



## Product Data Sheet

---

<b>Product Name:</b>	Hyp3-BK(1-8), Hyp3-Des-Arg9-Bradykinin	
<b>Catalog Number:</b>	AS-65644 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	Arg-Pro-Hyp-Gly-Phe-Ser-Pro-Phe (3-letter code) RPOGFSPF (1-letter code)	
<b>Molecular Weight:</b>	920.1	
<b>% Peak Area by HPLC</b>	≥95%	
<b>Appearance</b>	Lyophilized white powder	
<b>Peptide Reconstitution</b>	Hyp3-Des-Arg9-Bradykinin peptide is freely soluble in H <sub>2</sub> O	
<b>Storage:</b>	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.	
<b>Description:</b>	Hyp3-Des-Arg9-Bradykinin acts as an agonist for the type 1 (B1) bradykinin receptor. It is cleaved from kininogens, hydroxylated on the third amino acid, proline, of the bradykinin sequence by plasma kallikrein.	
<b>References:</b>	Ann-Maree Duncan. et al <i>J Physiol Regul Integr Comp Physiol.</i> 278(4): R897-904.(2000)	

*For Research Use Only*