

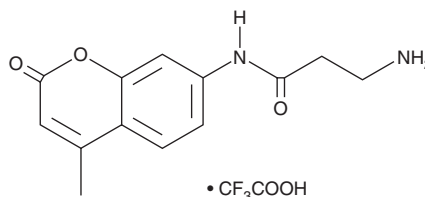
PRODUCT INFORMATION



H-β-Ala-AMC (trifluoroacetate salt)

Item No. 42913

CAS Registry No.: 201847-54-3
Formal Name: 3-amino-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-propanamide, 2,2,2-trifluoroacetate salt
Synonyms: H-β-Alanine-AMC, 7-N-β-Alanylamino-4-methylcoumarin
MF: C₁₃H₁₄N₂O₃ • CF₃COOH
FW: 360.3
Purity: ≥98%
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

H-β-Ala-AMC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the H-β-Ala-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. H-β-Ala-AMC (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Description

H-β-Ala-AMC is a fluorogenic substrate for *P. aeruginosa* β-alanyl aminopeptidase.¹ Upon enzymatic cleavage by β-alanyl aminopeptidase, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify β-alanyl aminopeptidase activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively. H-β-Ala-AMC has also been used in the synthesis of acetylspermidine-AMC, a fluorogenic substrate for histone deacetylase 10 (HDAC10).²

References

1. Váradi, L., Hibbs, D.E., Orega, S., *et al.* β-Alanyl aminopeptidase-activated fluorogenic probes for the rapid identification of *Pseudomonas aeruginosa* in clinical samples. *RSC Advances* **6**(64), (2016).
2. Herp, D., Ridinger, J., Robaa, D., *et al.* First fluorescent acetylspermidine deacetylation assay for HDAC10 identifies selective inhibitors with cellular target engagement. *Chembiochem* **23**(14), e202200180 (2022).

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, 03/11/2025

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