

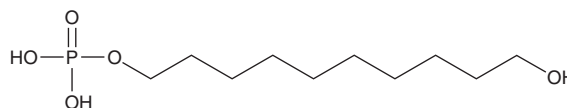
PRODUCT INFORMATION



10-(Phosphonoxy)-1-decanole

Item No. 9004346

CAS Registry No.: 235097-47-9
Formal Name: 1-(dihydrogen phosphate) 1,10-decanediol
Synonyms: 10-Hydroxydecyl dihydrogen phosphate, HDP
MF: C₁₀H₂₃O₅P
FW: 254.3
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

10-(Phosphonoxy)-1-decanole is a hydrolytic degradation product of 10-methacryloyloxydecyl dihydrogen phosphate-N-methylacryloyl glycine (MDP-NMGly) dental primers.¹

Reference

1. Ma, S. Development of a self-etching primer with higher shelf life and greater dentin bond stability. *Dent. Mater. J.* **29(1)**, 59-67 (2010).

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, 02/13/2024

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