

PRODUCT INFORMATION

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

Item No. 27131

Formal Name:	glycyl-L-tyrosyl-L-prolylglycyl-L-lysyl-L-phenylalanine, trifluoroacetate salt	
Synonyms:	GYPGKF, H-Gly-Tyr-Pro-Gly-Lys-Phe-OH	
MF:	$C_{33}H_{45}N_7O_8 \cdot XCF_3COOH$	H—Gly—Tyr—Pro—Gly—Lys—Phe—OH
FW:	667.8	• XCF_3COOH
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) (mouse) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PAR4 (1-6) (mouse) (trifluoroacetate salt) in water. The solubility of 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

PAR4 (1-6) is a synthetic peptide agonist of proteinase-activated receptor 4 (PAR4; $EC_{50} = 30 \mu M$ in *Xenopus* oocytes expressing human PAR4 receptors) 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.¹ It induces calcium release in wild-type and *PAR3*^{-/-} mouse platelets and in human platelets but has no activity in *Xenopus* oocytes expressing human PAR1. PAR4 (1-6) also triggers [³H]inositol phosphate release from KOLF cells expressing human PAR4.²

References

1. Kahn, M.L., Zheng, Y.-W., Huang, W., *et al.* A dual thrombin receptor system for platelet activation. *Nature* **394**(6694), 690-694 (1998).
2. 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).

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, 10/27/2022

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