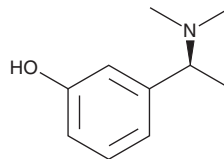


PRODUCT INFORMATION



NAP 226-90
Item No. 30124

CAS Registry No.: 139306-10-8
Formal Name: 3-[(1S)-1-(dimethylamino)ethyl]-phenol
MF: C₁₀H₁₅NO
FW: 165.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218 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

NAP 226-90 is supplied as a crystalline solid. A stock solution may be made by dissolving the NAP 226-90 in the solvent of choice, which should be purged with an inert gas. NAP 226-90 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NAP 226-90 in these solvents is approximately 30 mg/ml.

Description

NAP 226-90 is a metabolite of rivastigmine (Item No. 14270) that is formed by hydrolysis.¹

Reference

1. Polinsky, R.J. Clinical pharmacology of rivastigmine: A new-generation acetylcholinesterase inhibitor for the treatment of Alzheimer's disease. *Clin. Ther.* **20(4)**, 634-647 (1998).

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, 05/18/2020

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