

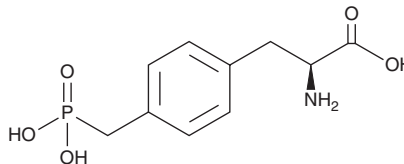
PRODUCT INFORMATION



4-Phosphonomethylphenylalanine

Item No. 45209

CAS Registry No.: 142434-81-9
Formal Name: 4-(phosphonomethyl)-L-phenylalanine
Synonym: *p*-(Phosphonomethyl)phenylalanine
MF: $C_{10}H_{14}NO_5P$
FW: 259.2
Purity: $\geq 95\%$
Supplied as: A 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

4-Phosphonomethylphenylalanine is supplied as a solid. Aqueous solutions of 4-phosphonomethylphenylalanine can be prepared by directly dissolving the solid in aqueous buffers. 4-Phosphonomethylphenylalanine is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

4-Phosphonomethylphenylalanine is a synthetic non-proteinogenic amino acid.¹ It is a phosphatase-stable phosphotyrosine mimetic that has been used in the synthesis of phosphorylated peptides.^{1,2}

References

1. Choi, W.J., Shi, Z.D., Worthy, K.M., et al. Application of azide-alkyne cycloaddition 'click chemistry' for the synthesis of Grb2 SH2 domain-binding macrocycles. *Bioorg. Med. Chem. Lett.* **16**(20), 5265-5269 (2006).
2. McNamara, D.J., Dobrusin, E.M., Zhu, G., et al. Inhibition of binding of phospholipase C γ SH2 domains to phosphorylated epidermal growth factor receptor by phosphorylated peptides. *Int. J. Peptide Protein Res.* **42**, 240-248 (1993).

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, 03/16/2026

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