

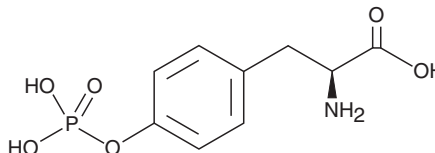
PRODUCT INFORMATION



O-Phospho-L-Tyrosine

Item No. 23064

CAS Registry No.: 21820-51-9
Formal Name: O-phosphono-L-Tyrosine
Synonyms: Phosphotyrosine, P-Tyr
MF: $C_9H_{12}NO_6P$
FW: 261.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 220 nm
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

O-Phospho-L-Tyrosine is supplied as a crystalline solid. Aqueous solutions of o-phospho-L-tyrosine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of o-phospho-L-tyrosine 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-Phospho-L-tyrosine is a phosphorylated form of L-tyrosine. Tyrosine phosphorylation was originally discovered in proteins containing tyrosines found to be modified by tumor viruses.¹ O-phospho-L-tyrosine is used in studies of tyrosine phosphorylation and has been used in affinity chromatography for the purification of IgG.²⁻⁴

References

1. Hunter, T. Phosphotyrosine - a new protein modification *Trends Biochem. Sci.* **7(7)**, 246-249 (1982).
2. Gmeiner, B., and Seelos, C. Phosphorylation of tyrosine prevents dityrosine formation *in vitro*. *FEBS. Lett.* **255(2)**, 395-397 (1989).
3. Afonso, A., Pereira, P., Queiroz, J.A., *et al.* Purification of pre-miR-29 by a new O-phospho-L-tyrosine affinity chromatographic strategy optimized using design of experiments. *J. Chromatogr. A.* **1343**, 119-27 (2014).
4. Pavan, G.L., Bresolin, I.T.L., Grespan, A., *et al.* Phosphorylated-tyrosine based pseudobioaffinity adsorbent for the purification of immunoglobulin G. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1052**, 10-18 (2017).

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, 10/24/2022

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