

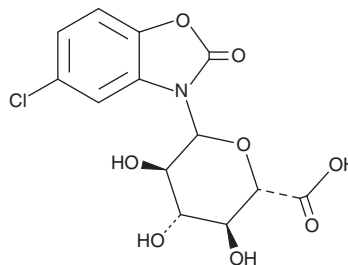
PRODUCT INFORMATION



Chlorzoxazone N-Glucuronide (lithium salt)

Item No. 19879

Formal Name: (2S,3S,4S,5R)-6-(5-chloro-2-oxobenzo[d]oxazol-3(2H)-yl)-3,4,5-trihydroxytetrahydro-2H-pyran-2-carboxylic acid
MF: C₁₃H₁₂ClNO₈
FW: 345.7
Purity: ≥95%
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

Chlorzoxazone N-glucuronide (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the chlorzoxazone N-glucuronide (lithium salt) in the solvent of choice, which should be purged with an inert gas. Chlorzoxazone N-glucuronide (lithium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of chlorzoxazone N-glucuronide (lithium salt) in these solvents is approximately 20, 10, and 16 mg/ml, respectively.

Description

Chlorzoxazone N-glucuronide is a phase II metabolite of the skeletal muscle relaxant chlorzoxazone (Item No. 18869).¹ Chlorzoxazone N-glucuronide is formed *via* glucuronidation of chlorzoxazone by the UDP-glucuronosyltransferase (UGT) isoform UGT1A9.

Reference

1. Quesnot, N., Bucher, S., Gade, C., *et al.* Production of chlorzoxazone glucuronides via cytochrome P4502E1 dependent and independent pathways in human hepatocytes. *Arch. Toxicol.* **92(10)**, 3077-3091 (2018).

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, 11/07/2022

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