

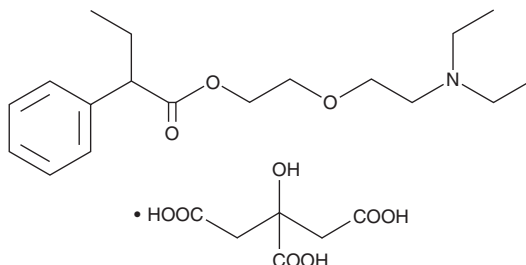
PRODUCT INFORMATION



Butamirate (citrate)

Item No. 42561

CAS Registry No.: 18109-81-4
Formal Name: α -ethyl-benzeneacetic acid, 2-[2-(diethylamino)ethoxy] ethyl ester, 2-hydroxy-1,2,3-propanetricarboxylate
MF: $C_{18}H_{29}NO_3 \cdot C_6H_8O_7$
FW: 499.6
Purity: $\geq 98\%$
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

Butamirate (citrate) is supplied as a solid. A stock solution may be made by dissolving the butamirate (citrate) in the solvent of choice, which should be purged with an inert gas. Butamirate (citrate) is sparingly soluble (1-10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of butamirate (citrate) can be prepared by directly dissolving the solid in aqueous buffers. Butamirate (citrate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Butamirate is an antitussive agent.¹ It reduces proliferation, STAT3 activity, migration, and invasion in U87 glioma tumorspheres. Butamirate (2 mg/kg per day) decreases tumor volume in a U87MG glioblastoma mouse xenograft model. Formulations containing butamirate have been used as cough suppressants.

Reference

1. Lee, S.-J., Yeom, S.-Y., Lee, J.-Y., *et al.* Application of the antitussive agents oxelaidin and butamirate as anti-glioma agents. *Sci. Rep.* **11**(1), 10145 (2021).

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, 01/22/2025

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