

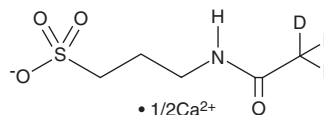
# PRODUCT INFORMATION



## Acamprosate-d<sub>3</sub> (calcium salt)

Item No. 26780

**CAS Registry No.:** 1225580-94-8  
**Formal Name:** 3-(acetyl-2,2,2-d<sub>3</sub>-amino)-1-propanesulfonic acid, hemicalcium salt (2:1)  
**Synonyms:** N-acetyl Homotaurinate-d<sub>3</sub>, N-acetyl Homotaurine-d<sub>3</sub>  
**MF:** C<sub>5</sub>H<sub>7</sub>D<sub>3</sub>NO<sub>4</sub>S • 1/2Ca  
**FW:** 203.3  
**Chemical Purity:** ≥98% (Acamprosate)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>3</sub>); ≤1% d<sub>0</sub>  
**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

Acamprosate-d<sub>3</sub> (calcium salt) is intended for use as an internal standard for the quantification of acamprosate (Item No. 23899) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

### Description

Acamprosate is an acetylated derivative of the GABA analog homotaurine (tramiprosate; Item No. 20701).<sup>1</sup> Despite its structural similarity to GABA, acamprosate does not act at GABA<sub>A</sub> receptors but does decrease paired-pulse inhibition of GABA<sub>A</sub> inhibitory post-synaptic currents (IPSCs) at short inter-stimulus intervals when used at a concentration of 300 μM, indicating that it may inhibit GABA<sub>B</sub> autoreceptor-mediated inhibition of GABA release.<sup>2-4</sup> It is an NMDA receptor modulator with antagonist or agonist effects depending on brain region, receptor subunit composition, and other factors.<sup>4</sup> Acamprosate (0.26 and 0.52 mmol/kg per day, i.p.) reduces voluntary intake of ethanol in rats, an effect that can be blocked by the GABA antagonist bicuculline (Item No. 11727). Formulations containing acamprosate have been used for the maintenance of alcohol abstinence.

### References

1. Boismare, F., Daoust, M., Moore, N., *et al. Pharmacol. Biochem. Behav.* **21(5)**, 787-789 (1984).
2. Reilly, M.T., Lobo, I.A., McCracken, L.M., *et al. Alcohol. Clin. Exp. Res.* **32(2)**, 188-196 (2008).
3. Berton, F., Francesconi, W.G., Madamba, S.G., *et al. Alcohol Clin. Exp. Res.* **22(1)**, 183-191 (1998).
4. Tomek, S.E., Lacrosse, A.L., Nemirovsky, N.E., *et al. Pharmaceuticals (Basel)* **6(2)**, 251-268 (2013).

#### 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, 12/21/2022

#### 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