



## Product Data Sheet

---

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

**Catalog Number:** AS-65487 (1 mg)

**Sequence:** Ser-parafluoroPhe-Aad-Leu-Arg-Asn-Pro-NH<sub>2</sub>

S-F(para-Fluoro)-Aad-LRNP-NH<sub>2</sub>

**Molecular Weight:** 892.9

**Peptide Purity:** Peak Area by HPLC ≥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:**

This peptide is a PAR-1 antagonist which inhibited SFLLRNP induced platelet aggregation with an IC<sub>50</sub> of 115 μM. Antagonists specific for the thrombin receptor can be promising compounds for treatment of thrombosis.

**References:**

Fujita, T. et al. *J Biochem* **126**, 174 (1999).

*For Research Use Only*