

PRODUCT INFORMATION



Parasin I (catfish) (trifluoroacetate salt)

Item No. 36724

Formal Name: L-lysylglycyl-L-arginylglycyl-L-lysyl-L-glutaminyglycylglycyl-L-lysyl-L-valyl-L-arginyl-L-alanyl-L-lysyl-L-alanyl-L-lysyl-L-threonyl-L-arginyl-L-seryl-L-serine, trifluoroacetate salt

Peptide Sequence: KGRGKQGKVRRAKAKTRSS-OH

MF: $C_{82}H_{154}N_{34}O_{24} \cdot XCF_3COOH$

FW: 2,000.3

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

H — Lys — Gly — Arg — Gly — Lys — Gln — Gly — Gly — Lys — Val —
Arg — Ala — Lys — Ala — Lys — Thr — Arg — Ser — Ser — OH
• XCF_3COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Description

Parasin I is an antimicrobial peptide fragment from the N-terminal region of histone H2A in catfish (*P. asotus*) that is involved in host defense.¹ It is formed from histone H2A by cathepsin D in the skin mucosa of wounded *P. asotus*. Parasin I is active against various Gram-positive and Gram-negative bacteria, as well as fungi (MICs = 1-2, 1-4, and 1-2 $\mu g/ml$, respectively).²

References

1. Cho, J.H., Park, I.Y., and Kim, S.C. Cathepsin D produces the potent antimicrobial peptide parasin I from histone H2A in scaleless fish skin. *Peptides: The wave of the future*. Lebl, M. Houghten, R.A., editors, Springer (2001).
2. Park, I.Y., Park, C.B., Kim, M.S., et al. Parasin I, an antimicrobial peptide derived from histone H2A in the catfish, *Parasilurus asotus*. *FEBS Lett.* **437**(3), 258-262 (1998).

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, 09/25/2023

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