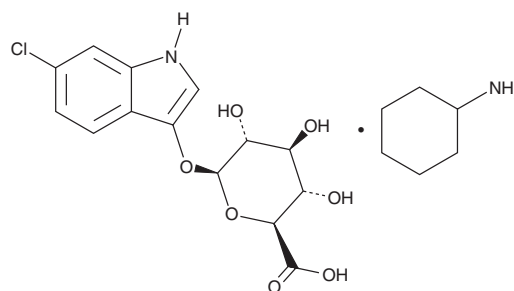


PRODUCT INFORMATION

6-Chloro-3-indolyl-β-D-Glucuronide (cyclohexylammonium salt)

Item No. 30131

CAS Registry No.: 138182-20-4
Formal Name: 6-chloro-1H-indol-3-yl, β-D-glucopyranosiduronic acid, compd. with cyclohexanamine
MF: C₁₄H₁₄ClNO₇ • C₆H₁₃N
FW: 442.9
Purity: ≥95%
UV/Vis.: λ_{max}: 231 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

6-Chloro-3-indolyl-β-D-glucuronide (cyclohexylammonium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-chloro-3-indolyl-β-D-glucuronide (cyclohexylammonium salt) in the solvent of choice, which should be purged with an inert gas. 6-Chloro-3-indolyl-β-D-glucuronide (cyclohexylammonium salt) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

6-Chloro-3-indolyl-β-D-glucuronide (cyclohexylammonium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 6-chloro-3-indolyl-β-D-glucuronide (cyclohexylammonium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 6-Chloro-3-indolyl-β-D-Glucuronide (cyclohexylammonium salt) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

6-Chloro-3-indolyl-β-D-glucuronide is a chromogenic substrate for β-glucuronidase.¹ Upon cleavage by β-glucuronidase, an insoluble salmon precipitate is formed that can be used in histochemical analysis of β-glucuronidase activity.

Reference

1. Wozniak, C.A. and Owens, L.D. Hydrolysis of glucuronide-based substrates mediated by tungsten, Cu²⁺, Fe²⁺, and Zn²⁺. *Physiologia Plantarum* **96(3)**, 484-490 (1996).

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, 05/12/2020

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