

PRODUCT INFORMATION



Hydrochlorothiazide Impurity C

Item No. 22384

CAS Registry No.: 402824-96-8
Formal Name: N-[[7-(aminosulfonyl)-6-chloro-2,3-dihydro-1,1-dioxido-4H-1,2,4-benzothiadiazin-4-yl]methyl]-6-chloro-3,4-dihydro-2H-1,2,4-benzothiadiazine-7-sulfonamide, 1,1-dioxide

MF: C₁₅H₁₆Cl₂N₆O₈S₄

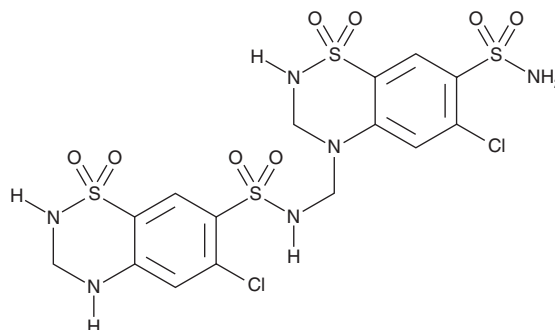
FW: 607.5

Purity: ≥90%

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

Hydrochlorothiazide impurity C is supplied as a solid. A stock solution may be made by dissolving the hydrochlorothiazide impurity C in the solvent of choice, which should be purged with an inert gas. Hydrochlorothiazide impurity C is soluble in the organic solvent DMSO.

Description

Hydrochlorothiazide impurity C is a potential impurity found in commercial hydrochlorothiazide preparations.¹ Hydrochlorothiazide (Item No. 21304) is a thiazide diuretic that inhibits the apical membrane sodium chloride cotransporter in the distal and collecting duct nephron segments.²⁻⁴

References

1. Fang, X., Bibart, R.T., Mayr, S., *et al.* Purification and identification of an impurity in bulk hydrochlorothiazide. *J. Pharm. Sci.* **90(11)**, 1800-1809 (2001).
2. 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).
3. 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).
4. 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, 01/03/2023

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