

# PRODUCT INFORMATION



## PAR3 (1-6) amide (mouse) (trifluoroacetate salt)

Item No. 27130

<b>Formal Name:</b>	L-seryl-L-phenylalanyl-L-asparaginylglycylglycyl-L-prolinamide, trifluoroacetate salt	
<b>Synonyms:</b>	H-Ser-Phe-Asn-Gly-Gly-Pro-NH <sub>2</sub> , SFNGGP amide, SFNGGP-NH <sub>2</sub> , SFN-NH <sub>2</sub>	
<b>MF:</b>	C <sub>25</sub> H <sub>36</sub> N <sub>8</sub> O <sub>8</sub> • XCF <sub>3</sub> COOH	H—Ser—Phe—Asn—Gly—Gly—Pro—NH <sub>2</sub>
<b>FW:</b>	576.6	• XCF <sub>3</sub> COOH
<b>Purity:</b>	≥95%	
<b>Supplied as:</b>	A solid	
<b>Storage:</b>	-20°C	
<b>Stability:</b>	≥4 years	

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

### Laboratory Procedures

PAR3 (1-6) amide (mouse) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PAR3 (1-6) amide (mouse) (trifluoroacetate salt) in water. The solubility of PAR3 (1-6) amide (mouse) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PAR3 (1-6) amide is a synthetic hexapeptide that corresponds to amino acid residues 1-6 of the amino terminal tethered ligand sequence of mouse proteinase-activated receptor 3 (PAR3) and residues 38-43 of the full-length sequence. PAR3 (1-6) amide activates PAR1 and PAR2, but not PAR3 or PAR4.<sup>1,2</sup> It does not induce thrombin signaling in COS-7 cells expressing PAR3, PAR4, or both receptors at concentrations less than or equal to 500 μM, however, it does activate PAR1 and PAR2 in a cross-desensitization assay in Jurkat T cells when used at concentrations of 500 and 1,000 μM.<sup>1,2</sup> PAR3 (1-6) increases calcium levels in Kirsten virus-transformed rat kidney (KNRK) cells expressing rat or human PAR2, with a higher potency for rat PAR2.

### References

1. Nakanishi-Matsui, M., Zheng, Y.W., Sulciner, D.J., *et al.* PAR3 is a cofactor for PAR4 activation by thrombin. *Nature* **404(6778)**, 609-613 (2000).
2. Hansen, K.K., Saifeddine, M., and Hollenberg, M.D. Tethered ligand-derived peptides of proteinase-activated receptor 3 (PAR3) activate PAR1 and PAR2 in Jurkat T cells. *Immunology* **112(2)**, 183-190 (2004).

#### 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