

PRODUCT INFORMATION

PAR3 (1-6) amide (human) (trifluoroacetate salt)

Item No. 28201

Formal Name:	L-threonyl-L-phenylalanyl-L-arginylglycyl-L-alanyl-L-prolinamide, trifluoroacetate salt	
Synonyms:	Proteinase-Activated Receptor 3, TFRGAP-NH ₂	
MF:	C ₂₉ H ₄₆ N ₁₀ O ₇ • XCF ₃ COOH	H—Thr—Phe—Arg—Gly—Ala—Pro—NH ₂
FW:	646.7	• XCF ₃ COOH
Purity:	≥95%	
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

PAR3 (1-6) amide (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PAR3 (1-6) amide (human) (trifluoroacetate salt) in water. The solubility of PAR3 (1-6) amide (human) (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 peptide agonist of proteinase-activated receptor 1 (PAR1) and PAR2 that corresponds to residues 1-6 of the amino terminal tethered ligand sequence of human PAR3 and residues 39-44 of the full-length human sequence. PAR3 (1-6) amide activates PAR1 and PAR2 in Jurkat and HEK cells, inducing calcium accumulation in Jurkat cells, which is reduced following desensitization with the PAR1- and PAR2-selective peptide agonist SFLR-NH₂.¹ It decreases the expression of CDC42 in wild-type, but not PAR3 knockdown, PANC-1 cells and reduces invasion of PANC-1 cells.²

References

1. 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).
2. Segal, L., Katz, L.S., Lupu-Meiri, M., et al. Proteinase-activated receptors differentially modulate *in vitro* invasion of human pancreatic adenocarcinoma PANC-1 cells in correlation with changes in the expression of CDC42 protein. *Pancreas* **43**(1), 103-108 (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, 12/05/2022

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