



## Product Data Sheet

---

<b>Product Name:</b>	Fibrinolysis Inhibiting Factor	
<b>Catalog Number:</b>	AS-24262 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Gly-Pro-Arg-Pro-OH (3-letter code) GPRP (1-letter code)	
<b>Molecular Weight:</b>	425.5	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Fibrinolysis peptide is freely soluble in water.

**Storage:** Fibrinolysis peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This peptide mimics the N terminus of the alpha-chain of fibrin. It binds to the "a" polymerization pocket in the alpha-chain more tightly than a peptide corresponding to the chain native sequence GPRV. The complex between GPRP and fibrinogen causes an increased resistance to degradation by plasmin, and the polymerization reaction is depressed by the addition of excess peptide. Fibrin assembly, as well as the acceleration of the factor XIIIa reaction, can be prevented by this homologue of the natural sequence of amino acids. Ref: Pratt, K. et al. *Proc. Natl. Acad. Sci. USA* **94**, 7176 (1997); Lorand, L. et al. *Proc. Natl. Acad. Sci. USA* **95**, 537 (1998).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
Fibrinogen Binding Inhibitor Peptide HHLGGAKQAGDV	AS-60700	1 mg
Fibrinopeptide A, human ADSGEGDFLAEGGGVR	AS-20735	1 mg
Fibrinopeptide B, human Pyr-GVNDNEEGFFSAR	AS-60697	1 mg

Gamma-Fibrinogen (377-395) AS-62128 1 mg  
YSMKETTMKIIPFNRLSIG

Gamma-Fibrinogen (377-395), scrambled AS-62129 1 mg  
KMMISYTFPIERTGLISNK

*For Research Use Only*