

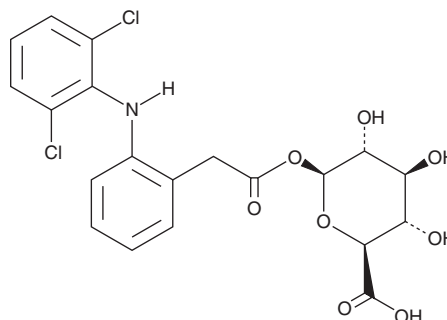
# PRODUCT INFORMATION



## Diclofenac Acyl-β-D-Glucuronide

Item No. 22455

**CAS Registry No.:** 64118-81-6  
**Formal Name:** β-D-glucopyranuronic acid, 1-[2-[(2,6-dichlorophenyl)amino]benzeneacetate]  
**Synonyms:** Diclofenac Glucuronide, Diclofenac-β-D-glucuronide  
**MF:** C<sub>20</sub>H<sub>19</sub>Cl<sub>2</sub>NO<sub>8</sub>  
**FW:** 472.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 277 nm  
**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

Diclofenac acyl-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the diclofenac acyl-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Diclofenac acyl-β-D-glucuronide is soluble in methanol and DMSO.

### Description

Diclofenac acyl-β-D-glucuronide is an active metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor diclofenac (Item Nos. 70680 | 22983).<sup>1</sup> It is formed from diclofenac via UDP-glucuronosyltransferase (UGT) isoform UGT2B7.<sup>2</sup> Diclofenac acyl-β-D-glucuronide inhibits the proliferation of HT-29 colorectal adenocarcinoma cells (IC<sub>50</sub> = 130 μM).<sup>3</sup>

### References

1. Kenny, J.R., Maggs, J.L., Meng, X., *et al.* Syntheses and characterization of the acyl glucuronide and hydroxy metabolites of diclofenac. *J. Med. Chem.* **47(11)**, 2816-2825 (2004).
2. King, C., Tang, W., Ngui, J., *et al.* Characterization of rat and human UDP-glucuronosyltransferases responsible for the in vitro glucuronidation of diclofenac. *Toxicol. Sci.* **61(1)**, 49-53 (2001).
3. Cannell, G.R., Vesey, D.A., and Dickinson, R.G. Inhibition of proliferation of HT-29 colon adenocarcinoma cells by carboxylate NSAIDs and their acyl glucuronides. *Life Sci.* **70(1)**, 37-48 (2001).

#### 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, 09/28/2022

#### 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