

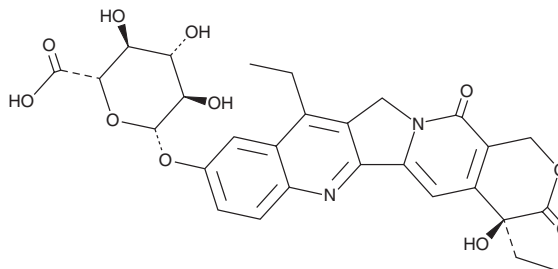
# PRODUCT INFORMATION



## SN-38 Glucuronide

Item No. 35642

**CAS Registry No.:** 121080-63-5  
**Formal Name:** (4S)-4,11-diethyl-3,4,12,14-tetrahydro-4-hydroxy-3,14-dioxo-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinolin-9-yl,  $\beta$ -D-glucopyranosiduronic acid  
**Synonym:** SN-38G  
**MF:** C<sub>28</sub>H<sub>28</sub>N<sub>2</sub>O<sub>11</sub>  
**FW:** 568.5  
**Purity:**  $\geq$ 95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

SN-38 Glucuronide is supplied as a solid. A stock solution may be made by dissolving the SN-38 glucuronide in the solvent of choice, which should be purged with an inert gas. SN-38 Glucuronide is slightly soluble in methanol and DMSO.

### Description

SN-38 Glucuronide is a phase II metabolite of irinotecan (Item No. 14180).<sup>1</sup> It is formed *via* glucuronidation of the irinotecan phase I metabolite SN-38 (Item No. 15632) by the UDP-glucuronosyltransferase (UGT) isoform UGT1A1.

### Reference

1. Hirose, K., Kozu, C., Yamashita, K., *et al.* Correlation between plasma concentration ratios of SN-38 glucuronide and SN-38 and neutropenia induction in patients with colorectal cancer and wild-type UGT1A1 gene. *Oncol. Lett.* **3**(3), 694-698 (2012).

**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, 09/26/2023

### 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