

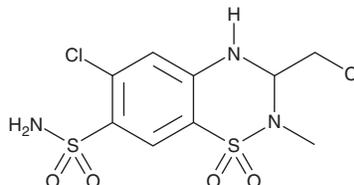
# PRODUCT INFORMATION



## Methyclothiazide

Item No. 23799

**CAS Registry No.:** 135-07-9  
**Formal Name:** 6-chloro-3-(chloromethyl)-3,4-dihydro-2-methyl-2H-1,2,4-benzothiadiazine-7-sulfonamide, 1,1-dioxide  
**Synonym:** NSC 110431  
**MF:** C<sub>9</sub>H<sub>11</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 360.2  
**Purity:** ≥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

Methyclothiazide is supplied as a solid. A stock solution may be made by dissolving the methyclothiazide in the solvent of choice, which should be purged with an inert gas. Methyclothiazide is slightly soluble in organic solvents such as DMSO and methanol.

### Description

Methyclothiazide is a thiazide diuretic.<sup>1</sup> It inhibits norepinephrine- (Item No. 16673) and vasopressin-induced contractions of aortic rings isolated from spontaneously hypertensive rats at a concentration of 100 μM. Methyclothiazide reduces salt-induced increases in blood pressure in salt-sensitive but not salt-resistant rats without affecting body weight or heart rate.<sup>2</sup> Formulations containing methyclothiazide have been used to treat hypertension.

### References

1. Colas, B., Slama, M., Masson, H., et al. Direct vascular actions of methyclothiazide and indapamide in aorta of spontaneously hypertensive rat. *Fundam. Clin. Pharmacol.* **14(4)**, 363-368 (2000).
2. Sasaki, S. and Buñag, R.D. Methyclothiazide attenuates salt-induced hypertension without affecting sympathetic responsiveness in Dahl rats. *J. Cardiovasc. Pharmacol.* **5(3)**, 378-383 (1983).

#### 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, 11/16/2022

#### 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