion, the SURFE<sup>2</sup>R 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/cddis.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

Nanion Technologies GmbH  
Ganghoferstr. 70A  
80339 Munich, Germany

phone +49 89 219 095-0  
fax +49 89 218997960  
www.nanion.de • info@nanion.de

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  $\text{MgCl}_2$ , 30mM HEPES, pH 7.4 with NMG. Activating buffer contained: 140 mM NaCl, 5 mM  $\text{MgCl}_2$ , 30 mM HEPES, pH 7.4 with NMG, x mM uridine/cytidine.

### SURFE<sup>2</sup>R sensor preparation

According to the Nanion standard procedure "SURFE<sup>2</sup>R Sensor Preparation". Sensors are prepared in resting buffer, membrane is diluted 1:10 with resting buffer.

### SURFE<sup>2</sup>R 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.