

# PRODUCT INFORMATION



## <sup>10</sup>Panx (trifluoroacetate salt)

Item No. 36867

**Formal Name:** L-tryptophyl-L-arginyl-L-glutamyl-L-alanyl-L-alanyl-L-phenylalanyl-L-valyl-L-α-aspartyl-L-seryl-L-tyrosine, trifluoroacetate salt

**Synonyms:** <sup>10</sup>Panx1, WRQAAFVDSY

**MF:** C<sub>58</sub>H<sub>79</sub>N<sub>15</sub>O<sub>16</sub> • XCF<sub>3</sub>COOH

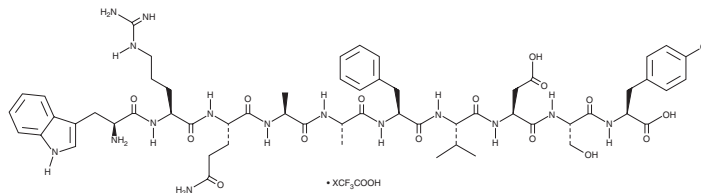
**FW:** 1,242.3

**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

<sup>10</sup>Panx (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the <sup>10</sup>Panx (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. <sup>10</sup>Panx (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of <sup>10</sup>Panx (trifluoroacetate salt) in these solvents is approximately 2 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of <sup>10</sup>Panx (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of <sup>10</sup>Panx (trifluoroacetate salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

<sup>10</sup>Panx is a peptide mimetic of the pannexin 1 (PANX1) ectodomain and inhibitor of PANX1, an innexin family gap domain protein and component of membrane channels.<sup>1</sup> It inhibits purinergic P2X<sub>7</sub> receptor activation-induced dye uptake, constitutive dye uptake, and hemichannel currents, as well as ATP-induced caspase activation and IL-1β release in HEK293 cells when used at a concentration of 200 μM. <sup>10</sup>Panx inhibits pannexin channel currents and connexin 46 channel currents in *Xenopus* oocytes expressing the human channels.<sup>2</sup>

### References

1. Pelegrin, P. and Surprenant, A. Pannexin-1 mediates large pore formation and interleukin-1β release by the ATP-gated P2X<sub>7</sub> receptor. *EMBO J.* **25**(21), 5071-5082 (2006).
2. Wang, J., Ma, M., Locovei, S., et al. Modulation of membrane channel currents by gap junction protein mimetic peptides: Size matters. *Am. J. Physiol. Cell Physiol.* **293**(3), C1112-C1119 (2007).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)