

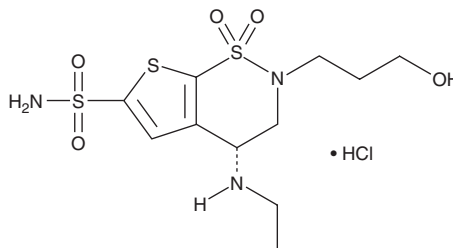
# PRODUCT INFORMATION



## O-desmethyl Brinzolamide (hydrochloride)

Item No. 34276

**CAS Registry No.:** 2967475-56-3  
**Formal Name:** (4R)-(ethylamino)-3,4-dihydro-2-(3-hydroxypropyl)-2H-thieno[3,2-e]-1,2-thiazine-6-sulfonamide, 1,1-dioxide, monohydrochloride  
**MF:**  $C_{11}H_{19}N_3O_5S_3 \cdot HCl$   
**FW:** 405.9  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 256 nm  
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

O-desmethyl Brinzolamide is an active metabolite of the carbonic anhydrase (CA) inhibitor brinzolamide (Item No. 11660).<sup>1,2</sup> It inhibits CAII and CAIV ( $IC_{50}$ s = 0.136 and 165 nM, respectively).<sup>1</sup>

### References

1. Huang, Q., Rui, E.Y., Cobbs, M., *et al.* Design, synthesis, and evaluation of NO-donor containing carbonic anhydrase inhibitors to lower intraocular pressure. *J. Med. Chem.* **58**(6), 2821-2833 (2015).
2. Lo Faro, A.F., Tini, A., Gottardi, M., *et al.* Development and validation of a fast ultra-high-performance liquid chromatography tandem mass spectrometry method for determining carbonic anhydrase inhibitors and their metabolites in urine and hair. *Drug Test Anal.* **13**(8), 1552-1560 (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, 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