

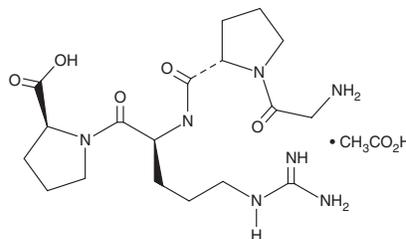
PRODUCT INFORMATION



H-Gly-Pro-Arg-Pro-OH (acetate)

Item No. 21163

CAS Registry No.: 157009-81-9
Formal Name: glycyl-L-prolyl-L-arginyl-L-proline, monoacetate
Synonym: GPRP
MF: C₁₈H₃₁N₇O₅ • C₂H₄O₂
FW: 485.5
Purity: ≥95%
UV/Vis.: λ_{max}: 208, 279 nm
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-Gly-Pro-Arg-Pro-OH (GPRP) (acetate) is supplied as a crystalline solid. Aqueous solutions of GPRP acetate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GPRP acetate in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GPRP is a peptide corresponding partly to the amino terminus of the A α chain and B β chain of fibrin. It inhibits the spontaneous polymerization of fibrin monomers and also inhibits ADP-induced platelet aggregation (IC₅₀ = 3 mM).¹⁻³ GPRP binds to the fibrinogen D domain, inhibiting the ability of fibrinogen to bind to the platelet membrane glycoprotein GPIIb/IIIa.^{1,2}

References

1. Laudano, A.P. and Doolittle, R.F. Synthetic peptide derivatives that bind to fibrinogen and prevent the polymerization of fibrin monomers. *Proc. Natl. Acad. Sci. USA* **75(7)**, 3085-3089 (1978).
2. Plow, E.F. and Marguerie, G. Inhibition of fibrinogen binding to human platelets by the tetrapeptide glycyl-L-prolyl-L-arginyl-L-proline. *Proc. Natl. Acad. Sci. USA* **79(12)**, 3711-3715 (1982).
3. Adelman, B., Gennings, C., Stony, J., et al. Synergistic inhibition of platelet aggregation by fibrinogen-related peptides. *Circ. Res.* **67(4)**, 941-947 (1990).

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, 12/05/2022

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