

PRODUCT INFORMATION

Glycerophosphate (sodium salt hydrate)

Item No. 14405

Formal Name: 1,2,3-propanetriol, 2-(dihydrogen phosphate), 1,2,3-propanetriol, (2R)1-(dihydrogen phosphate), disodium salt, pentahydrate

Synonym: Glycerol Phosphate

MF: $C_3H_7O_6P \cdot 2Na [5H_2O]$

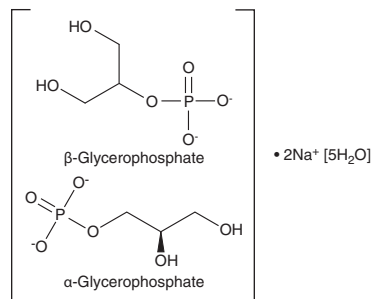
FW: 306.1

Purity: $\geq 95\%$ (mixture of α and β)

Supplied as: A crystalline solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Glycerophosphate is a mixture of the endogenous metabolite α -glycerophosphate and the alkaline phosphatase (ALP) inhibitor β -glycerophosphate (Item No. 38367).

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, 05/12/2023

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