



Product Data Sheet

Product Name:	Des-Arg9-[Leu8]-Bradykinin	
Catalog Number:	AS-65643 (5 mg)	Lot Number: See label on vial
Sequence:	Arg-Pro-Pro-Gly-Phe-Ser-Pro-Leu (3-letter code) RPPGFSPL (1-letter code)	
Molecular Weight:	870.1	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	Des-Arg9-[Leu8]-Bradykinin peptide 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:	Des-Arg9-[Leu8]-Bradykinin is a specific antagonist for the type 1 (B1) bradykinin receptor.	
References:	Ann-Maree Duncan. et al <i>J Physiol Regul Integr Comp Physiol.</i> 278(4):R897-904.(2000) Joan M. Annu. Rev. Pharmacol. Toxicol. 31 :129—62(1991)	

For Research Use Only