

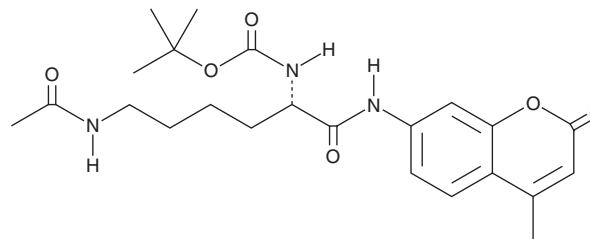
PRODUCT INFORMATION



Boc-Lys(Ac)-AMC

Item No. 29682

CAS Registry No.: 233691-67-3
Formal Name: N-[(1S)-5-(acetylamino)-1-[[[4-methyl-2-oxo-2H-1-benzopyran-7-yl]amino]carbonyl]pentyl]-carbamic acid, 1,1-dimethylethyl ester
MF: C₂₃H₃₁N₃O₆
FW: 445.5
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 229, 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-Lys(Ac)-AMC is supplied as a crystalline solid. A stock solution may be made by dissolving the Boc-Lys(Ac)-AMC in the solvent of choice, which should be purged with an inert gas. Boc-Lys(Ac)-AMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Boc-Lys(Ac)-AMC in these solvents is approximately 3, 25, and 30 mg/ml, respectively.

Description

Boc-Lys(Ac)-AMC is a fluorescent substrate for histone deacetylases (HDACs).¹ Following deacetylation by an HDAC, Boc-Lys-AMC is cleaved by trypsin and 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify HDAC activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Wegener, D., Wirsching, F., Riester, D., *et al.* A fluorogenic histone deacetylase assay well suited for high-throughput activity screening. *Chem. Biol.* **10**(1), 61-68 (2003).

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, 12/09/2022

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