



Product Data Sheet

Product Name: Protease-Activated Receptor-1, PAR-1 Agonist, amide

Catalog Number: AS-65479 (1 mg)

Sequence: Ala-parafluoroPhe-Arg-Cha-Cit-Tyr-NH₂

AF(para-Fluoro)R-Cha-Cit-Y-NH₂

Molecular Weight: 883

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 PAR-1 selective agonist displaying a high level of specificity to PAR-1 over PAR-2. The specificity of peptide was evaluated in cell-based calcium signaling assay using HEK293 cells. PAR-1 selective agonists can be used to study PAR-1 activation in vivo.

References:

Kawabata, A. et al. *J Pharmacol Exp Ther* **288**, 358 (1999).

For Research Use Only