

Product Data Sheet

Product Name: Hyp3-KBK(1-8), Hyp4-Des-Arg10-Kallidin

Catalog Number: AS-65648 (5 mg) Lot Number: See label on vial

Sequence: Lys-Arg-Pro-Hyp-Gly-Phe-Ser-Pro-Phe (3-letter code)

KRPOGFSPF (1-letter code)

Molecular Weight: 1048.3

% Peak Area by HPLC ≥95%

Appearance Lyophilized white powder

Peptide Reconstitution Hyp4-Des-Arg10-Kallidin peptide is freely soluble in H₂O

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: Hyp4-Des-Arg10-Kallidin 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 tissue kallikrein.

References: Ann-Maree Duncan. et al J Physiol Regul Integr Comp Physiol. 278(4): R897-904. (2000)

For Research Use Only