

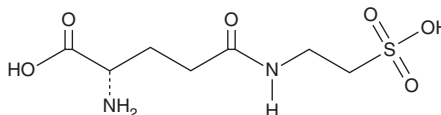
PRODUCT INFORMATION



Glutaurine

Item No. 39586

CAS Registry No.: 56488-60-9
Formal Name: N-(2-sulfoethyl)-L-glutamine
Synonyms: γ -Glutamyl Taurine, γ -L-Glutamyl Taurine
MF: C₇H₁₄N₂O₆S
FW: 254.3
Purity: $\geq 90\%$
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

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

Description

Glutaurine is a dipeptide composed of L-glutamic acid (Item No. 30377) and taurine (Item No. 27031) that has been found in rat brain.¹ It is formed from glutathione and taurine by γ -glutamyl transpeptidase. Glutaurine enhances potassium-induced release of L-glutamate and inhibits kainate-induced formation of cGMP in isolated mouse cerebral cortical slices when used at concentrations of 0.1 and 1 mM, respectively.² It decreases plasma levels of triiodothyronine (T3) and increases plasma thyroid-stimulating hormone (TSH) levels in rats when administered at a dose of 600 μ g/kg.³ Glutaurine (50 μ g/animal, p.o.) prevents electroconvulsive shock-induced amnesia in rats.⁴

References

1. Török, K., Varga, V., Somogyi, J., *et al.* Formation of γ -glutamyl-*taurine* in the rat brain. *Neurosci. Lett.* **27(2)**, 145-149 (1981).
2. Varga, V., Janáky, R., Marnela, K.-M., *et al.* Interactions of γ -L-glutamyltaurine with excitatory aminoacidergic neurotransmission. *Neurochem. Res.* **19(3)**, 243-248 (1994).
3. Baskin, S.I., Bartuska, D., Thampi, N., *et al.* The effect of glutaurine on thyroid hormones in the rat. *Neuropeptides* **9(1)**, 45-50 (1987).
4. Balazs, M. and Telegdy, G. Effects of glutaurine treatment on electroshock-induced amnesia. Antiamnesic action of glutaurine. *Neuropeptides* **12(2)**, 55-58 (1988).

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, 08/18/2023

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