

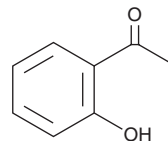
PRODUCT INFORMATION



2'-Hydroxyacetophenone

Item No. 37418

CAS Registry No.: 118-93-4
Formal Name: 1-(2-hydroxyphenyl)-ethanone
Synonym: 1-(2-Hydroxyphenyl)ethenone, NSC 9263, NSC 16933, NSC 44452, NSC 83568
MF: C₈H₈O₂
FW: 136.2
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 251 nm
Supplied as: A neat liquid
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'-Hydroxyacetophenone is supplied as a neat liquid. A stock solution may be made by dissolving the 2'-hydroxyacetophenone in the solvent of choice, which should be purged with an inert gas. 2'-Hydroxyacetophenone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2'-hydroxyacetophenone in these solvents is approximately 11, 20, and 16 mg/ml, respectively.

Description

2'-Hydroxyacetophenone is an acetophenone synthetic building block.¹ It has been used in the synthesis of hydroxychalcone derivatives with inhibitory activity towards COX-2.²

References

1. Albayrak, C., Gümrükçüoğlu, I., Odabaşoğlu, M., *et al.* Synthesis, spectroscopic, and molecular structure characterizations of some azo derivatives of 2-hydroxyacetophenone. *J. Mol. Struct.* **932(1-3)**, 43-54 (2009).
2. Tran, T.-D., Park, H., Kim, H.P., *et al.* Inhibitory activity of prostaglandin E₂ production by the synthetic 2'-hydroxychalcone analogues: Synthesis and SAR study. *Bioorg. Med. Chem. Lett.* **19(6)**, 1650-1653 (2009).

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, 03/20/2023

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