

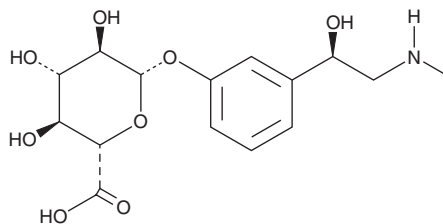
PRODUCT INFORMATION



Phenylephrine Glucuronide

Item No. 26647

CAS Registry No.: 2021255-73-0
Formal Name: (2S,3S,4S,5R,6S)-3,4,5-trihydroxy-6-(3-((R)-1-hydroxy-2-(methylamino)ethyl)phenoxy)tetrahydro-2H-pyran-2-carboxylic acid
MF: C₁₅H₂₁NO₈
FW: 343.3
Purity: ≥95%
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

Phenylephrine glucuronide is supplied as a solid. A stock solution may be made by dissolving the phenylephrine glucuronide in water. We do not recommend storing the aqueous solution for more than one day.

Description

Phenylephrine glucuronide is a metabolite of the α_{1A} -adrenergic receptor agonist phenylephrine (Item Nos. 18619 | 17205).¹

Reference

1. Karaman, R., Karaman, D., and Zeiadeh, I. Computationally-designed phenylephrine prodrugs - a model for enhancing bioavailability. *Mol. Phys.* **111**(21), 3249-3264 (2013).

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, 12/19/2022

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