

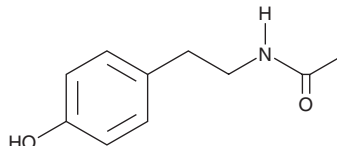
PRODUCT INFORMATION



N-Acetyltyramine

Item No. 9001373

CAS Registry No.: 1202-66-0
Formal Name: N-[2-(4-hydroxyphenyl)ethyl]-acetamide
MF: $C_{10}H_{13}NO_2$
FW: 179.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

N-acetyltyramine is supplied as a solid. A stock solution may be made by dissolving the N-acetyltyramine in the solvent of choice, which should be purged with an inert gas. N-acetyltyramine is soluble in methanol, ethanol, DMSO, and dimethyl formamide.

Description

N-acetyltyramine is a metabolite of the biogenic amine tyramine (Item No. 18601).¹ It enhances cytotoxicity of doxorubicin (Item No. 15007) in resistant P388 murine leukemia cells with an IC_{50} value of 0.13 $\mu g/ml$ compared with an IC_{50} value of 0.48 $\mu g/ml$ for doxorubicin alone.

Reference

1. Kunimoto, S., Xu, C.Z., Naganawa, H., *et al.* Reversal of resistance by N-acetyltyramine or N-acetyl-2-phenylethylamine in doxorubicin-resistant leukemia P388 cells. *J. Antibiot. (Tokyo)* **40(11)**, 1651-1652 (1987).

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, 06/20/2018

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