

PRODUCT INFORMATION



PapRIV (trifluoroacetate salt)

Item No. 37453

Formal Name: (S)-4-((S)-2-((S)-1-(L-seryl-L-aspartyl-L-leucyl)pyrrolidine-2-carboxamido)-3-phenylpropanamido)-5-(((S)-1-carboxy-2-(1H-imidazol-4-yl)ethyl)amino)-5-oxopentanoic acid

Peptide Sequence: SDLPFEH-OH

MF: C₃₈H₅₃N₉O₁₃ • XCF₃COOH

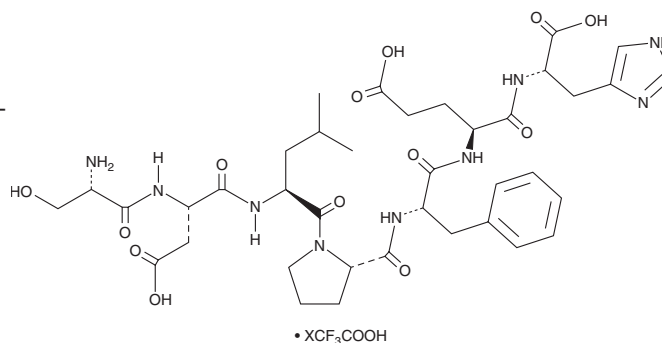
FW: 843.9

Purity: ≥98%

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

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

Description

PapRIV is a quorum-sensing heptapeptide originally isolated from *B. cereus*.¹ It is translated as a 48-amino acid polypeptide, which is secreted and extracellularly processed by NprB proteases to form the active heptapeptide.² PapRIV (1-25 μM) induces IL-6 and TNF-α production by, and NF-κB translocation in, BV-2 microglia cells.¹

References

1. Janssens, Y., Debonne, N., De Spiegeleer, A., *et al.* PapRIV, a BV-2 microglial cell activating quorum sensing peptide. *Sci. Rep.* **11(1)**, 10723 (2021).
2. Pomerantsev, A.P., Pomerantseva, O.M., Camp, A.S., *et al.* PapR peptide maturation: role of the NprB protease in *Bacillus cereus* 569 PlcR/PapR global gene regulation. *FEMS Immunol. Med. Microbiol.* **55(3)**, 361-377 (2009).

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, 10/04/2023

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