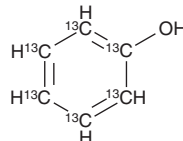


PRODUCT INFORMATION



Phenol-¹³C₆
Item No. 30526

CAS Registry No.: 89059-34-7
Formal Name: Phenol-1,2,3,4,5,6-¹³C₆
MF: [¹³C]₆H₆O
FW: 100.1
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 273 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

Phenol-¹³C₆ is supplied as a crystalline solid. A stock solution may be made by dissolving the phenol-¹³C₆ in the solvent of choice, which should be purged with an inert gas. Phenol-¹³C₆ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of phenol-¹³C₆ in these solvents is approximately 30 mg/ml.

Phenol-¹³C₆ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, phenol-¹³C₆ should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Phenol-¹³C₆ has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Phenol-¹³C₆ is a building block.¹ It has been used in the synthesis of ¹³C₆-labeled metabolites of (±)-asenapine (Item No. 9000496).

Reference

1. Kuethe, J.T. Synthesis of stable isotope-labeled metabolites of asenapine. *J. Labelled Comp. Radiopharm.* 55(5), 180-185 (2012).

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, 05/01/2020

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