

# PRODUCT INFORMATION



## Bz-FVR-AMC (trifluoroacetate salt)

Item No. 24845

**Formal Name:** N-((S)-1-(((S)-1-(((S)-5-guanidino-1-((4-methyl-2-oxo-2H-chromen-7-yl)amino)-1-oxopentan-2-yl)amino)-3-methyl-1-oxobutan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)benzamide, trifluoroacetate salt

**Synonyms:** Benzoyl-Phe-Val-Arg-AMC, Thrombin Fluorogenic Substrate III

**MF:** C<sub>37</sub>H<sub>43</sub>N<sub>7</sub>O<sub>6</sub> • XCF<sub>3</sub>COOH

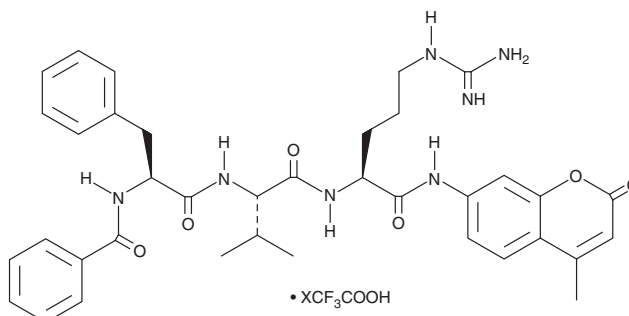
**FW:** 681.8

**Purity:** ≥95%

**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

Bz-FVR-AMC (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Bz-FVR-AMC (trifluoroacetate salt) in water. The solubility of Bz-FVR-AMC (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Bz-FVR-AMC is a fluorogenic peptide substrate for bovine cathepsins S and H, *Xenopus* endopeptidase XSECP1, and esterase E-II from *B. gabonica* venom, among others.<sup>1-3</sup> Upon enzymatic cleavage, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify enzyme activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

### References

1. Xin, X.-Q., Gunsekera, B., and Mason, R.W. The specificity and elastolytic activities of bovine cathepsins S and H. *Arch. Biochem. Biophys.* **299**(2), 334-339 (1992).
2. Darby, N.J., Lackey, D.B., and Smyth, D.G. Purification of a cysteine endopeptidase which is secreted with bioactive peptides from the epidermal glands of *Xenopus laevis*. *Eur. J. Biochem.* **195**(1), 65-70 (1991).
3. Cloete, D., Viljoen, C.C., and Schabort, J.C. Deviation from Michaelis-Menten kinetics for esterase E-II from the venom of *Bitis gabonica*, using synthetic arginine amides as substrates. *Toxicon* **21**(6), 857-869 (1983).

#### 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, 09/30/2022

#### 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