



Product Data Sheet

Product Name: Protease-Activated Receptor-1, PAR-1 Antagonist

Catalog Number: AS-65486 (1 mg)

Sequence: H-Tyr-Phe-Leu-Leu-Arg-Asn-Pro-OH

YFLLRNP-OH

Molecular Weight: 922.1

Peptide Purity: Peak Area by HPLC $\geq 95\%$

Storage: This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description:

The peptide YFLLRNP is an antagonist to thrombin and SFLLRNP (thrombin receptor agonist peptide) in human platelets. YFLLRNP can induce partial activation of human platelets, visible by shape change; but it is unable to fully activate the platelets.

References:

Rasmussen, UB. et al. *J Biol Chem* **268**, 14322 (1993).

For Research Use Only