



Product Data Sheet

Product Name: P2 (53-78), Peripheral Myelin Protein P2 (53-78), bovine

Catalog Number: AS-65472 (1 mg) **Lot Number:** See label on vial

Sequence: H-Thr-Glu-Ser-Pro-Phe-Lys-Asn-Thr-Glu-Ile-Ser-Phe-
Lys-Leu-Gly-Gln-Glu-Phe-Glu-Glu-Thr-Thr-Ala-Asp-Asn-Arg-OH

TESPFKNTEISFKLGQEFEEETTADNR

Molecular Weight: 3019.9

Peptide Purity: Peak Area by HPLC \geq 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 bovine peripheral myelin P2 protein amino acid residues 53-78. It is neurotogenic, inducing experimental autoimmune neuritis (EAN) in Lewis rats.

Related Products:

P0 (180-199), Peripheral Myelin Protein P0 (180-199), mouse, Cat# AS-65470

References:

Hahn, AF. et al. *Acta Neuropathol* **82**, 60 (1991).

For Research Use Only