PRODUCT INFORMATION



β-hydroxy Pyruvic Acid (lithium salt hydrate)

Item No. 21856

Formal Name:	3-hydroxy-2-oxo-propanoic acid, monolithium salt, hydrate	
Synonym:	3-hydroxy Pyruvic Acid	0
MF:	$C_3H_4O_4 \bullet Li [XH_2O]$	
FW:	110.0	-O TO OH
Purity:	≥95%	0
Supplied as:	A crystalline solid	• Li ⁺ [XH ₂ O]
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

β-hydroxy Pyruvic acid (lithium salt hydrate) is supplied as a crystalline solid. Aqueous solutions of β -hydroxy pyruvic acid (lithium salt hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of β -hydroxy pyruvic acid (lithium 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

 β -hydroxy Pyruvic acid is a metabolite in the process of photorespiration in plants.¹ In peroxisomes, hydroxypyruvate is reduced to glycerate by NADH.^{2,3}

References

- 1. Stafford, H.A., Magaldi, A., and Vennesland, B. The enzymatic reduction of hydroxypyruvic acid to D-glyceric acid in higher plants. J. Biol. Chem. 207(2), 621-629 (1954).
- 2. Douce, R. and Heldt, H.-W. Photorespiration in Photosynthesis: Physiology and Metabolism. Springer Netherlands, 9, 115-136 (2000).
- 3. Givan, C.V. and Kleczkowski, L.A. The enzymic reduction of glyoxylate and hydroxypyruvate in leaves of higher plants. Plant Physiol. 100(2), 552-556 (1992).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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.

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, 12/22/2022

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