

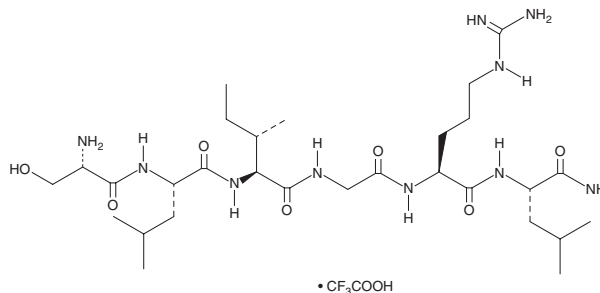
# PRODUCT INFORMATION



## SLIGRL-NH<sub>2</sub> (trifluoroacetate salt)

Item No. 16723

**Formal Name:** L-seryl-L-leucyl-L-isoleucylglycyl-L-arginyl-L-Leucinamide, trifluoroacetate salt  
**MF:** C<sub>29</sub>H<sub>56</sub>N<sub>10</sub>O<sub>7</sub> • XCF<sub>3</sub>COOH  
**FW:** 656.8  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

SLIGRL-NH<sub>2</sub> (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the SLIGRL-NH<sub>2</sub> (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SLIGRL-NH<sub>2</sub> (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SLIGRL-NH<sub>2</sub> (trifluoroacetate salt) is approximately 5 mg/ml in ethanol and approximately 20 mg/ml in DMSO and DMF.

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 SLIGRL-NH<sub>2</sub> can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of SLIGRL-NH<sub>2</sub> in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Proteinase-activated receptor 2 (PAR2) is a G protein-coupled receptor that is cleaved by serine proteases, resulting in self-activation of the receptor by a tethered ligand.<sup>1,2</sup> The ligand corresponds to residues 39-44 (SLIGRL in mouse PAR2).<sup>1,2</sup> SLIGRL-NH<sub>2</sub> is a recombinant peptide that activates PAR2 (EC<sub>50</sub> = ~5 μM), without requiring receptor cleavage.<sup>1,2</sup> This peptide does not activate PAR1. Through its effects on PAR2, SLIGRL-NH<sub>2</sub> stimulates gastric and intestinal smooth muscle contraction and induces thermal hyperalgesia in mice.<sup>3,4</sup> SLIGRL-NH<sub>2</sub> is used to explore signaling through PAR2 in cells and in animals.<sup>5,6</sup>

### References

1. Nystedt, S., Emilsson, K., Wahlestedt, C., *et al.* Molecular cloning of a potential proteinase activated receptor. *Proc. Natl. Acad. Sci. USA* **97**(20), 9208-9212 (1994).
2. Nystedt, S., Larsson, A.-K., Aberg, H., *et al.* The mouse proteinase-activated receptor-2 cDNA and gene. Molecular cloning and functional expression. *J. Biol. Chem.* **270**(11), 5950-5955 (1995).
3. Kawabata, A., Kuroda, R., Nagata, N., *et al.* *In vivo* evidence that protease-activated receptors 1 and 2 modulate gastrointestinal transit in the mouse. *Br. J. Pharmacol.* **133**(8), 1213-1218 (2001).
4. Liu, Q., Weng, H.-J., Patel, K.N., *et al.* The distinct roles of two GPCRs, MrgprC11 and PAR2, in itch and hyperalgesia. *Sci. Signal.* **4**(181), 1-6 (2011).
5. Sriwai, W., Mahavadi, S., Al-Shboul, O., *et al.* Distinctive G protein-dependent signaling by protease-activated receptor 2 (PAR2) in smooth muscle: Feedback inhibition of RhoA by cAMP-independent PKA. *PLoS One* **8**(6), 1-12 (2013).
6. Elmariah, S.B., Reddy, V.B., and Lerner, E.A. Cathepsin S signals via PAR2 and generates a novel tethered ligand receptor agonist. *PLoS One* **9**(6), 1-9 (2014).

#### 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, 11/14/2022

#### 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