

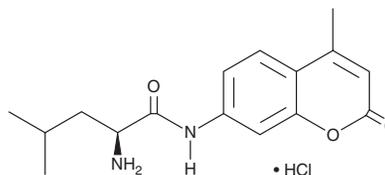
PRODUCT INFORMATION



L-Leu-AMC (hydrochloride)

Item No. 34267

CAS Registry No.:	62480-44-8
Formal Name:	(2S)-2-amino-4-methyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-pentanamide, monohydrochloride
Synonyms:	Leu-MCA, L-Leucine-7-amido-4-methylcoumarin, 7-L-Leucyl-4-methylcoumarinylamide
MF:	$C_{16}H_{20}N_2O_3 \cdot HCl$
FW:	342.8
Purity:	≥98%
Ex./Em. Max:	340-360/440-460 nm
UV/Vis.:	λ_{max} : 214, 296, 326 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

L-Leu-AMC is supplied as a solid. A stock solution may be made by dissolving the L-Leu-AMC in the solvent of choice, which should be purged with an inert gas. L-Leu-AMC is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of L-Leu-AMC in these solvents is approximately 25 mg/ml.

L-Leu-AMC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, L-Leu-AMC should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. L-Leu-AMC has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

L-Leu-AMC is a fluorogenic substrate for leucine aminopeptidase.¹ Upon enzymatic cleavage by leucine aminopeptidase, AMC is released and its fluorescence can be used to quantify leucine aminopeptidase activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Izquierdo, M., Lin, D., O'Neill, S., *et al.* Development of a high-throughput screening assay to identify inhibitors of the major M17-leucyl aminopeptidase from *Trypanosoma cruzi* using rapidfire mass spectrometry. *SLAS Discov.* **25(9)**, 1064-1071 (2020).

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, 12/08/2022

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