

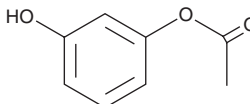
PRODUCT INFORMATION



Resorcinol Monoacetate (technical grade)

Item No. 39019

CAS Registry No.: 102-29-4
Formal Name: 1-acetate, 1,3-benzenediol
Synonyms: Acetylresorcinol, NSC 40511
MF: $C_8H_8O_3$
FW: 152.1
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Resorcinol monoacetate (technical grade) is supplied as a neat oil. A stock solution may be made by dissolving the resorcinol monoacetate (technical grade) in the solvent of choice, which should be purged with an inert gas. Resorcinol monoacetate (technical grade) is soluble in DMSO.

Description

Resorcinol monoacetate is a phenol acetate with antibacterial activity.¹ It increases collagen cross-linking and prevents *B. subtilis* growth on sponge surfaces when attached to collagen sponges at 2 or 5% (w/w). Resorcinol monoacetate has been used as a starting material in the synthesis of $\alpha V\beta 3$ and $\alpha V\beta 5$ integrin dual inhibitors.²

References

1. Gunasekaran, S. and Chvapil, M. Collagen based biomaterials: An ideal way of increasing their resistance to infection. *Biomater. Artif. Cells Artif. Organs* **16**(4), 771-783 (1988).
2. Marugán, J.J., Manthey, C., Anaclerio, B., et al. Design, synthesis, and biological evaluation of novel potent and selective $\alpha V\beta 3/\alpha V\beta 5$ integrin dual inhibitors with improved bioavailability. Selection of the molecular core. *J. Med. Chem.* **48**(4), 926-934 (2005).

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, 05/28/2024

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