

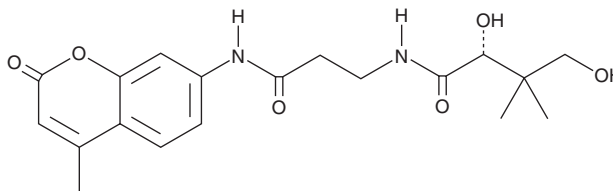
PRODUCT INFORMATION



Pantothenate-AMC

Item No. 9004463

CAS Registry No.: 1224691-12-6
Formal Name: (2R)-2,4-dihydroxy-3,3-dimethyl-N-[3-[(4-methyl-2-oxo-2H-1-benzopyran-7-yl)amino]-3-oxopropyl]-butanamide
Synonym: Pantothenate-7-amino-4-methylcoumarin
MF: C₁₉H₂₄N₂O₆
FW: 376.4
Purity: ≥95%
Ex./Em. Max: 340-360/440-460nm
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

Pantothenate-AMC is supplied as a solid. A stock solution may be made by dissolving the pantothenate-AMC in the solvent of choice, which should be purged with an inert gas. Pantothenate-AMC is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile. It is also slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Pantothenate-AMC is a fluorogenic substrate for vanin-1, an enzyme involved in the recycling of L-pantothenic acid.¹ Upon enzymatic cleavage by vanin-1, 7-amino-7-methylcoumarin (AMC) is released and its fluorescence can be used to quantify vanin-1 activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Ruan, B.H., Cole, D.C., Wu, P., *et al.* A fluorescent assay suitable for inhibitor screening and vanin tissue quantification. *Anal. Biochem.* **399(2)**, 284-292 (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, 07/24/2024

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