



Product Data Sheet

Product Name:	Des-Arg9-Bradykinin	
Catalog Number:	AS-65642 (5 mg)	Lot Number: See label on vial
Sequence:	Arg-Pro-Pro-Gly-Phe-Ser-Pro-Phe (3-letter code) RPPGFSPF (1-letter code)	
Molecular Weight:	904.1	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	Des-Arg9-Bradykinin is freely soluble in H ₂ O	
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:	Kininogen cleaved by plasma kallikrein forms bradykinin. Des-Arg9-Bradykinin is a carboxypeptidase metabolite of bradykinin and acts as an agonist for the type 1 (B1) bradykinin receptor.	
References:	Cyr, M. et al. <i>Am J Physiol Heart Circ Physiol</i> . 281, 275 (2001)	

For Research Use Only