

PRODUCT INFORMATION

Boc-Val-Pro-Arg-AMC (trifluoroacetate salt)

Item No. 45711

Formal Name: *tert*-butyl ((S)-1-((S)-2-(((S)-5-guanidino-1-((4-methyl-2-oxo-2H-chromen-7-yl)amino)-1-oxopentan-2-yl)carbamoyl)pyrrolidin-1-yl)-3-methyl-1-oxobutan-2-yl)carbamate, 2,2,2-trifluoroacetic acid

Synonym: Boc-VPR-AMC

MF: C₃₁H₄₅N₇O₇ • CF₃COOH

FW: 741.8

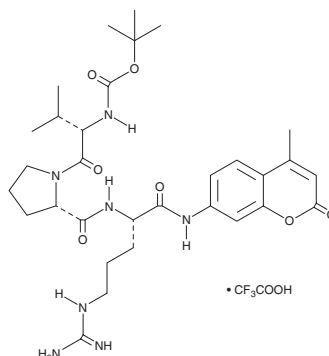
Purity: ≥95%

Ex./Em. Max: 340-360/440-460 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

Boc-Val-Pro-Arg-AMC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Boc-Val-Pro-Arg-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Boc-Val-Pro-Arg-AMC (trifluoroacetate salt) is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Boc-Val-Pro-Arg-AMC is a fluorogenic substrate for serine proteases.^{1,2} Upon enzymatic cleavage by serine proteases, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify serine protease activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

References

1. Michael, I.P., Sotiropoulou, G., Pampalakis, G., *et al.* Biochemical and enzymatic characterization of human kallikrein 5 (hK5), a novel serine protease potentially involved in cancer progression. *J. Biol. Chem.* **280**(15), 14628-14635 (2005).
2. Crossen, J., Shankar, K.N., and Diamond, S.L. Investigating thrombin-loaded fibrin in whole blood clot microfluidic assay via fluorogenic peptide. *Biophys. J.* **122**(4), 697-712 (2023).

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, 06/16/2026

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