

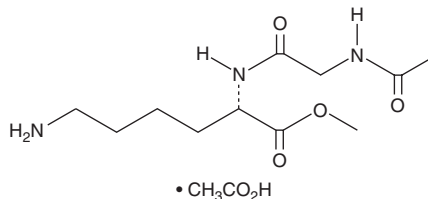
PRODUCT INFORMATION



Ac-Gly-Lys-OMe (acetate)

Item No. 30568

CAS Registry No.: 14752-92-2
Formal Name: N-acetylglycyl-L-lysine, methyl ester, monoacetate
Synonym: α -N-Acetylglycyl-L-Lysine
MF: $C_{11}H_{21}N_3O_4 \cdot C_2H_4O_2$
FW: 319.4
Purity: $\geq 95\%$
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

Ac-Gly-Lys-OMe (acetate) is supplied as a crystalline solid. Aqueous solutions of Ac-Gly-Lys-OMe (acetate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Ac-Gly-Lys-OMe (acetate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-Gly-Lys-OMe is a synthetic peptide urokinase substrate.^{1,2} It has been used to characterize the effects of small molecule inhibitors and metal ions on urokinase activity in cell-free assays.

References

1. Walton, P.L. The hydrolysis of α -N-acetylglycyl-L-lysine methyl ester by urokinase. *Biochim. Biophys. Acta* **132(1)**, 104-114 (1967).
2. Sumi, H., Takada, Y., and Takada, A. The effects of metal ions on esterase activities of urokinase. *Thromb. Haemost.* **39(1)**, 46-52 (1978).

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, 11/03/2022

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