

PRODUCT INFORMATION



Indolicidin (trifluoroacetate salt)

Item No. 35845

Formal Name:	L-isoleucyl-L-leucyl-L-prolyl-L-tryptophyl-L-lysyl-L-tryptophyl-L-prolyl-L-tryptophyl-L-tryptophyl-L-prolyl-L-tryptophyl-L-arginyl-L-argininamide, trifluoroacetate salt	H—Ile—Leu—Pro—Trp—Lys—Trp—Pro—Trp—Trp—Pro—Trp—Arg—Arg—NH ₂
Synonym:	ILPWKWPWWPWRR-NH ₂	• XCF ₃ COOH
MF:	C ₁₀₀ H ₁₃₂ N ₂₆ O ₁₃ • XCF ₃ COOH	
FW:	1,906.3	
Purity:	≥98%	
UV/Vis.:	λ _{max} : 272 nm	
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

Indolicidin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the indolicidin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Indolicidin (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of indolicidin (trifluoroacetate salt) in these solvents is approximately 2, 5, and 10 mg/ml, respectively.

Description

Indolicidin is an antimicrobial peptide.¹ It is active against multidrug-resistant isolates of enteroaggregative *E. coli* (MDR-EAEC; MICs = 32 μM for all). Indolicidin is also active against *S. cerevisiae*, *T. beigellii*, and *C. albicans* (MICs = 5-10, 2.5-5, and 5-10 μM, respectively) and reduces viral replication of HIV-1 in MT-2 cells when used at concentrations ranging from 67 to 100 μg/ml.^{2,3} *In vivo*, indolicidin increases survival of MDR-EAEC-infected *G. mellonella* larvae.¹

References

1. Vergis, J., Malik, S.S., Pathak, R., et al. Antimicrobial efficacy of indolicidin against multi-drug resistant enteroaggregative *Escherichia coli* in a *Galleria mellonella* model. *Front. Microbiol.* **10**, 2723 (2019).
2. Lee, D.G., Kim, H.K., Kim, S.A., et al. Fungicidal effect of indolicidin and its interaction with phospholipid membranes. *Biochem. Res. Commun.* **305(2)**, 305-310 (2003).
3. Robinson, W.E., Jr., McDougall, B., Tran, D., et al. Anti-HIV-1 activity of indolicidin, an antimicrobial peptide from neutrophils. *J. Leukoc. Biol.* **63(1)**, 94-100 (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, 12/15/2022

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