

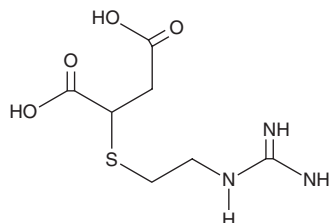
PRODUCT INFORMATION



2-Guanidinoethylmercaptosuccinic Acid

Item No. 19354

CAS Registry No.: 77482-44-1
Formal Name: 2-[[2-[(aminoiminomethyl)amino]ethyl]thio]-butanedioic acid
Synonym: GEMSA
MF: C₇H₁₃N₃O₄S
FW: 235.3
Purity: ≥98%
Supplied as: A 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

2-Guanidinoethylmercaptosuccinic acid (GEMSA) is supplied as a solid. A stock solution may be made by dissolving the GEMSA in water. We do not recommend storing the aqueous solution for more than one day.

Description

Enkephalin convertase, also known as carboxypeptidase E, removes C-terminal residues during the processing of propeptides, including enkephalin and proinsulin. GEMSA is a potent inhibitor of enkephalin convertase ($K_i = 8.8$ nM) that is functional *in vivo*.^{1,2} It is a peptidomimetic inhibitor, occupying the specificity pocket of the enzyme.³ GEMSA can also inhibit other carboxypeptidases, including thrombin-activatable fibrinolysis inhibitor.^{4,5}

References

1. Przewlocka, B., Dziedzicka, M., Silberring, J., *et al.* Analgesic and convulsant effects of guanidinoethylmercaptosuccinic acid (GEMSA)-a potent enkephalin convertase inhibitor. *Neuropeptides* **8(4)**, 359-365 (1986).
2. Bommer, M., Nikolarakis, K., Noble, E.P., *et al.* *In vivo* modulation of rat hypothalamic opioid peptide content by intracerebroventricular injection of guanidinoethylmercaptosuccinic acid (GEMSA): Possible physiological role of enkephalin convertase. *Brain Res.* **492(1-2)**, 305-313 (1989).
3. Aloy, P., Companys, V., Vendrell, J., *et al.* The crystal structure of the inhibitor-complexed carboxypeptidase D domain II and the modeling of regulatory carboxypeptidases. *J. Biol. Chem.* **276(19)**, 16177-16184 (2001).
4. Bajzar, L., Manuel, R., and Nesheim, M.E. Purification and characterization of TAFI, a thrombin-activable fibrinolysis inhibitor. *J. Biol. Chem.* **270(24)**, 14477-14484 (1995).
5. Ore, F.D., Ajandouz, E.H., Giardina, T., *et al.* The membrane-bound basic carboxypeptidase from hog intestinal mucosa. *Biochim. Biophys. Acta.* **1421(2)**, 234-248 (2016).

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, 10/06/2022

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