

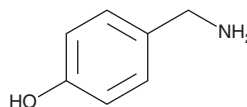
PRODUCT INFORMATION



4-HOBA

Item No. 25398

CAS Registry No.: 696-60-6
Formal Name: 4-(aminomethyl)-phenol
Synonyms: 4-(Aminomethyl)phenol,
4-Hydroxybenzylamine,
NSC 125720
MF: C₇H₉NO
FW: 123.2
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 277 nm
Supplied as: A crystalline 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

4-HOBA is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-HOBA in the solvent of choice, which should be purged with an inert gas. 4-HOBA is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 4-HOBA in these solvents is approximately 30 mg/ml.

4-HOBA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-HOBA should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4-HOBA has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-HOBA is an isomer of the isoketal scavenger 2-HOBA (Item No. 25357) that has reduced efficacy as an isoketal scavenger.¹ Unlike 2-HOBA, 4-HOBA has no effect on hypertension induced by angiotensin II (Item No. 17150) in mice.¹ 4-HOBA has been used as a negative control for the activity of 2-HOBA and related compounds in a mouse model of hypertension.¹

Reference

1. Kirabo, A., Fontana, V., de Faria, A.P.C., *et al.* DC isoketal-modified proteins activate T cells and promote hypertension. *J. Clin. Invest.* **124**(10), 4642-4656 (2014).

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, 11/03/2022

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