

PRODUCT INFORMATION



PR 26 (trifluoroacetate salt)

Item No. 37460

Synonym: Proline-arginine-rich 26
Peptide Sequence: RRRPRPPYLPRPRPPFFPPRLPPRI-OH H-Arg-Arg-Arg-Pro-Arg-Pro-Pro-Tyr-Leu-Pro-Arg-Pro-Arg-Pro-Pro-Pro-Phe-Phe-Pro-Pro-Arg-Leu-Pro-Pro-Arg-Ile-OH
MF: C₁₅₃H₂₄₂N₅₀O₂₈ • XCF₃COOH
FW: 3,229.9
Purity: ≥95%
Supplied as: A solid • XCF₃COOH
Storage: -20°C
Stability: ≥4 years

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

Laboratory Procedures

PR 26 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PR 26 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas.

PR 26 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Proline-arginine-rich 26 (PR 26) is a synthetic peptide fragment of the antibiotic peptide PR 39 peptide (Item No. 36717).¹ It is active against *E. coli*, *S. typhimurium*, *S. choleraesuis*, and *S. suis* (MICs = 2, 1, 2, and 1 µg/ml, respectively). PR 26 (3 µM) inhibits *S. typhimurium* infection of rat IEC-6 intestinal cells. It increases phagocytosis of *S. choleraesuis* in isolated porcine neutrophils when used at a concentration of 2 µM. PR 26 (1 µM) induces chemotaxis in isolated porcine neutrophils.²

References

1. Shi, J., Ross, C.R., Chengappa, M.M., *et al.* Antibacterial activity of a synthetic peptide (PR-26) derived from PR-39, a proline-arginine-rich neutrophil antimicrobial peptide. *Antimicrob Agents Chemother.* **40**(1), 115-121 (1996).
2. Huang, H.J., Ross, C.R., and Blecha, F. Chemoattractant properties of PR-39, a neutrophil antibacterial peptide. *J. Leukoc. Biol.* **61**(5), 624-629 (1997).

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, 03/19/2025

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