

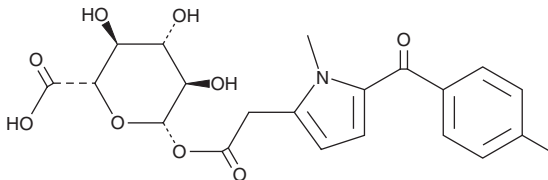
# PRODUCT INFORMATION



## Tolmetin $\beta$ -D-Glucuronide

Item No. 22161

CAS Registry No.: 71595-19-2  
Formal Name: 1-[1-methyl-5-(4-methylbenzoyl)-1H-pyrrole-2-acetate]  $\beta$ -D-glucopyranuronic acid  
MF:  $C_{21}H_{23}NO_9$   
FW: 433.4  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Tolmetin  $\beta$ -D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the tolmetin  $\beta$ -D-glucuronide in the solvent of choice, which should be purged with an inert gas. Tolmetin  $\beta$ -D-glucuronide is slightly soluble in DMSO and methanol (warmed).

### Description

Tolmetin  $\beta$ -D-glucuronide is a major phase II metabolite of tolmetin (Item No. 18195).<sup>1</sup> Tolmetin  $\beta$ -D-glucuronide has been classified into a 'warning' category in terms of its potential to cause idiosyncratic drug toxicity, based on its half-life in several *in vitro* systems.<sup>2</sup>

### References

1. Hyneck, M.L., Smith, P.C., Unseld, E., *et al.* High-performance liquid chromatographic determination of tolmetin, tolmetin glucuronide and its isomeric conjugates in plasma and urine. *J. Chromatogr.* **420**(2), 349-356 (1987).
2. Sawamura, R., Okudaira, N., Watanabe, K., *et al.* Predictability of idiosyncratic drug toxicity risk for carboxylic acid-containing drugs based on the chemical stability of acyl glucuronide. *Drug. Metab. Dispos.* **38**(10), 1857-1864 (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, 12/16/2022

#### 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