

PRODUCT INFORMATION



GGACK (hydrochloride)

Item No. 34618

Formal Name: L- α -glutamyl-N-[(1S)-4-[(aminoiminomethyl)amino]-1-(2-chloroacetyl)butyl]-glycinamide, monohydrochloride

Synonyms: Glu-Gly-Arg-Chloromethyl Ketone, Glu-Gly-Arg-CH₂Cl

MF: C₁₄H₂₅ClN₆O₅ • HCl

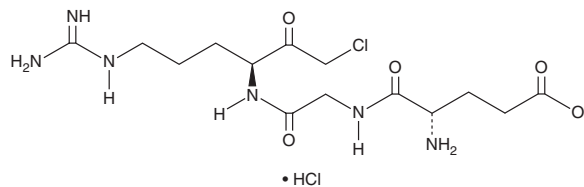
FW: 429.3

Purity: \geq 95%

Supplied as: A 10 mg/ml solution in water

Storage: -20°C

Stability: \geq 4 years



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

Description

GGACK is an inhibitor of urokinase-type plasminogen activator (uPA) and factor Xa (K_i s = 1.8 and 8.7 μ M, respectively).^{1,2}

References

1. Kettner, C. and Shaw, E. The susceptibility of urokinase to affinity labeling by peptides of arginine chloromethyl ketone. *Biochim. Biophys. Acta* **569**(1), 31-40 (1979).
2. Kettner, C. and Shaw, E. The selective affinity labeling of factor Xa by peptides of arginine chloromethyl ketone. *Thromb. Res.* **22**(5-6), 645-652 (1981).

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, 03/25/2026

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