

PRODUCT INFORMATION



Kinetensin (human) (trifluoroacetate salt)

Item No. 33138

Formal Name: L-isoleucyl-L-alanyl-L-arginyl-L-arginyl-L-histidyl-L-prolyl-L-tyrosyl-L-phenylalanyl-L-leucine, trifluoroacetate salt

Synonyms: IARRHPYFL, Ile-Ala-Arg-Arg-His-Pro-Tyr-Phe-Leu

MF: C₅₆H₈₅N₁₇O₁₁ • XCF₃COOH

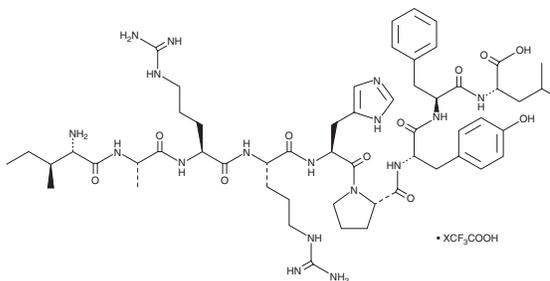
FW: 1,172.4

Purity: ≥95%

Supplied as: A crystalline 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

Kinetensin (human) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the kinetensin (human) (trifluoroacetate salt) in water. The solubility of kinetensin (human) (trifluoroacetate salt) in water is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Kinetensin is a nonapeptide that has been found in pepsin-treated human plasma.¹ It binds to neurotensin receptors in HT-29 cells and mouse brain ($K_{iS} = 1,396$ and 149.6 nM, respectively, in radioligand binding assays) but does not induce the accumulation of intracellular calcium in HT-29 cells ($EC_{50} = >10,000$ nM).² It induces histamine release in isolated rat peritoneal mast cells with an EC_{50} value of approximately 10 μ M.³ Intradermal administration of kinetensin induces vascular permeability in rats.

References

1. Mogard, M.H., Kobayashi, R., Chen, C.F., *et al.* The amino acid sequence of kinetensin, a novel peptide isolated from pepsin-treated human plasma: Homology with human serum albumin, neurotensin and angiotensin. *Biochem. Biophys. Res. Commun.* **136(3)**, 983-988 (1986).
2. Akunne, H.C., Demattos, S.B., Whetzel, S.Z., *et al.* Agonist properties of a stable hexapeptide analog of neurotensin, N^oMeArg-Lys-Pro-Trp-tLeu-Leu (NT1). *Biochem. Pharmacol.* **49(8)**, 1147-1154 (1995).
3. Sydbom, A., Ware, J., and Mogard, M.H. Stimulation of histamine release by the peptide kinetensin. *Agents Actions* **27(1-2)**, 68-71 (1989).

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, 08/01/2023

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