

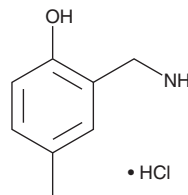
PRODUCT INFORMATION



5-methyl-2-HOBA (hydrochloride)

Item No. 25351

CAS Registry No.: 2044714-53-4
Formal Name: 2-(aminomethyl)-4-methyl-phenol, monohydrochloride
Synonyms: 5-Me-2-HOBA, 5-methyl-2-Hydroxybenzylamine
MF: C₈H₁₁NO • HCl
FW: 173.6
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 286 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

5-methyl-2-HOBA (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-methyl-2-HOBA (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 5-methyl-2-HOBA (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-methyl-2-HOBA (hydrochloride) in these solvents is approximately 0.25, 2.5, and 0.3 mg/ml, respectively.

5-methyl-2-HOBA (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-methyl-2-HOBA (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 5-methyl-2-HOBA (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

5-methyl-2-HOBA is an isoketal scavenger.¹ It reduces angiotensin II-induced increases in systolic blood pressure in mice.

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, 11/21/2022

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