



Product Data Sheet

Product Name: MBP (69–85), guinea pig
Catalog Number: AS-62081 (1 mg) **Lot Number:** See label on vial
Sequence: H-Tyr-Gly-Ser-Leu-Pro-Gln-Lys-Ser-Gln-Arg-Ser-Gln-Asp-Glu-Asn-Pro-Val-OH (3-letter sequence)
YGSLPQKSQRSQDENPV (1-letter code)

Molecular Weight: 1933.1

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: This is a myelin basic protein encephalitogenic epitope used to induce experimental autoimmune encephalomyelitis (EAE) in rats. Ref: Grabie, N. et al. *Int Immunol* **11**, 899 (1999); Xiao, B. et al. *Clin Exp Immunol*. **125**, 300 (2001).

For Research Use Only