

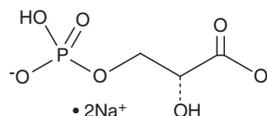
PRODUCT INFORMATION



D-(-)-3-Phosphoglyceric Acid (sodium salt)

Item No. 20123

CAS Registry No.: 80731-10-8
Formal Name: 2R-hydroxy-3-(phosphonoxy)-
propanoic acid, disodium salt
MF: C₃H₅O₇P • 2Na
FW: 230.0
Purity: ≥90%
Supplied as: A crystalline 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

D-(-)-3-Phosphoglyceric acid (sodium salt) is supplied as a crystalline solid. Aqueous solutions of D-(-)-3-phosphoglyceric acid (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of D-(-)-3-phosphoglyceric acid (sodium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

D-(-)-3-Phosphoglyceric acid is an intermediate in several biological pathways, most notably glycolysis/gluconeogenesis and the biosynthesis of serine, glycine, and threonine.¹ It also has numerous roles as an intermediate in biosynthetic pathways in plants, eukaryotes, and prokaryotes.

Reference

1. G. Zubay, Principles of Biochemistry, Second Edition. Brown (1995).

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, 05/26/2023

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