

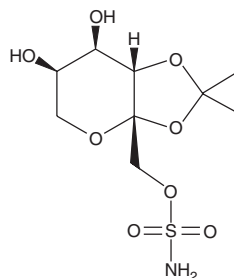
PRODUCT INFORMATION



5-Desisopropylidene Topiramate

Item No. 35071

CAS Registry No.: 106881-41-8
Formal Name: 2,3-O-(1-methylethylidene)-β-D-fructopyranose 1-sulfamate
Synonym: 4,5-Desisopropylidene TPM
MF: C₉H₁₇NO₈S
FW: 299.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

4,5-Desisopropylidene topiramate is supplied as a solid. A stock solution may be made by dissolving the 4,5-desisopropylidene topiramate in the solvent of choice, which should be purged with an inert gas. 4,5-Desisopropylidene topiramate is slightly soluble in methanol (sonicated). It is also slightly soluble in water (heated). We do not recommend storing the aqueous solution for more than one day.

Description

4,5-Desisopropylidene topiramate is an inactive metabolite of the anticonvulsant topiramate.^{1,2}

References

1. Maryanoff, B.E., Nortey, S.O., Gardocki, J.F., *et al.* Anticonvulsant O-alkyl sulfamates. 2,3:4,5-Bis-O-(1-methylethylidene)-β-D-fructopyranose sulfamate and related compounds. *J. Med. Chem.* **30(5)**, 880-887 (1987).
2. Milosheska, D. A novel LC-MS/MS method for the simultaneous quantification of topiramate and its main metabolites in human plasma. *J. Pharm. Biomed. Anal.* **138**, 180-188 (2017).

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, 03/31/2022

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