

PRODUCT INFORMATION



4-hydroxy Amphetamine Glucuronide (trifluoroacetate salt)

Item No. 9002853

Formal Name: 4-(2-aminopropyl)phenyl-β-D-glucopyranosiduronic acid, trifluoroacetate salt

Synonyms: 4-HA Glucuronide, p-hydroxy Amphetamine, para-hydroxy Amphetamine, POHAG

MF: C₁₅H₂₁NO₇ • CF₃COOH

FW: 441.4

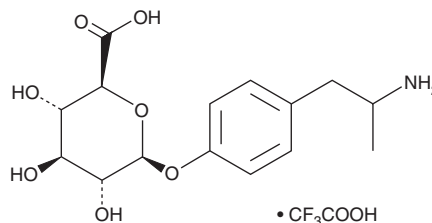
Purity: ≥98%

UV/Vis.: λ_{max}: 221, 272, 278 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 1 year from the QC date provided on the Certificate of Analysis, when stored properly



Description

4-hydroxy Amphetamine glucuronide (Item No. 9002853) is an analytical reference standard that is structurally categorized as an amphetamine metabolite. It is the glucuronide of 4-hydroxy Amphetamine (Item No. 9001830), a bioactive metabolite of amphetamine. 4-hydroxy Amphetamine glucuronide is the primary metabolite of amphetamine found in bile of rats.¹ It may be found in the urine of female rats and guinea pigs but is not found in human or male rat urine.² This product is intended for research and forensic applications.

References

1. Caldwell, J., Dring, L.G., and Williams, R.T. Biliary excretion of amphetamine and methamphetamine in the rat. *Biochem. J.* **129**(1), 25-29 (1972).
2. Sever, P.S., Dring, L.G., and Williams, R.T. Urinary metabolites of p-hydroxyamphetamine in man, rat and guinea-pig. *Xenobiotica* **6**(6), 345-353 (1976).

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, 12/13/2016

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