

Product Data Sheet

Product Name: Kallikrein-2 inhibitor, cyclic (KLK2b ht4)

AS-65652 (1 mg) Catalog Number: Lot Number: See label on vial

Sequence: Cyclo-Ala-Ala-Arg-Phe-Lys-Val-Trp-Trp-Ala-Ala-Asn (3-letter code)

Cyclo-AARFKVWWAAN (1-letter code)

Molecular Weight: 1301.6

% Peak Area by HPLC ≥95%

Appearance Lyophilized white powder

Peptide Reconstitution KLK2b ht4 inhibitor peptide is freely soluble in H2O

This peptide is shipped at ambient temperature. Upon receipt, store Storage:

lyophilized peptide at -20°C or lower. Reconstituted peptide can be

aliquoted and stored at -20°C or lower.

Description: KLK2b ht4, 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 ht4 cyclic peptide shows higher stability and biological activity

compared with linear peptide inhibitors.

References:

Miikka Pakkala et al., *J Pept Sci.* 13(5):348-53. (2007)

For Research Use Only