

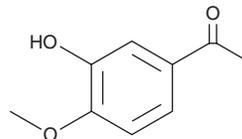
PRODUCT INFORMATION



3-Hydroxy-4-methoxyacetophenone

Item No. 36801

CAS Registry No.: 6100-74-9
Formal Name: 1-(3-hydroxy-4-methoxyphenyl)-ethanone
Synonyms: Acetoisovanillone, NSC 30050
MF: C₉H₁₀O₃
FW: 166.2
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 272 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

3-Hydroxy-4-methoxyacetophenone is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy-4-methoxyacetophenone in the solvent of choice, which should be purged with an inert gas. 3-Hydroxy-4-methoxyacetophenone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-hydroxy-4-methoxyacetophenone in ethanol is approximately 10 mg/ml and approximately 2 mg/ml in DMSO and DMF.

Description

3-Hydroxy-4-methoxyacetophenone is a building block.^{1,2} It has been used in the synthesis of anti-inflammatory and antiviral agents.

References

1. Yu, Y.-Y., Li, X.-Q., Hu, W.-P., *et al.* Self-developed NF-κB inhibitor 270 protects against LPS-induced acute kidney injury and lung injury through improving inflammation. *Biomed. Pharmacother.* **147**, 112615 (2022).
2. Liu, J., Ma, L., Song, C., *et al.* Anti-HIV effects of baculiferins are regulated by the potential target protein DARS. *ACS Chem. Biol.* **16(8)**, 1377-1389 (2021).

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, 12/21/2022

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