

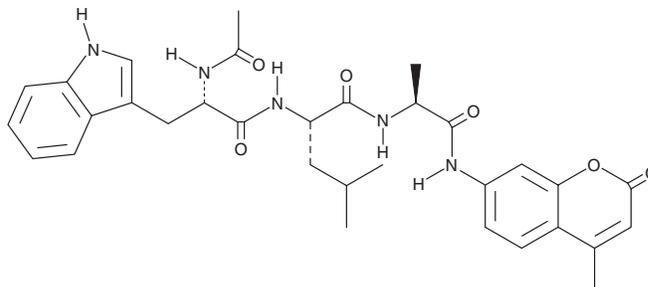
PRODUCT INFORMATION



Ac-WLA-AMC

Item No. 26639

CAS Registry No.: 1104011-59-7
Formal Name: N-acetyl-L-tryptophyl-L-leucyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-alaninamide
Synonyms: Acetyl-Trp-Leu-Ala-7-amino-4-Methylcoumarin, Ac-Trp-Leu-Ala-AMC
MF: C₃₂H₃₇N₅O₆
FW: 587.7
Purity: ≥95%
Ex./Em. Max: 351/430 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

Ac-WLA-AMC is supplied as a solid. A stock solution may be made by dissolving the Ac-WLA-AMC in the solvent of choice, which should be purged with an inert gas. Ac-WLA-AMC is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Ac-WLA-AMC is a fluorogenic substrate for the β 5c subunit of the 20S proteasome.¹ Upon cleavage, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify the activity of the β 5c subunit of the 20S proteasome. AMC displays excitation/emission maxima of 351/430 nm, respectively.

Reference

1. Blackburn, C., Gigstad, K.M., Hales, P., *et al.* Characterization of a new series of non-covalent proteasome inhibitors with exquisite potency and selectivity for the 20S β 5-subunit. *Biochem. J.* **430(3)**, 461-476 (2010).

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, 11/29/2022

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