

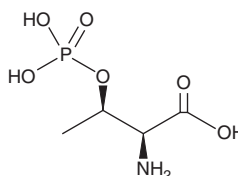
# PRODUCT INFORMATION



## O-Phospho-L-Threonine

Item No. 35057

CAS Registry No.: 1114-81-4  
Synonym: O-Phosphothreonine  
MF:  $C_4H_{10}NO_6P$   
FW: 199.1  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

O-Phospho-L-threonine is supplied as a solid. Aqueous solutions of O-phospho-L-threonine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of O-phospho-L-threonine in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

O-Phospho-L-threonine is a phosphorylated form of the essential amino acid L-threonine (Item No. 30194). It has been used as a substrate in the characterization of phosphatases.<sup>1</sup>

### Reference

1. Polya, G.M. and Wettenhall, R.E. Rapid purification and N-terminal sequencing of a potato tuber cyclic nucleotide binding phosphatase. *Biochim. Biophys. Acta* **1159**(2), 179-184 (1992).

#### 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, 12/08/2022

#### 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