

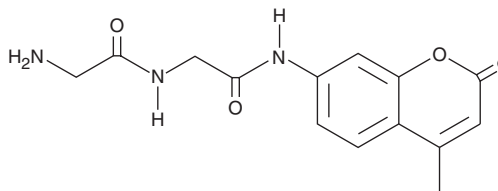
PRODUCT INFORMATION



Gly-Gly-AMC

Item No. 85140

CAS Registry No.: 208645-74-3
Formal Name: glycyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-glycinamide
MF: C₁₄H₁₅N₃O₄
FW: 289.3
Purity: ≥98%
UV/Vis.: λ_{max}: 327 nm
Supplied as: A crystalline 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

Gly-Gly-AMC is supplied as a crystalline solid. A stock solution may be made by dissolving the Gly-Gly-AMC in the solvent of choice, which should be purged with an inert gas. Gly-Gly-AMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Gly-Gly-AMC in ethanol and DMF is approximately 1 mg/ml and approximately 20 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Gly-Gly-AMC can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Gly-Gly-AMC in PBS, pH 7.2, is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Gly-Gly-AMC is a fluorogenic peptide substrate. It has been used to assess the activity of bacterial proteases from *P. aeruginosa* and *S. aureus*.¹ Activity can be quantified by fluorescent detection of free AMC (also known as 7-amino-4-methylcoumarin), which is excited at 340-360 nm and emits at 440-460 nm.

Reference

1. Wildeboer, D., Jeganathan, F., Price, R.G., *et al.* Characterization of bacterial proteases with a panel of fluorescent peptide substrates. *Anal. Biochem.* **384**(2), 321-328 (2009).

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, 10/31/2022

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