

Product Data Sheet

Product Name:	(Arg)9	
Catalog Number:	AS-61204 (1 mg)	Lot Number: See label on vial
Sequence:	H-Arg-Arg-Arg-Arg-Arg-Arg-Arg-Arg-OH (3-letter code) RRRRRRRR (1-letter code)	
Molecular Weight:	1423.7	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	
Peptide Reconstitution:	(Arg)9 peptide is freely soluble in w	ater.

Storage: Arg)9 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.

Related Products:

Name (Arg)9 biotin labeled Biotin-LC-RRRRRRRRR-NH2	Cat # AS-64078	Size 1 mg
(Arg)9, FAM-labeled FAM-RRRRRRRR	AS-61207	1 mg
(Arg)9, TAMRA-labeled TAMRA-RRRRRRRR	AS-61208	1 mg

For Research Use Only