

PRODUCT INFORMATION



Z-VRPR-FMK (trifluoroacetate salt)

Item No. 19617

Formal Name: N-[(phenylmethoxy)carbonyl]-L-valyl-L-arginyl-N-
[[1S]-4-[(aminoiminomethyl)amino]-1-(2-fluoroacetyl)
butyl]-L-prolinamide, 2,2,2-trifluoroacetate

Synonym: Z-Val-Arg-Pro-DL-Arg-FMK

MF: C₃₁H₄₉FN₁₀O₆ • CF₃COOH

FW: 790.8

Purity: ≥90%

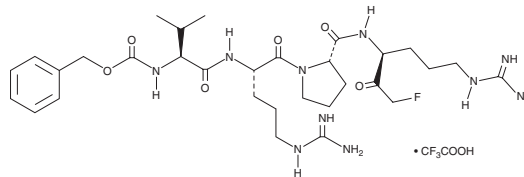
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Avoid freeze/thaw cycles. Hygroscopic. Protect from moisture.

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



Laboratory Procedures

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

Description

Z-VRPR-FMK is a selective, irreversible inhibitor of the paracaspase mucosa-associated lymphoid tissue lymphoma translocation protein 1 (MALT1) an endopeptidase that targets BCL10 and reduces NF-κB activation in lymphocytes, at 12.5 to 75 μM.¹⁻³ Other MALT1 substrates that may be affected by Z-VRPR-FMK are TNFAIP3 and CYLD.³

References

1. Rebeaud, F., Hailfinger, S., Posevitz-Fejfar, A., *et al.* The proteolytic activity of the paracaspase MALT1 is key in T cell activation. *Nat. Immunol.* **9(3)**, 272-281 (2008).
2. Hachmann, J., Snipas, S.J., van Raam, B.J., *et al.* Mechanism and specificity of the human paracaspase MALT1. *Biochem. J.* **443(1)**, 287-295 (2012).
3. Fernandez-Jimenez, N., Castellanos-Rubio, A., Plaza-Izurietta, L., *et al.* Coregulation and modulation of NFκB-related genes in celiac disease: Uncovered aspects of gut mucosal inflammation. *Hum. Mol. Genet.* **23(5)**, 1298-1310 (2014).

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, 11/29/2022

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