

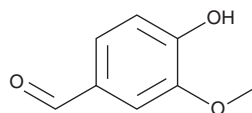
PRODUCT INFORMATION



Vanillin

Item No. 36422

CAS Registry No.: 121-33-5
Formal Name: 4-hydroxy-3-methoxy-benzaldehyde
Synonyms: 3-methoxy-4-Hydroxybenzaldehyde, NSC 15351, NSC 403658, NSC 48383
MF: C₈H₈O₃
FW: 152.1
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 279, 308 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Vanillin (Item No. 36422) is an analytical reference standard categorized as a precursor in the synthesis of 3,4-MDMA (Item Nos. ISO60190 | 13971).¹ Vanillin is used as a flavoring additive in vaping mixtures for electronic cigarettes.² This product is intended for research and forensic applications.

References

1. Gallagher, R.T., Shimmon, R., and McDonagh, A.M. Synthesis and impurity profiling of MDMA prepared from commonly available starting materials. *Forensic Sci. Int.* **223(1-3)**, 306-313 (2012).
2. Erythropel, H.C., Jabba, S.V., DeWinter, T.M., et al. Formation of flavorant-propylene glycol adducts with novel toxicological properties in chemically unstable e-cigarette liquids. *Nicotine Tob. Res.* **21(9)**, 1248-1258 (2019).

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, 07/07/2023

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