



Product Data Sheet

Product Name:	Penetratin-Arg		
Catalog Number:	AS-65464 (1 mg)	Lot Number:	See label on vial
Sequence:	H-Arg-Gln-Ile-Arg-Ile-Trp-Phe-Gln-Asn-Arg-Arg-Met-Arg-Trp-Arg-Arg-OH RQIRIWFQNRMRWRR		
Molecular Weight:	2358.8		
Peptide Purity:	Peak Area by HPLC \geq 95%		
Description:	<p>Penetratin-Arg is a cell-penetrating peptide (CPP) that is derived from the 3rd helix of Drosophila Antennapedia homeodomain protein. Penetratin-Arg is the same as Penetratin except that Lysine residues were substituted with Arginines. It forms an α-helix structure in lipid environment and can permeate cell membrane at low micromolar concentration without significantly affecting membrane. CPP can be conjugated with large molecules and used as a drug delivery vehicle.</p>		
Storage:	<p>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.</p>		
References:	<ol style="list-style-type: none">1. Nasrollahi, SA. et al. <i>Chem Biol Drug Des</i> 08, 639-646 (2012).2. Ye, J. et al. <i>J Am Chem Soc</i> 132 (3), 1-21 (2010).		

For Research Use Only