

PRODUCT INFORMATION



(Ala1)-PAR4 (1-6) (mouse) (trifluoroacetate salt)

Item No. 27132

Formal Name: L-alanyl-L-tyrosyl-L-prolylglycyl-L-lysyl-L-phenylalanine, trifluoroacetate salt

Synonyms: H-Ala-Tyr-Pro-Gly-Lys-Phe-OH, AYPGKF

MF: $C_{34}H_{47}N_7O_8 \cdot XCF_3COOH$ H—Ala—Tyr—Pro—Gly—Lys—Phe—OH

FW: 681.8 • XCF₃COOH

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

(Ala1)-PAR4 (1-6) (mouse) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the (Ala1)-PAR4 (1-6) (mouse) (trifluoroacetate salt) in water. The solubility of (Ala1)-PAR4 (1-6) (mouse) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(Ala1)-PAR4 (1-6) 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 mouse PAR4 and residues 60-65 of the full-length sequence with an alanine residue substituted for glycine at position 1/60.¹ It is 10-fold more potent than PAR4 (1-6) (Item No. 27131), triggering [³H]inositol phosphate release from KOLF cells expressing human PAR4 at levels comparable to 30 nM thrombin (Item No. 13188) when used at a concentration of 500 μM. (Ala1)-PAR4 (1-6) induces aggregation of human platelets.²

References

1. Faruqi, T.R., Weiss, E.J., Shapiro, M.J., *et al.* Structure-function analysis of protease-activated receptor 4 tethered ligand peptides. *J. Biol. Chem.* **275(26)**, 19728-19734 (2000).
2. Lidington, E.A., Steinberg, R., Kinderlerer, A.R., *et al.* A role for proteinase-activated receptor 2 and PKC-ε in thrombin-mediated induction of decay-accelerating factor on human endothelial cells. *Am. J. Physiol. Cell Physiol.* **289(6)**, 1437-1447 (2005).

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, 12/01/2022

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