

Product Data Sheet

Product Name: P0 (180-199), Peripheral Myelin P0 Protein (180-199), mouse

Catalog Number: AS-65470 (1 mg) Lot Number: See label on vial

Sequence: H-Ser-Ser-Lys-Arg-Gly-Arg-Gln-Thr-Pro-Val-Leu-Tyr-

Ala-Met-Leu-Asp-His-Ser-Arg-Ser-OH

SSKRGRQTPVLYAMLDHSRS

Molecular Weight: 2290.6

Peptide Purity: Peak Area by HPLC ≥95%

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:

This peptide is derived from murine peripheral myelin protein P0 amino acid residues 180-199. It is neuritogenic inducing experimental autoimmune neuritis (EAN) in murine.

Related Products:

P0 (106-125), Peripheral Myelin Protein P0 (106-125), mouse, Cat# AS-65469 P2 (53-78), Peripheral Myelin Protein P2 (53-78), bovine, Cat# AS-65472

References:

Zhang, HL. et al. PLoS One 7, e38157 (2012).

For Research Use Only