

PRODUCT INFORMATION



N-desethyl Brinzolamide (oxalate)

Item No. 34275

CAS Registry No.: 2775292-24-3

Formal Name: (4R)-4-amino-3,4-dihydro-2-(3-methoxypropyl)-2H-thieno[3,2-e]-1,2-thiazine-6-sulfonamide, 1,1-dioxide, monoethanedioate

Synonym: AL 8520

MF: $C_{10}H_{17}N_3O_5S_3 \cdot C_2H_2O_4$

FW: 445.5

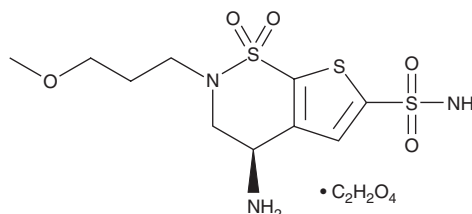
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 255 nm

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

N-desethyl Brinzolamide (oxalate) is supplied as a solid. A stock solution may be made by dissolving the N-desethyl brinzolamide (oxalate) in the solvent of choice, which should be purged with an inert gas. N-desethyl Brinzolamide (oxalate) is soluble in DMSO.

Description

N-desethyl Brinzolamide is an active metabolite of the carbonic anhydrase (CA) inhibitor brinzolamide (Item No. 11660).¹ It inhibits CAII and CAIV activities (IC_{50} s = 1.28 and 128 nM, respectively).

Reference

1. Kim, C.-Y., Whittington, D.A., Chang, J.S., *et al.* Structural aspects of isozyme selectivity in the binding of inhibitors to carbonic anhydrases II and IV. *J. Med. Chem.* **45**(4), 888-893 (2002).

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, 09/14/2023

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