



## Product Data Sheet

---

Product Name:	Hyp4-Kallidin	
Catalog Number:	AS-65646 (5 mg)	Lot Number: See label on vial
Sequence:	Lys-Arg-Pro-Hyp-Gly-Phe-Ser-Pro-Phe-Arg (3-letter code) KRPOGFSPFR (1-letter code)	
Molecular Weight:	1204.4	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	peptide is freely soluble in H <sub>2</sub> 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 tissue kallikrein forms kallidin (Lys-bradykinin). A portion of kininogen is hydroxylated on the third amino acid of the bradykinin sequence, proline, leading to the formation of hydroxylated kinin peptides. Hydroxylated kinins have similar biological activity as non-hydroxylated kinins. Kallidin is one of the main kinins in mammals. It acts as an agonist for the type 2 (B <sub>2</sub> ) bradykinin receptor and is a potent inflammatory mediator produced during acute and chronic inflammation.	
References:	Altamura, M. et al <i>Regul. Pept.</i> 80, 13 (1999) Siebeck, M. et al. <i>Immunopharmacol.</i> 40, 49 (1998) Wiernas, T. et al. <i>Br. J. Pharmacol.</i> 121, 649 (1997) Mindroiu T, et al. <i>Biochem Biophys Res Commun.</i> 29;152(2):519-26. (1988)	

*For Research Use Only*