

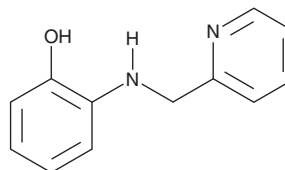
PRODUCT INFORMATION



2-PMAP

Item No. 21021

CAS Registry No.: 102212-26-0
Formal Name: 2-[(2-pyridinylmethyl)amino]-phenol
MF: C₁₂H₁₂N₂O
FW: 200.2
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 293 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

2-PMAP is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-PMAP in the solvent of choice. 2-PMAP is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of 2-PMAP in these solvents is approximately 30 mg/ml.

2-PMAP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-PMAP should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-PMAP has a solubility of approximately 0.33 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

2-PMAP is an inhibitor of the production of amyloid precursor protein (APP) and amyloid-β (Aβ) secretion.¹ It reduces levels of APP and its intracellular C-terminal fragments α-CTF and β-CTF in CHO APP751SW cells when used at a concentration of 10 μM. 2-PMAP also reduces secretion of Aβ (1-40) and Aβ (1-42) by 88.6 and 84.9%, respectively, in CHO APP751SW cells when used at a concentration of 50 μM without inducing cellular toxicity. *In vivo*, 2-PMAP reduces brain levels of full-length APP, α-CTF and β-CTF, and soluble Aβ (1-40) and Aβ (1-42) in the APP_{SW}PS1_{dE9} transgenic mouse model of Alzheimer's disease. It also improves novel object memory as well as spatial memory in the radial arm maze in APP_{SW}PS1_{dE9} transgenic mice.

Reference

1. Asuni, A.A., Guridi, M., Pankiewicz, J.E., *et al.* Modulation of APP expression reduces Aβ deposition in a mouse model. *Ann. Neurol.* **75**(5), 684-699 (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, 12/22/2022

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