

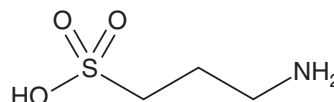
PRODUCT INFORMATION



Tramiprosate

Item No. 20701

CAS Registry No.: 3687-18-1
Formal Name: 3-amino-1-propanesulfonic acid
Synonyms: Homotaurine, NC 531, NC 758, NSC 77071, 3-Sulfopropylamine
MF: C₃H₉NO₃S
FW: 139.2
Purity: ≥95%
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Tramiprosate is supplied as a crystalline solid. Aqueous solutions of tramiprosate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of tramiprosate in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Tramiprosate is a sulfated glycosaminoglycan mimetic that was reported to bind soluble A β and inhibit the interaction of A β with endogenous glycosaminoglycans, thereby preventing β -sheet formation.¹ Conversely, tramiprosate was also reported to promote tau polymerization into fibrillar aggregates.² It is also an analog of GABA that can act as a partial agonist of GABA_A receptors, as well as an antagonist at GABA_B receptors.³

References

1. Gervais, F., Chalifour, R., Garceau, D., *et al.* Glycosaminoglycan mimetics: A therapeutic approach to cerebral amyloid angiopathy. *Amyloid* **8(Suppl. 1)**, 28-35 (2001).
2. Santa-Maria, I., Hernandez, F., Del Rio, J., *et al.* Tramiprosate, a drug of potential interest for the treatment of Alzheimer's disease, promotes an abnormal aggregation of tau. *Molecular Neurodegeneration* **2**, 17-29 (2007).
3. Giotti, A., Luzzi, S.S. and Zilletti, L. Homotaurine: A GABA_B antagonist in guinea-pig ileum. *Br. J. Pharmac.* **79**, 855-862 (1983).

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, 11/08/2022

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