

## **Product Data Sheet**

Product Name: Hyp3- Bradykinin

Catalog Number: AS-65641 (5 mg) Lot Number: See label on vial

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

RPOGFSPFR (1-letter code)

Molecular Weight: 1076.3

% Peak Area by HPLC ≥95%

Appearance Lyophilized white powder

Peptide Reconstitution Hyp3- Bradykinin 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°C or lower. Reconstituted peptide can be aliquoted and stored at

-20 °C or lower.

Description: Hyp3- Bradykinin is a natural analog of bradykinin, where the third amino acid,

proline, is modified to (2S,4R)-4-Hydroxyproline. It has similar biological activity to Bradykinin and acts as an agonist for the type 2(B2) bradykinin receptor. Recent research studies reported that bradykinin is also linked to Alzheimer's disease (AD) pathology. The level of plasma bradykinin in AD patients has been shown to be elevated significantly. It interacts with the fibrillar form of  $A\beta$  and co-localizes with

Aβ plaques in the post-mortem human AD brain.

References: Pradeep K. Singh et al. Neurobiology of Disease. 139:104833. (2020)

For Research Use Only