

Product Name: MOG (35-55), mouse, rat

 Catalog Number:
 AS-60130-1 (1 mg)
 Lot Number: See label on vial

 AS-60130-5 (5 mg)
 AS-60130-10 (10 mg)

Sequence: H-Met-Glu-Val-Gly-Trp-Tyr-Arg-Ser-Pro-Phe-Ser-Arg-Val-Val-His-Leu-Tyr-Arg-Asn-Gly-Lys-OH (3-letter code) MEVGWYRSPFSRVVHLYRNGK (1-letter code)

Molecular Weight: 2582.0

% 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 peptide at -20° C or lower. Reconstituted peptide can be aliquoted and stored at -20° C or lower.

Description: This peptide* induced anti-MOG(35-55) autoantibody production in C57BL/6 mice with 100% incident (10/10), providing quantitative support for the suitability of the MOG peptide for *in vivo* anti-MOG(35-55) autoantibody study. Cysteine containing MOG (35-55) peptide induced polyclonal anti-MOG(35-55) antibody production in rabbits with high specificity, high affinity, and high titer indicating its high purity, as well as suitability for peptide-coated ELISA and Western blot assay. *Tested for lot# 61918 & 71369

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

MOG (35-55) peptides (MEVGWYRSPFSRVVHLYRNGK) were purchased from AnaSpec. EAE was induced in WT and different mutant mice 8 to 12 weeks of age by immunization of mice s.c. with 200 µl of an emulsion containing 200 µg MOG (35-55) peptides in 100 µl PBS and 100 µl CFA. CFA was prepared by mixing of incomplete Freund's adjuvant (DIFCO Laboratories) with 8 mg/ml of *Mycobacterium tuberculosis* H37RA (desiccated; DIFCO Laboratories) – <u>Guo, B. et al. *J Clin Invest*</u> **118**, 1680 (2008).

The MOG (35-55) peptide (MEVGWYRSPFSRVVHLYRNGK) was obtained from Anaspec (San Jose, CA), and purified to >98% purity by reverse phase HPLC analysis. Mice were injected subcutaneously (two 100 μ L injections into adjacent areas in one hind limb) with an emulsion of 300 μ g MOG (35–55) dissolved in 100 μ L PBS, mixed with 100 μ L CFA containing 500 μ g of *Mycobacterium tuberculosis* (Difco, Detroit, MI) –<u>Sharp, A. et al. *J Neuroinflamm* 5, 33 (2008).</u>

MOG peptides (35-55), (40-54), or (44-54) which have previously been utilized in the MOG

EAE disease model, and in which CD8 T cell responses have been implicated. The peptides were purchased or synthesized by Anaspec Inc.,San Jose, CA. Briefly, peptides were emulsified (1.5 mg/ml) in complete freund's adjuvant (CFA; Sigma, St Louis, MO) containing 200ng/ml *Mycobacterium tuberculosis* (Difco, Detroit, MI). Emulsion was then deposited

© AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555 Tel: (800) 452-5530 | (510) 791-9560 | service@anaspec.com | www.anaspec.com **Published Citations:**

Polak, P. et al. *Brain* 10.1093/brain/awq362 (2011).Simonini, M. et al. *ASN Neuro* 10.1042/AN20090033 (2010).Li, Q. et al. *Am J Pathol* **175**, 661 (2009).Li, Q. et al. Mol Immunol **46**, 2885 (2009).Guo, B. et al. *J Clin Invest* **118**, 1680 (2008).Sharp, A. et al. *J Neuroinflamm* **5**, 33 (2008).Frausto, R. et al. *J Neuroimmuno* **192**, 124 (2007).

Related Products:

Name MOG (35 - 55), human (MEVGWYRPPFSRVVHLYRNGK)	Cat # AS-63918	Size 1 mg
Recombinant mouse MOG (1 - 125)	AS-55150-100 AS-55150-500 AS-55150-1000	100 ug 500 ug 1000 ug
Recombinant human MOG (1 - 125)	AS-55158-100 AS-55158-500 AS-55158-1000	100 ug 500 ug 1000 ug
Recombinant rat MOG (1 - 125)	AS-55152-100 AS-55152-500 AS-55152-1000	100 ug 500 ug 1000 ug

SensoLyte® Anti-Mouse MOG (1-125) IgG Quantitative ELISA Kit	AS-55156	1 kit
SensoLyte® Anti-Rat MOG (1-125) IgG Quantitative ELISA Kit	AS-55157	1 kit
SensoLyte® Anti-MOG(35-55) IgG Quantitative ELISA Kit (mouse/rat)	AS-54465	1 kit

For Research Use Only