

Product Data Sheet

Product Name: Protease-Activated Receptor-2, PAR-2 Agonist, amide

Catalog Number: AS-65490 (1 mg)

Sequence: 2-Furoyl-Leu-Ile-Gly-Arg-Leu-Orn-NH₂

2-Furoyl-LIGRLO-NH₂

Molecular Weight: 778

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:

The peptide 2-furoyl-LIGRLO-NH₂ showed higher potency and receptor selectivity for *in vitro* assays than other PAR-2-activating peptides. 2-Furoyl-LIGRLO-NH2 peptide was equally or more potent than SLIGRL-NH₂ for increasing intracellular calcium in cultured human and rat PAR-2-expressing cells and significantly more potent than SLIGRL-NH₂ in assays of tissue PAR-2 activity.

References:

McGuire, JJ. et al. J Pharmacol Exp Ther 309, 1124 (2004).

For Research Use Only