



Product Data Sheet

Product Name:	[Pyr1]-Apelin-13	
Catalog Number:	AS-60833 (1 mg)	Lot Number: See label on vial
Sequence:	Pyr-Arg-Pro-Arg-Leu-Ser-His-Lys-Gly-Pro-Met-Pro-Phe-OH (3-letter code) Pyr-RPRLSHKGPMPF-OH (1-letter code)	
Molecular Weight:	1533.8	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: [Pyr1]-Apelin-13 peptide is freely soluble in water.

Storage: [Pyr1]-Apelin-13 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: [125I]-(Pyr1)Apelin-13 binds with high affinity and selectivity to human cardiac tissue. Ref: Katugampola, SD. et al. *Brit. J. Pharma.* **132**, 1255 (2001); Tatemoto, K. et al. *Biochem. Biophys. Res. Commun.* **251**, 471 (1998).

Related Products:

Name	Cat #	Size
[Ala13]-Apelin-13 QRPRLSHKGPMPA	AS-61078	1 mg
Apelin 12 RPRLSHKGPMPF	AS-60836	1 mg
Apelin-16, human, bovine FRRQRPRLSHKGPMPF	AS-63699	1 mg
Apelin-16, human, bovine FRRQRPRLSHKGPMPF	AS-63699	1 mg

For Research Use Only