

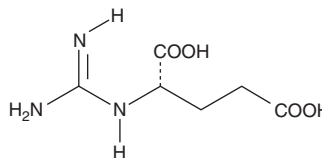
PRODUCT INFORMATION



α -Guanidinoglutaric Acid

Item No. 80340

CAS Registry No.: 73477-53-9
Formal Name: N-(aminoiminomethyl)-L-glutamic acid
Synonyms: GGA
MF: C₆H₁₁N₃O₄
FW: 189.2
Purity: ≥98%
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

α -Guanidinoglutaric acid is supplied as a crystalline solid. Aqueous solutions of α -guanidinoglutaric acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of α -guanidinoglutaric acid in PBS (pH 7.2) is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day. Due to the highly charged characteristics of GGA at neutral pH, the compound will not be cell membrane permeable.

Description

α -Guanidinoglutaric Acid (GGA) is a non-N^G-substituted inhibitor of nitric oxide synthase (NOS).¹ GGA is a linear mixed-type inhibitor of nNOS having a K_i of 2.69 μ M.¹ It is normally present in small quantities in the cerebral cortex of cats with increased levels following seizure.²

References

1. Yokoi, I., Kabuto, H., Habu, H., *et al.* α -Guanidinoglutaric acid, an endogenous convulsant, as a novel nitric oxide synthase inhibitor. *J. Neurochem.* **63(4)**, 1565-1567 (1994).
2. Mori, A., Akagi, M., Katayama, Y., *et al.* α -Guanidinoglutaric acid in cobalt-induced epileptogenic cerebral cortex of cats. *J. Neurochem.* **35(3)**, 603-605 (1980).

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, 12/07/2022

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