

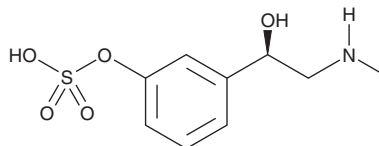
PRODUCT INFORMATION



Phenylephrine-3-O-Sulfate

Item No. 26620

CAS Registry No.: 1242184-39-9
Formal Name: α R-[(methylamino)methyl]-3-(sulfooxy)-benzenemethanol
MF: $C_9H_{13}NO_5S$
FW: 247.3
Purity: $\geq 85\%$
Supplied as: An oil
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Phenylephrine-3-O-sulfate is supplied as an oil. A stock solution may be made by dissolving the phenylephrine-3-O-sulfate in the solvent of choice, which should be purged with an inert gas. Phenylephrine-3-O-sulfate is soluble in the organic solvent DMSO.

Description

Phenylephrine-3-O-sulfate is metabolite of the α_{1A} -adrenergic receptor agonist phenylephrine (Item Nos. 18619 | 17205) formed *via* sulfation.¹

Reference

1. Shah, H.N., Halquist, M.T., and Gerk, P.M. Direct detection of phenylephrine 3-O-sulfate in LS180 human intestinal cells using a novel hydrophilic interaction liquid chromatography (HILIC) assay. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1040**, 67-72 (2017).

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, 11/29/2022

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