

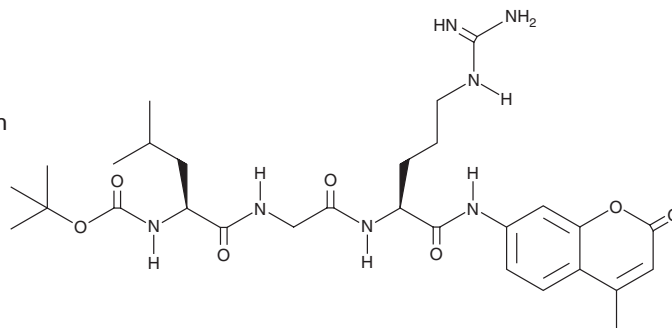
PRODUCT INFORMATION



Boc-Leu-Gly-Arg-AMC

Item No. 45716

CAS Registry No.: 65147-09-3
Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-L-leucylglycyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-argininamide
Synonym: Boc-Leu-Gly-Arg-7-amino-4-methylcoumarin
MF: C₂₉H₄₃N₇O₇
FW: 601.7
Purity: ≥90%
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-Leu-Gly-Arg-AMC is supplied as a solid. A stock solution may be made by dissolving the Boc-Leu-Gly-Arg-AMC in the solvent of choice, which should be purged with an inert gas. Boc-Leu-Gly-Arg-AMC is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Boc-Leu-Gly-Arg-AMC is a fluorogenic substrate for complement convertases, human airway trypsin-like protease (HAT), and urokinase-type plasminogen activator (uPA).¹⁻³ Upon enzymatic cleavage by these proteases, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify complement convertase, HAT, or uPA activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

References

1. Kölln, J., Spillner, E., Andrä, J., *et al.* Complement inactivation by recombinant human C3 derivatives. *J. Immunol.* **173**(9), 5540-5545 (2004).
2. Böttcher-Friebertshäuser, E., Freuer, C., Sielaff, F., *et al.* Cleavage of influenza virus hemagglutinin by airway proteases TMPRSS2 and HAT differs in subcellular localization and susceptibility to protease inhibitors. *J. Virol.* **84**(11), 5605-5614 (2010).
3. Lee, S.-L., Dickson, R.B., and Lin, C.-Y. Activation of hepatocyte growth factor and urokinase/plasminogen activator by matriptase, an epithelial membrane serine protease. *J. Biol. Chem.* **275**(47), 36720-36725 (2000).

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, 07/20/2026

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