

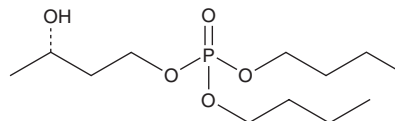
PRODUCT INFORMATION



(S)-Dibutyl 3-Hydroxybutyl Phosphate

Item No. 14787

CAS Registry No.: 2108602-99-7
Formal Name: (S)-dibutyl (3-hydroxybutyl) phosphate
Synonym: (S)-TBP-OH
MF: C₁₂H₂₇O₅P
FW: 282.3
Purity: ≥95%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

(S)-Dibutyl 3-hydroxybutyl phosphate ((S)-TBP-OH) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (S)-TBP-OH in DMF is approximately 25 mg/ml and approximately 15 mg/ml in ethanol and DMSO.

(S)-TBP-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methanolic solution of (S)-TBP-OH should be diluted with the aqueous buffer of choice. The solubility of (S)-TBP-OH in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dibutyl 3-hydroxybutyl phosphate (TBP-OH; Item No. 9001825) is a compound produced from the metabolism of the organophosphorus solvent, tributyl phosphate (TBP). Incubation of goldfish liver microsomes in the presence of NADPH has been shown to convert TBP into TBP-OH and dibutyl phosphate.¹ TBP-OH has also been produced during radiolysis of TBP.² (S)-TBP-OH is an optically active isomer of TBP-OH.

References

1. Sasaki, K., Suzuki, T., Takeda, M., *et al.* Metabolism of phosphoric acid triesters by goldfish, *Carassius auratus*. *Eisei Kagaku* **31(6)**, 397-404 (1985).
2. Lesage, D., Virelizier, H., Jankowski, C.K., *et al.* Identification of minor products obtained during radiolysis of tributylphosphate (TBP). *Spectroscopy* **13(4)**, 275-290 (1997).

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, 07/25/2022

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