

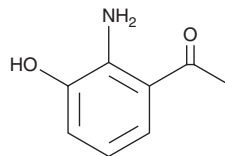
PRODUCT INFORMATION



1-(2-Amino-3-hydroxyphenyl)ethanone

Item No. 36416

CAS Registry No.: 4502-10-7
Formal Name: 1-(2-amino-3-hydroxyphenyl)-ethanone
MF: C₈H₉NO₂
FW: 151.2
Purity: ≥98%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A 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

1-(2-Amino-3-hydroxyphenyl)ethanone is supplied as a solid. A stock solution may be made by dissolving the 1-(2-amino-3-hydroxyphenyl)ethanone in the solvent of choice, which should be purged with an inert gas. 1-(2-Amino-3-hydroxyphenyl)ethanone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1-(2-amino-3-hydroxyphenyl)ethanone in DMSO is approximately 1 mg/ml and approximately 2 mg/ml in DMF. 1-(2-Amino-3-hydroxyphenyl)ethanone is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1-(2-amino-3-hydroxyphenyl)ethanone can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1-(2-amino-3-hydroxyphenyl)ethanone in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1-(2-Amino-3-hydroxyphenyl)ethanone is an active metabolite of the quorum sensing-regulated molecule 2-amino acetophenone.¹ It increases survival in a mouse model of *P. aeruginosa*-induced post-burn infection when administered at a dose of 7.6 mg/kg.

Reference

1. Bandyopadhaya, A., Kesarwani, M., Que, Y.-A., *et al.* The quorum sensing volatile molecule 2-amino acetophenone modulates host immune responses in a manner that promotes life with unwanted guests. *PLoS Pathog.* **8(11)**, e1003024 (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, 02/09/2023

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