

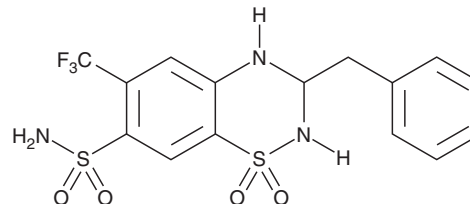
# PRODUCT INFORMATION



## Bendroflumethiazide

Item No. 21311

**CAS Registry No.:** 73-48-3  
**Formal Name:** 3,4-dihydro-3-(phenylmethyl)-6-(trifluoromethyl)-2H-1,2,4-benzothiadiazine-7-sulfonamide-1,1-dioxide  
**Synonym:** Bendrofluazide  
**MF:** C<sub>15</sub>H<sub>14</sub>F<sub>3</sub>N<sub>3</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 421.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 271, 324 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Bendroflumethiazide (Item No. 21311) is an analytical reference standard categorized as a diuretic.<sup>1,2</sup> Diuretics, including bendroflumethiazide, have been abused as performance-enhancing drugs and masking agents in sports doping.<sup>3</sup> This product is intended for research and forensic applications.

### References

1. Majid, D.S.A. and Navar, L.G. Blockade of distal nephron sodium transport attenuates pressure natriuresis in dogs. *Hypertension* **23(6 Pt 2)**, 1040-1045 (1994).
2. Vongpatanasin, W. Hydrochlorothiazide (HCTZ) is not the most useful nor versatile thiazide diuretic. *Curr. Opin. Cardiol.* **30(4)**, 361-365 (2015).
3. Cadwallader, A.B., de la Torre, X., Tieri, A., et al. The abuse of diuretics as performance-enhancing drugs and masking agents in sport doping: Pharmacology, toxicology and analysis. *Br. J. Pharmacol.* **161(1)**, 1-16 (2010).

#### 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, 07/06/2023

#### 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