

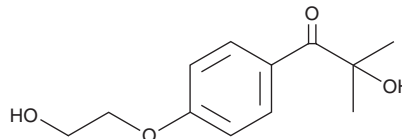
PRODUCT INFORMATION



2-Hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone

Item No. 36428

CAS Registry No.: 106797-53-9
Formal Name: 2-hydroxy-1-[4-(2-hydroxyethoxy)phenyl]-2-methyl-1-propanone
MF: C₁₂H₁₆O₄
FW: 224.3
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 275 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

2-Hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone in the solvent of choice, which should be purged with an inert gas. 2-Hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone in these solvents is approximately 30 mg/ml.

Description

2-Hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone is a photoinitiator.^{1,2} Upon UV activation, 2-hydroxy-4'-(2-hydroxyethoxy)-2-methylpropiophenone dissociates into benzoyl and ketyl radicals, which induce polymerization. It has been used in the synthesis of hyperbranched polymers for various industrial applications, as well as in the synthesis of noble metal nanoparticles.

References

1. Choi, G., Oh, Y., Jeong, S., *et al.* Synthesis of renewable, recyclable, degradable thermosets endowed with highly branched polymeric structures and reinforced with carbon fibers. *Macromol.* **56(6)**, 2526-2535 (2023).
2. Liu, M., Li, M.-D., Xue, J., *et al.* Time-resolved spectroscopic and density functional theory study of the photochemistry of Irgacure-2959 in an aqueous solution. *J. Phys. Chem. A* **118(38)**, 8701-8707 (2013).

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, 04/12/2023

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