

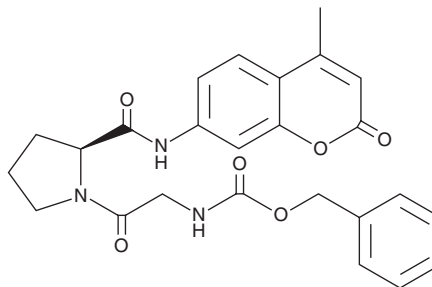
PRODUCT INFORMATION



Cbz-Gly-Pro-AMC

Item No. 42320

CAS Registry No.: 68542-93-8
Formal Name: N-[(phenylmethoxy)carbonyl]glycyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-prolinamide
Synonym: Z-Gly-Pro-MeCouNH
MF: C₂₅H₂₅N₃O₆
FW: 463.5
Purity: ≥98%
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

Cbz-Gly-Pro-AMC is supplied as a solid. A stock solution may be made by dissolving the Cbz-Gly-Pro-AMC in the solvent of choice, which should be purged with an inert gas. Cbz-Gly-Pro-AMC is soluble (≥10 mg/ml) in DMSO and ethanol.

Description

Cbz-Gly-Pro-AMC is a fluorogenic substrate for prolyl endopeptidase.¹ Upon enzymatic cleavage by prolyl endopeptidase, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify prolyl endopeptidase activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Yoshimoto, T., Ogita, K., Walter, R., *et al.* Post-proline cleaving enzyme. Synthesis of a new fluorogenic substrate and distribution of the endopeptidase in rat tissues and body fluids of man. *Biochim. Biophys. Acta* **569(2)**, 184-192 (1979).

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, 12/10/2024

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