

PRODUCT INFORMATION



Scyliorhinin I (dogfish) (trifluoroacetate salt)

Item No. 43081

Formal Name: L-alanyl-L-lysyl-L-phenylalanyl-L- α -aspartyl-L-lysyl-L-phenylalanyl-L-tyrosylglycyl-L-leucyl-L-methioninamide, trifluoroacetate salt

Synonym: Ala-Lys-Phe-Asp-Lys-Phe-Tyr-Gly-Leu-Met-NH₂

Peptide Sequence: AKFDKIFYGLM-NH₂

MF: C₅₉H₈₇N₁₃O₁₃S • XCF₃COOH

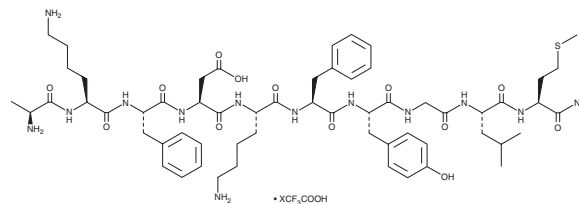
FW: 1,218.5

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

Scyliorhinin I (dogfish) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the scyliorhinin I (dogfish) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Scyliorhinin I (dogfish) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of scyliorhinin I (dogfish) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Scyliorhinin I (dogfish) (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Scyliorhinin I is a neurokinin-1 (NK₁) and NK₂ receptor peptide agonist that has been found in dogfish intestines.^{1,2} It binds to membrane preparations of rat submandibular gland and hamster urinary bladder (K_ds = 0.9 and 2 nM, respectively), tissues that endogenously express high levels of NK₁ and NK₂ receptors, respectively. It increases intracellular levels of inositol-1,4,5-trisphosphate (IP₃) in KNRK rat kidney cells expressing the NK₁ receptor when used at a concentration of 1 μM.² Scyliorhinin I induces contractions in isolated and perfused carp intestinal bulb smooth muscle strips (EC₅₀ = 15 nM).³ Intrathecal or intracerebroventricular administration of scyliorhinin I induces reciprocal hindlimb neck-scratching behavior in mice (ED₅₀s = 0.07 and 0.05 nmol/animal, respectively).⁴

References

1. Buck, S.H. and Krstenansky, J.L. The dogfish peptides scyliorhinin I and scyliorhinin II bind with differential selectivity to mammalian tachykinin receptors. *Eur. J. Pharmacol.* **144**(1), 109-111 (1987).
2. Vigna, S.R. The role of the amino-terminal domain of tachykinins in neurokinin-1 receptor signaling and desensitization. *Neuropeptides* **37**(1), 30-35 (2003).
3. Kitazawa, T. Excitatory responses to scyliorhinins I and II in smooth muscle strips isolated from the carp intestinal bulb (*Cyprinus carpio*). *Naunyn Schmiedebergs Arch Pharmacol.* **343**(5), 525-531 (1991).
4. Raffa, R.B., Martinez, R.P., and Connelly, C.D. Scyliorhinin-I and -II induce reciprocal hindlimb scratching in mice: Differentiation of spinal and supraspinal neurokinin receptors in vivo. *Neurosci. Lett.* **158**(1), 87-91 (1993).

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, 05/08/2025

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