

PRODUCT INFORMATION



Boc-LRR-AMC (trifluoroacetate salt)

Item No. 26642

Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-L-leucyl-L-arginyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-argininamide, trifluoroacetate salt

Synonyms: Boc-Leu-Arg-Arg-AMC, Boc-Leu-Arg-Arg-7-amido-4-Methylcoumarin

MF: C₃₃H₅₂N₁₀O₇ • XCF₃COOH

FW: 700.8

Purity: ≥90%

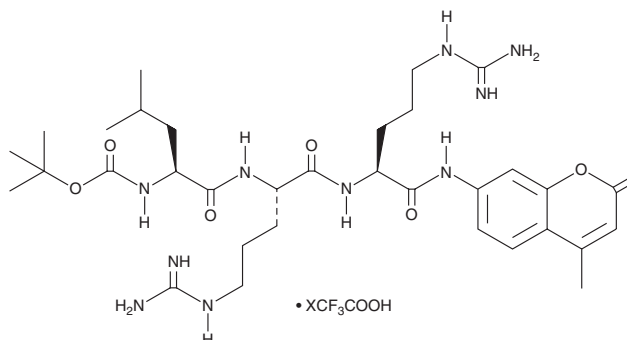
UV/Vis.: λ_{max}: 328 nm

Ex./Em. Max: 340-360/440-460 nm, respectively

Supplied as: A crystalline 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-LRR-AMC (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Boc-LRR-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Boc-LRR-AMC (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Boc-LRR-AMC (trifluoroacetate salt) in these solvents is approximately 10, 12, and 14 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Boc-LRR-AMC (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Boc-LRR-AMC (trifluoroacetate salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Boc-LRR-AMC is a fluorogenic substrate for the trypsin-like activity of the 26S proteasome or 20S proteolytic core.^{1,2} Upon enzymatic cleavage by the 26S proteasome or 20S proteolytic core, amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify 26S proteasome or 20S proteolytic core trypsin-like activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

References

1. Prudnikov, I.M. and Smirnov, A.N. Short peptide tools for monitoring caspase and proteasome activities in embryonal and adult rat brain lysates: An approach for the differential identification of proteases. *J. Biochem.* **151(3)**, 299-316 (2012).
2. Aki, M., Shimbara, N., Takashina, M., *et al.* Interferon-γ induces different subunit organizations and functional diversity of proteasomes. *J. Biochem.* **115(2)**, 257-269 (1994).

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, 04/11/2023

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