

PRODUCT INFORMATION



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

Item No. 27128

Formal Name: glycyl-L-tyrosyl-L-prolylglycyl-L-lysyl-L-phenylalaninamide, trifluoroacetate salt
MF: $C_{33}H_{46}N_8O_7 \cdot XCF_3COOH$
FW: 666.8
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years

$H-Gly-Tyr-Pro-Gly-Lys-Phe-NH_2$
 $\cdot XCF_3COOH$

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

Laboratory Procedures

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

PAR4 (1-6) amide is a synthetic peptide agonist of proteinase-activated receptor 4 (PAR4) and an amide form of PAR4 (1-6) (Item No. 27131) that corresponds to residues 1-6 of the amino terminal tethered ligand sequence of mouse PAR4 and residues 60-65 of the full-length sequence. It induces aggregation of rat platelets with an EC_{50} value of $40 \mu M$.¹ PAR4 (1-6) amide ($100 \mu M$) induces smooth muscle relaxation and prostaglandin E_2 (PGE_2 ; Item No. 14010) release in isolated mouse tracheal segments, effects that are inhibited by indomethacin (Item No. 70270).²

References

- Hollenberg, M.D. and Saifeddine, M. Proteinase-activated receptor 4 (PAR4): Activation and inhibition of rat platelet aggregation by PAR4-derived peptides. *Can. J. Physiol. Pharmacol.* **79(5)**, 439-442 (2001).
- Lan, R.S., Knight, D.A., Stewart, G.A., et al. Role of PGE_2 in protease-activated receptor-1, -2 and -4 mediated relaxation in the mouse isolated trachea. *Br. J. Pharmacol.* **132(1)**, 93-100 (2001).

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