



Product Data Sheet

Product Name:	Kallikrein-2 inhibitor, cyclic (KLK2b ht2)	
Catalog Number:	AS-65651 (1 mg)	Lot Number: See label on vial
Sequence:	Cyclo-Ala-Arg-Phe-Lys-Val-Trp-Trp-Ala-Asn (3-letter code) Cyclo-ARFKVWWAN (1-letter code)	
Molecular Weight:	1159.4	
% Peak Area by HPLC	≥95%	
Appearance	Lyophilized white powder	
Peptide Reconstitution	KLK2b ht2 inhibitor 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:	<p>KLK2b ht2, a head to tail cyclic peptide, is an inhibitor of human glandular kallikrein enzyme, kallikrein-2 (KLK2). This enzyme belongs to kallikreins, serine protease family, which comprises of 15 members (KLK1-KLK15). KLK2 is increased in tumor tissues and in serum of patients with prostate cancer and is considered as a biomarker for this type of cancer. Inhibition of KLK2 activity might be a promising therapy for prostate cancer. KLK2b ht2 cyclic peptide shows higher stability and biological activity compared with linear peptide inhibitors.</p>	
References:	Miikka Pakkala et al., <i>J Pept Sci.</i> 13(5):348-53. (2007)	

For Research Use Only