

Product Data Sheet

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

AS-60217-5 (5 mg) Lot Number: See label on vial Catalog Number:

H-Ser-Leu-Ile-Gly-Lys-Val-NH2 (3-letter code) Sequence:

SLIGKV-NH2 (1-letter code)

Molecular Weight: 615.8

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 µl to 1 mg PAR peptide. This peptide is also soluble in 1%NH₄OH.

Storage: PAR 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 protease-activated receptor activating peptide (PAR-2-AP) corresponds to the PAR2 tethered ligand. Ref: Dulon et al. Am. J. Resp. Cell Mol. Biol. 28, 339 (2003); Kim, M. et al. Cell Biochem. Funct. 20, 339 (2002); Vesey, D. et al. Kidney International 67, 1315 (2005); Hollenberg, M. Can. J. Physiol. Pharmacol. 75, 832 (1997); Nishikawa, H. et al. J. Pharmacol. Exp. Ther. 312, 324 (2005).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Amidated PAR-activating peptides (PAR1 = AP-1 NH₂-TFLLRN-NH₂; PAR2 = AP-2 NH₂-SLIGKV-NH₂) were from Anaspec, Inc. Transfection of DU-145 cells with PAR1 and PAR2 SMARTpool (75 nmol/L) siRNAs (Dharmacon) or Silencer GAPDH siRNA (Ambion) with DharmaFECT1 (0.5 µL) was done with a modification of the double-transfection method. The cells were transfected for 5 h, the growth medium replaced, and the cells were treated with a second transfection after 24 h. The cells were reincubated (24 h) then shifted to serum-free medium prior to the ERK signaling assay-Mize, GJ. et al. Mole. Cancer Res. 6, 1043 (2008).

Published Citations:

Ahn, H. et al. Mol. Pharmacol. 51, 350 (1997). Rauch, BH. et al. Amer. Heart Assoc. 94, 340 (2004). Black, PC. et al. *The Prostate* **67**, 743 (2007). Mize, GJ. et al. Mole. Cancer Res. 6, 1043 (2008). Mize, GJ. et al. Protein Expression & Purification 57, 280 (2008). Wilson, TJ. et al. Cancer Res. 69, 3188 (2009).

Related Products:

Name

	Cat #	Size
Protease-Activated Receptor-1, PAR-1 Agonist, amide (TFLLRNPNDK-NH2)	AS-62936	1 mg
Protease-Activated Receptor-1, PAR-1 Agonist, amide (TFLLRN-NH2)	AS-62937	5 mg
Protease-Activated Receptor-1, PAR-1 Agonist (TFLLRN)	AS-61530	1 mg
Protease-Activated Receptor-3, PAR-3 Agonist, amide (SFNGGP-NH2)	AS-62938	1 mg
Protease-Activated Receptor-3 (1-6), PAR-3 (1-6), human (TFRGAP-NH2)	AS-62657	1 mg
Protease-Activated Receptor-4, PAR-4 Agonist, amide (AYPGKF-NH2)	AS-60218-1 AS-60218-5	1 mg 5 mg
Protease-Activated Receptor-4, PAR-4 Agonist, amide, murine (GYPGKF-NH2)	AS-60778	1 mg

For Research Use Only