

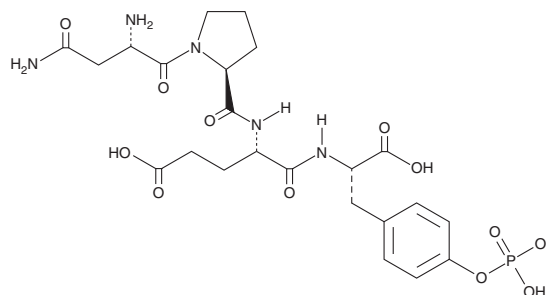
PRODUCT INFORMATION



H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH

Item No. 44048

CAS Registry No.: 290810-63-8
Formal Name: L-asparaginy-L-prolyl-L-α-glutamyl-L-tyrosine, dihydrogen phosphate (ester)
Synonym: Asn-Pro-Glu-Tyr(PO₃H₂)-OH
Peptide Sequence: NPEX-OH (X = Phosphorylated Tyrosine)
MF: C₂₃H₃₂N₅O₁₂P
FW: 601.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH is supplied as a crystalline solid. A stock solution may be made by dissolving the H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH in the solvent of choice, which should be purged with an inert gas. H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH is slightly soluble (0.1-1 mg/ml) in methanol.

H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

H-Asn-Pro-Glu-Tyr(PO₃H₂)-OH is a phosphorylated peptide derivative of the tetramine motif Asn-Pro-Glu-Tyr (NPEY), which is found in the insulin receptor β (InsRβ) submembrane domain and is phosphorylated at Tyr⁹⁷² upon insulin binding.¹

Reference

1. Petersen, M.C.H. and Shulman, G.I. Mechanisms of insulin action and insulin resistance. *Physiol. Rev.* **98**(4), 2133-2223 (2018).

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, 11/04/2025

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