

## **Product Data Sheet**

Product Name: Protease-Activated Receptor-4, PAR-4 Antagonist, amide

Catalog Number: AS-65493 (1 mg)

Sequence: trans-Cinnamoyl-Tyr-Pro-Gly-Lys-Phe-NH<sub>2</sub>

trans-Cinnamoyl-YPGKF-NH<sub>2</sub>

Molecular Weight: 739.8

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 selective PAR4 antagonist which inhibits thrombin- and PAR-4 agonist induced rat platelet aggregation.

## References:

Hollenberg, MD. and Saifeddine, M. Can J Phys Pharm 79, 439 (2001).

For Research Use Only