

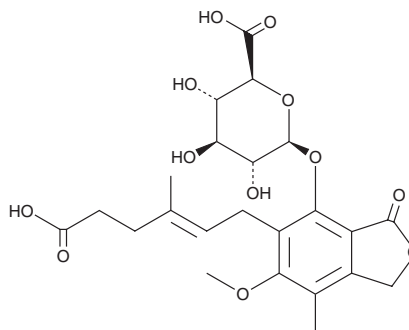
# PRODUCT INFORMATION



## Mycophenolic Acid $\beta$ -D-Glucuronide

Item No. 19078

**CAS Registry No.:** 31528-44-6  
**Formal Name:** 5-[(2E)-5-carboxy-3-methyl-2-penten-1-yl]-1,3-dihydro-6-methoxy-7-methyl-3-oxo-4-isobenzofuranyl  $\beta$ -D-glucopyranosiduronic acid  
**Synonyms:** MPAG, MPA-phenyl-glucuronide, Mycophenolic Acid Glucosiduronate  
**MF:**  $C_{23}H_{28}O_{12}$   
**FW:** 496.5  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 214, 250, 292 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Mycophenolic acid  $\beta$ -D-glucuronide (MPAG) is supplied as a crystalline solid. A stock solution may be made by dissolving the MPAG in the solvent of choice, which should be purged with an inert gas. MPAG is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MPAG in these solvents is approximately 10 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MPAG can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of MPAG in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Mycophenolic acid  $\beta$ -D-glucuronide (MPAG) is a metabolite of the immunosuppressant mycophenolic acid (Item No. 21716).<sup>1,2</sup> It is formed from mycophenolic acid by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A9, UGT1A8, UGT1A7, and UGT1A10 in insect cell-derived supersomes expressing the human enzymes.<sup>2</sup> Unlike mycophenolic acid, MPAG does not inhibit IMP dehydrogenase (IMPDH) type II in a cell-free assay.<sup>1</sup> However, it does reduce tumor growth in an Ehrlich murine spontaneous adenocarcinoma model when administered at a dose of 6 mg/animal.<sup>3</sup>

### References

1. Gensburger, O., Picard, N., and Marquet, P. Effect of mycophenolate acyl-glucuronide on human recombinant type 2 inosine monophosphate dehydrogenase. *Clin. Chem.* **55(5)**, 986-993 (2009).
2. Picard, N., Ratanasavanh, D., Prémaud, A., et al. Identification of the UDP-glucuronosyltransferase isoforms involved in mycophenolic acid phase II metabolism. *Drug Metab. Dispos.* **33(10)**, 139-146 (2005).
3. Ando, K., Suzuki, S., and Arita, M. Synthesis of mycophenolic acid  $\beta$ -D-glucuronide and its antitumor activity. *J. Antibiot. (Tokyo)* **23(8)**, 408-413 (1970).

#### 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, 02/15/2024

#### 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