

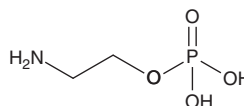
# PRODUCT INFORMATION



## O-Phosphoethanolamine

Item No. 34636

CAS Registry No.: 1071-23-4  
Formal Name: 2-amino-ethanol 1-(dihydrogen phosphate)  
Synonyms: O-Phosphorylethanolamine, NSC 254167  
MF:  $C_2H_8NO_4P$   
FW: 141.1  
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-Phosphoethanolamine is supplied as a solid. Aqueous solutions of O-phosphoethanolamine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of O-phosphoethanolamine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

O-Phosphoethanolamine is an endogenous phospholipid metabolite.<sup>1</sup> It is also a precursor in the biosynthesis of phosphatidylethanolamines in the endoplasmic reticulum. O-Phosphoethanolamine is also incorporated into lipid A, a component of LPS.<sup>2</sup> It decreases disease severity in a mouse model of cerulein-induced acute pancreatitis.<sup>3</sup> Levels of O-phosphoethanolamine are decreased in postmortem amygdala from patients with Parkinson's disease.<sup>4</sup>

### References

1. Podo, F. Tumour phospholipid metabolism. *NMR Biomed.* **12**(7), 413-349 (1999).
2. Samantha, A. and Vrielink, A. Lipid A phosphoethanolamine transferase: Regulation, structure and immune response. *J. Mol. Biol.* **432**(18), 5184-5196 (2020).
3. Sakai, A., Nishiumi, S., Shiomi, Y., et al. Metabolomic analysis to discover candidate therapeutic agents against acute pancreatitis. *Arch. Biochem. Biophys.* **522**(2), 107-120 (2012).
4. Kish, S.J., Rajput, A., Gilbert, J., et al. Elevated  $\gamma$ -aminobutyric acid level in striatal but not extrastriatal brain regions in Parkinson's disease: Correlation with striatal dopamine loss. *Ann. Neurol.* **20**(1), 26-31 (1986).

#### 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, 10/14/2022

#### 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