## ANASPEC

## Product Data Sheet

Product Name:	Des-Arg10-Kallidin	
Catalog Number:	AS-65647 (5 mg)	Lot Number: See label on vial
Sequence:	Lys-Arg-Pro-Pro-Gly-Phe-Ser-Pro-Phe (3-letter code) KRPPGFSPF (1-letter code)	
Molecular Weight:	1032.2	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	Des-Arg10-Kallidin peptide is freely soluble in H <sub>2</sub> O	
Storage:	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at $-20^{\circ}$ C or lower.	
Description:	Des-Arg10-Kallidin is a metabolite of kallic type 1 (B1) bradykinin receptor. It is induce myocardial ischemia and inflammation.	
References:	Ibeth Guevara-Lora. et al <i>Acta Biochim</i> . Pol 56(3):515-22. (2009) Ann-Maree Duncan. et al <i>J Physiol Regul Integr Comp Physiol</i> . 278(4):R897-904.(2000)	

For Research Use Only