

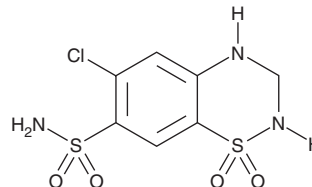
PRODUCT INFORMATION



Hydrochlorothiazide

Item No. 22671

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 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

Hydrochlorothiazide (Item No. 22671) is an analytical reference standard categorized as a thiazide diuretic. Diuretics, including hydrochlorothiazide, have been misused in sports for weight reduction and as masking agents.¹ This product is intended for use in analytical forensic applications. A general research tool is also available as hydrochlorothiazide (Item No. 21304).

Reference

1. Cooper, S.F., Masse, R., and Dugal, R. Comprehensive screening procedure for diuretics in urine by high-performance liquid chromatography. *J. Chromatogr. A.* **489(1)**, 65-88 (1989).

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, 06/16/2023

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