

Product Data Sheet

Product Name:PrP (106-126)Catalog Number:AS-61160 (1 mg)Lot Number: See label on vialSequence:H-Lys-Thr-Asn-Met-Lys-His-Met-Ala-Gly-Ala-Ala-Ala-Ala-Ala-Ala-Ala-Val-Val-
Gly-Gly-Leu-Gly-OH (3-letter code)
KTNMKHMAGAAAAGAVVGGLG (1-letter code)Molecular Weight:1912.3Peptide Purity:>95%Appearance:Lyophilized white powder

Peptide Reconstitution: Initially try to dissolve a small amount of PrP (106-126) peptide in water. If peptide does not dissolve, PrP (106-126) peptide is freely soluble in DMSO.

Storage: PrP (106-126) 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: This peptide increases membrane microviscosity, intracellular Ca2+ concentration and cell migration in circulating leucocytes, and oxygen production in monocytes and neutrophils. It also induces apoptotic cell death and impairment of L-type voltage-sensitive calcium channel activity in the GH3 cell lines. PrP (106–126) stimulates leucocyte migration in a dose-dependent manner. Ref: Florio, T. et al. *J. Neurosci. Res.***54**, 341 (1998).

For Research Use Only

©AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555 Tel: (800)-452-5530 | service@anaspec.com | www.anaspec.com