

PRODUCT INFORMATION



β-Glucuronidase Inhibitor

Item No. 41228

CAS Registry No.: 484006-66-8

Formal Name: N-[(1,2-dihydro-6,8-dimethyl-2-oxo-3-quinolinyl)methyl]-N'-(4-ethoxyphenyl)-N-(2-hydroxyethyl)-thiourea

MF: C₂₃H₂₇N₃O₃S

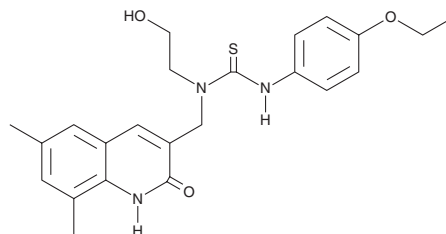
FW: 425.5

Purity: ≥98%

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

β-Glucuronidase inhibitor is supplied as a solid. A stock solution may be made by dissolving the β-glucuronidase inhibitor in the solvent of choice, which should be purged with an inert gas. β-Glucuronidase inhibitor is sparingly soluble (1-10 mg/ml) in DMSO.

Description

β-Glucuronidase inhibitor is an inhibitor of bacterial β-glucuronidase (IC₅₀ = 283 nM for the *E. coli* enzyme).¹ It inhibits wild-type but not Δ360-376 *E. coli* β-glucuronidase, a mutant lacking the bacteria-specific loop, when used at a concentration of 100 μM. *In vivo*, β-glucuronidase inhibitor (50 mg/kg) reduces colonic epithelial damage and the number of bloody diarrheal events induced by irinotecan (CPT-11) in a mouse model of chemotherapy-induced intestinal toxicity.

Reference

1. Wallace, B.D., Wang, H., Lane, K.T., *et al.* Alleviating cancer drug toxicity by inhibiting a bacterial enzyme. *Science* **330**(6005), 831-835 (2010).

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, 07/26/2024

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