

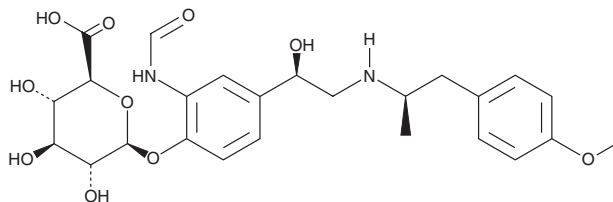
PRODUCT INFORMATION



Formoterol O-β-D-Glucuronide

Item No. 27805

CAS Registry No.: 615551-59-2
Formal Name: 2-(formylamino)-4-[(1R)-1-hydroxy-2-[[[(1R)-2-(4-methoxyphenyl)-1-methylethyl]amino]ethyl]phenyl, β-D-glucopyranosiduronic acid
MF: C₂₅H₃₂N₂O₁₀
FW: 520.5
Purity: ≥95% (mixture of diastereomers)
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Formoterol O-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the formoterol O-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Formoterol O-β-D-glucuronide is slightly soluble in methanol and DMSO.

Description

Formoterol O-β-D-glucuronide is a metabolite of formoterol (Item No. 15584).¹

Reference

1. Rosenberg, J., Larsson, P., Tegnér, K., et al. Mass balance and metabolism of [³H]Formoterol in healthy men after combined i.v. and oral administration-mimicking inhalation. *Drug Metab. Dispos.* **27(10)**, 1104-1116 (1999).

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, 11/10/2022

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