

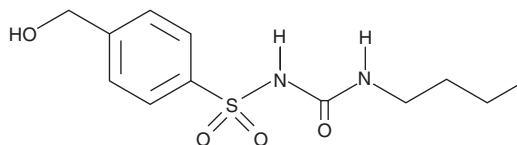
PRODUCT INFORMATION



4-hydroxy Tolbutamide

Item No. 10008520

CAS Registry No.: 5719-85-7
Formal Name: N-[(butylamino)carbonyl]-4-(hydroxymethyl)-benzenesulfonamide
MF: $C_{12}H_{18}N_2O_4S$
FW: 286.4
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 229 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.

Description

4-hydroxy Tolbutamide is a metabolite of the potassium channel antagonist tolbutamide (Item No. 19888).¹ It is formed from tolbutamide by the cytochrome P450 (CYP) isoform CYP2C9 in human liver microsomes and COS-7 cell lysates expressing the human enzyme. 4-hydroxy Tolbutamide has been used as a marker of CYP2C9 activity in rat hepatocytes.²

References

1. Miners, J.O. and Birkett, D.J. Use of tolbutamide as a substrate probe for human hepatic cytochrome P450 2C9. *Methods Enzymol.* **272**, 139-145 (1996).
2. Nguyen, V., Espiritu, M., and Elbarbry, F. Development and validation of a sensitive and specific LC-MS/MS cocktail assay for CYP450 enzymes: Application to study the effect of catechin on rat hepatic CYP activity. *Biomed. Chromatogr.* **34(3)**, e4789 (2019).

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, 01/22/2024

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