

# PRODUCT INFORMATION

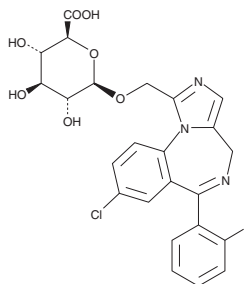


## 1 $\alpha$ -Hydroxymidazolam $\beta$ -D-Glucuronide

Item No. 10386

**CAS Registry No.:** 81256-81-7  
**Formal Name:** [8-chloro-6-(2-fluorophenyl)-4H-imidazo[1,5-a][1,4]benzodiazepin-1-yl]methyl- $\beta$ -D-glucopyranosiduronic acid

**MF:** C<sub>24</sub>H<sub>21</sub>ClFN<sub>3</sub>O<sub>7</sub>  
**FW:** 517.9  
**Purity:**  $\geq$ 98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years



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

### Laboratory Procedures

1 $\alpha$ -Hydroxymidazolam  $\beta$ -D-Glucuronide is supplied as a crystalline solid.

### Description

1 $\alpha$ -Hydroxymidazolam  $\beta$ -D-Glucuronide is the glucuronidated conjugate of the midazolam metabolite, 1'-hydroxymidazolam (Item No. 10385), that is mainly excreted in urine. O- and N-glucuronidation of 1'-hydroxymidazolam is mediated by the UDP-glucuronosyltransferases UGT2B4/2B7 and 1A4, respectively.<sup>1</sup>

### Reference

1. Seo, K.-A., Bae, S.K., Choi, Y.-K., *et al.* Metabolism of 1'- and 4-hydroxymidazolam by glucuronide conjugation is largely mediated by UDP-glucuronosyltransferases 1A4, 2B4, and 2B7. *Drug Metab. Dispos.* **38(11)**, 2007-2013 (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, 08/01/2023

#### 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