# **PRODUCT** INFORMATION



p-Cresol Glucuronide

Item No. 35113

CAS Registry No.: Formal Name:	17680-99-8 4-methylphenyl β-D-	
Synonyms:	4-Cresol Glucuronide, <i>p</i> -Tolyl-β-D-glucuronide	00 00
MF:	C <sub>13</sub> H <sub>16</sub> O <sub>7</sub>	
FW:	284.3	
Purity:	≥90%	HO* VH
UV/Vis.:	λ <sub>max</sub> : 220 nm	ÓН
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	
Information concerns the product experifications. Patch experific analytical results are provided on each contificate of an		

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

## Laboratory Procedures

p-Cresol glucuronide is supplied as a solid. A stock solution may be made by dissolving the p-cresol glucuronide in the solvent of choice, which should be purged with an inert gas. p-Cresol glucuronide is soluble in ethanol.

## Description

p-Cresol glucuronide is a metabolite of p-cresol, a uremic toxin formed by bacterial fermentation of proteins in the large intestine.<sup>1</sup> It is formed from p-cresol primarily by the UDP-glucuronosyltransferase (UGT) isoform UGT1A6, but also by UGT1A9, in human liver and kidney microsomes. p-Cresol glucuronide accumulates in feces, blood, and urine of patients with chronic kidney disease (CKD).<sup>2</sup>

## References

- 1. Rong, Y. and Kiang, T.K.L. Characterizations of human UDP-glucuronosyltransferase enzymes in the conjugation of *p*-cresol. Toxicol. Sci. 176(2), 285-296 (2020).
- 2. Gryp, T., De Paepe, K., Vanholder, R., et al. Gut microbiota generation of protein-bound uremic toxins and related metabolites is not altered at different stages of chronic kidney disease. Kidney Int. 97(6), 1230-1242 (2020).

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.

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, 10/21/2022

## 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