

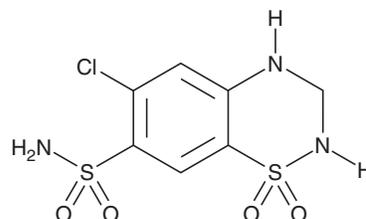
PRODUCT INFORMATION



Hydrochlorothiazide

Item No. 21304

CAS Registry No.: 58-93-5
Formal Name: 6-chloro-3,4-dihydro-2H-1,2,4-benzothiadiazine-7-sulfonamide, 1,1-dioxide
Synonyms: HCTZ, NSC 53477, SU 5879
MF: C₇H₈ClN₃O₄S₂
FW: 297.7
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 270, 315 nm
Supplied as: A crystalline 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

Hydrochlorothiazide (HCTZ) is supplied as a crystalline solid. A stock solution may be made by dissolving the HCTZ in the solvent of choice, which should be purged with an inert gas. HCTZ is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of HCTZ in these solvents is approximately 20 and 30 mg/ml, respectively.

HCTZ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HCTZ should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HCTZ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HCTZ is a thiazide diuretic that inhibits the apical membrane sodium chloride cotransporter in the distal and collecting duct nephron segments.¹⁻³

References

1. Odland, B., and Lönnerholm, G. Renal tubular secretion and effects of chlorothiazide, hydrochlorothiazide and clopamide: A study in the avian kidney. *Acta. Pharmacol. Toxicol (Copenh)* **51(3)**, 187-197 (1982).
2. 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).
3. Vongpatanasin, W. Hydrochlorothiazide (HCTZ) is not the most useful nor versatile thiazide diuretic. *Curr. Opin. Cardiol.* **30(4)**, 361-365 (2015).

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, 01/31/2024

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