

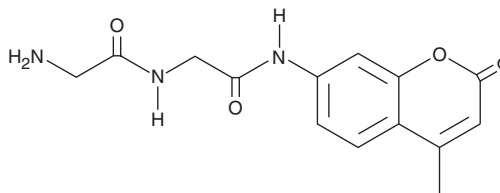
# PRODUCT INFORMATION



## 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:** C<sub>14</sub>H<sub>15</sub>N<sub>3</sub>O<sub>4</sub>  
**FW:** 289.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 327 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 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. Gly-Gly-AMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. 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*.<sup>1</sup> 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, 08/23/2017

#### 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