

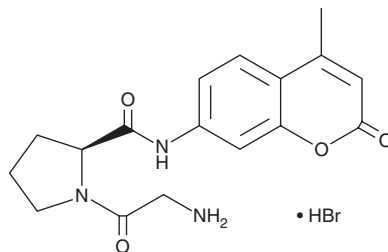
PRODUCT INFORMATION



GP-AMC (hydrobromide)

Item No. 34458

CAS Registry No.: 115035-46-6
Formal Name: glycyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-prolinamide, monohydrobromide
Synonyms: Gly-Pro-AMC, Gly-Pro-7-amido-4-methylcoumarin
MF: $C_{17}H_{19}N_3O_4 \cdot HBr$
FW: 410.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 328 nm
Ex./Em. Max 340-360 nm/440-460 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GP-AMC (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the GP-AMC (hydrobromide) in the solvent of choice, which should be purged with an inert gas. GP-AMC (hydrobromide) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GP-AMC (hydrobromide) in these solvents is approximately 1 and 10 mg/ml, respectively.

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 GP-AMC (hydrobromide) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of GP-AMC (hydrobromide) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GP-AMC is a fluorogenic substrate for dipeptidyl peptidase 4 (DPP-4).^{1,2} Upon enzymatic cleavage by DPP-4, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify DPP-4 activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

References

1. Matheeussen, V., Lambeir, A.-M., Jungraithmayr, W., *et al.* Method comparison of dipeptidyl peptidase IV activity assays and their application in biological samples containing reversible inhibitors. *Clin. Chim. Acta* **413**(3-4), 456-462 (2012).
2. Lammi, C., Bollati, C., Ferruzza, S., *et al.* Soybean- and lupin-derived peptides inhibit DPP-IV activity on in situ human intestinal Caco-2 cells and ex vivo human serum. *Nutrients* **10**(8), 1082 (2018).

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, 12/13/2022

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