

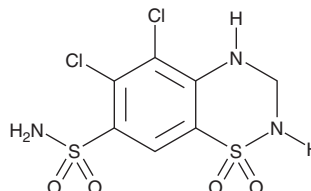
PRODUCT INFORMATION



5-chloro Hydrochlorothiazide

Item No. 17571

CAS Registry No.: 5233-42-1
Formal Name: 5,6-dichloro-3,4-dihydro-2H-1,2,4-benzothiadiazine-7-sulfonamide, 1,1-dioxide
MF: $C_7H_7Cl_2N_3O_4S_2$
FW: 332.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 230, 272 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

5-chloro Hydrochlorothiazide is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-chloro hydrochlorothiazide in the solvent of choice, which should be purged with an inert gas. 5-chloro Hydrochlorothiazide is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of 5-chloro hydrochlorothiazide in these solvents is approximately 20 and 30 mg/ml, respectively.

5-chloro Hydrochlorothiazide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-chloro hydrochlorothiazide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 5-chloro Hydrochlorothiazide 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

5-chloro Hydrochlorothiazide is a derivative of hydrochlorothiazide, which is a diuretic and antihypertensive agent that increases renal excretion of sodium, potassium, chloride, and bicarbonate ions by inhibiting tubular reabsorptive mechanisms.^{1,2}

References

1. Laragh, J.H. The mode of action and use of chlorothiazide and related compounds. *Circulation* **26**, 121-132 (1962).
2. Nath, M.R. and Saira, M. Synthesis, isolation, and characterization of hydrochlorothiazide dimer impurity. *J. Pharm. Sci.* **5(Suppl 3)**, 867-871 (2013).

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, 11/10/2022

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