

PRODUCT INFORMATION



PAR4 (1-6) amide (human) (trifluoroacetate salt)

Item No. 27127

Formal Name:	glycyl-L-tyrosyl-L-prolylglycyl-L-glutaminyll-L-valinamide, trifluoroacetate salt	
Synonyms:	GQV-NH ₂ , GYPGQV-NH ₂ , H-Gly-Tyr-Pro-Gly-Gln-Val-NH ₂	
MF:	C ₂₈ H ₄₂ N ₈ O ₈ • XCF ₃ COOH	H—Gly—Tyr—Pro—Gly—Gln—Val—NH ₂
FW:	618.7	• 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

PAR4 (1-6) amide (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PAR4 (1-6) amide (human) (trifluoroacetate salt) in water. The solubility of PAR4 (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

PAR4 (1-6) amide is a synthetic peptide agonist of proteinase-activated receptor 4 (PAR4) that corresponds to residues 1-6 of the amino terminal tethered ligand sequence of human PAR4 and residues 48-53 of the full-length sequence.¹ It induces contractions in isolated rat gastric longitudinal muscle tissue and endothelium-dependent relaxation of precontracted rat aortic rings (EC₅₀S = 300-400 μM).

References

1. Hollenberg, M.D., Saifedine, M., Al-Ani, B., *et al.* Proteinase-activated receptor 4 (PAR4): Action of PAR4-activating peptides in vascular and gastric tissue and lack of cross-reactivity with PAR1 and PAR2. *Can. J. Physiol. Pharmacol.* **77**(6), 458-464 (1999).

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