



Product Data Sheet

Product Name:	D-Arg-[Hyp3, Thi5,8,D-Phe7]-Bradykinin	
Catalog Number:	AS-65645 (5 mg)	Lot Number: See label on vial
Sequence:	D-Arg-Arg-Pro-Hyp-Gly-Thi-Ser-D-Phe-Thi-Arg (3-letter code) rRPOGXsfXR (1-letter code)	
Molecular Weight:	1294.5	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	D-Arg-[Hyp3, Thi5,8,D-Phe7]-Bradykinin peptide is 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:	D-Arg-[Hyp3, Thi5,8,D-Phe7]-Bradykinin is a potent, synthetic antagonist for the type 2(B2) bradykinin receptor.	
References:	J. W. Dear et al. <i>Br. J. Pharmacol.</i> 119(5): 1054–1062. (1996) Ulf H. Lerner et al. <i>Principles of Bone Biology (Third Edition)</i> . (2008)	

For Research Use Only