

PRODUCT INFORMATION



5-deoxy-D-Ribose

Item No. 30103

CAS Registry No.: 13039-75-3

MF: $C_5H_{10}O_4$

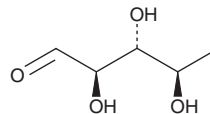
FW: 134.1

Purity: $\geq 95\%$

Supplied as: A liquid

Storage: $-20^\circ C$

Stability: ≥ 2 years



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

Laboratory Procedures

5-deoxy-D-Ribose is supplied as a solid. A stock solution may be made by dissolving the 5-deoxy-D-ribose in the solvent of choice, which should be purged with an inert gas. 5-deoxy-D-Ribose is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-deoxy-D-ribose in these solvents is approximately 15 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 5-deoxy-D-ribose can be prepared by directly dissolving the liquid in aqueous buffers. The solubility of 5-deoxy-D-ribose in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-deoxy-D-Ribose is a building block that has been used in the synthesis of carcinolytic compounds.¹

Reference

1. Shunk, C.H., Lavigne, J.B., and Folkers, K. Studies on carcinolytic compounds. V. 6,7-Dimethyl-9-[1'-(5-desoxy-D-ribityl)]-isoalloxazine. *J. Am. Chem. Soc.* **77(8)**, 2210-2212 (1955).

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, 05/12/2020

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