

PRODUCT INFORMATION



Hemokinin 1 (human) (trifluoroacetate salt)

Item No. 35458

Formal Name: L-threonylglycyl-L-lysyl-L-alanyl-L-seryl-L-glutamyl-L-phenylalanyl-L-phenylalanylglycyl-L-leucyl-L-methioninamide, trifluoroacetate salt

Synonyms: HK-1, Tachykinin-4 (57-67), Thr-Gly-Lys-Ala-Ser-Gln-Phe-Phe-Gly-Leu-Met-NH₂

Peptide Sequence: TGKASQFFGLM-NH₂

MF: C₅₄H₈₄N₁₄O₁₄S • XCF₃COOH

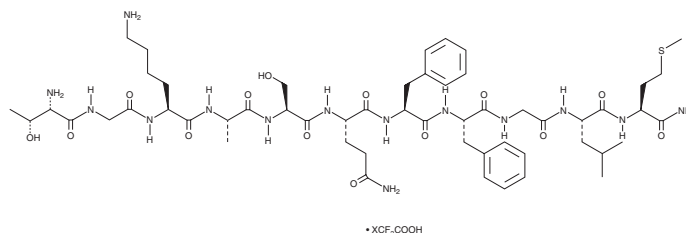
FW: 1,185.4

Purity: ≥98%

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

Hemokinin 1 (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the hemokinin 1 (human) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Hemokinin 1 is a tachykinin peptide neurokinin-1 (NK₁) receptor agonist.¹⁻⁵ It induces migration of A375 and B16/F10 melanoma cells when used at a concentration of 1 μM, an effect that can be reversed by the NK₁ receptor antagonist L-732,138 (Item No. 30894).¹ Hemokinin 1 (0.001-0.1 μM) promotes the proliferation, migration, and tube formation of human umbilical vein endothelial cells (HUVECs).² It induces the production of prostaglandin E₂ (PGE₂; Item No. 14010) in human colon mucosal and muscle explants.³ Intracerebroventricular administration of hemokinin 1 increases the latency to withdrawal in the tail-flick test in mice in a dose-dependent manner.⁴ It decreases mean arterial pressure (MAP) and heart rate in anesthetized rats when administered at doses ranging from 1 to 30 nmol/kg.⁵

References

1. Zhang, Y., Li, X., Li, J., et al. *Peptides* **83**, 8-15 (2016).
2. Song, H., Yin, W., Zeng, Q., et al. *Int. J. Biochem. Cell Biol.* **44**(9), 1410-1421 (2012).
3. Dai, L., Perera, D.S., King, D.W., et al. *J. Pharmacol. Exp. Ther.* **340**(1), 27-36 (2012).
4. Fu, C.Y., Zhao, Y.L., Dong, L., et al. *Brain Behav. Immun.* **22**(6), 850-860 (2008).
5. Kong, Z.-Q., Fu, C.-Y., Chen, Q., et al. *Eur. J. Pharmacol.* **590**(1-3), 310-316 (2008).

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, 02/09/2026

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