

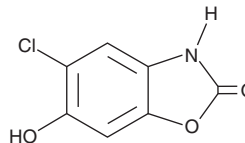
PRODUCT INFORMATION



6-hydroxy Chlorzoxazone

Item No. 10009029

CAS Registry No.: 1750-45-4
Formal Name: 5-chloro-6-hydroxy-2(3H)-benzoxazolone
Synonym: NSC 24955
MF: C₇H₄ClNO₃
FW: 185.6
Purity: ≥95%
UV/Vis.: λ_{max}: 301 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

6-hydroxy Chlorzoxazone is supplied as a solid. A stock solution may be made by dissolving the 6-hydroxy chlorzoxazone in the solvent of choice, which should be purged with an inert gas. 6-hydroxy Chlorzoxazone is slightly soluble in methanol and DMSO.

Description

6-hydroxy Chlorzoxazone is a metabolite of chlorzoxazone (Item No. 18869).¹ It is formed by metabolism of chlorzoxazone by the cytochrome P450 (CYP) isoform CYP2E1 in human hepatic microsomes.

Reference

1. Court, M.H., Von Moltke, L.L., Shader, R.I., *et al.* Biotransformation of chlorzoxazone by hepatic microsomes from humans and ten other mammalian species. *Biopharm. Drug Dispos.* **18(3)**, 213-226 (1997).

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, 04/18/2019

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