



Product Data Sheet

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) – [Guo, B. et al. J Clin Invest 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 µL injections into adjacent areas in one hind limb) with an emulsion of 300 µg MOG (35–55) dissolved in 100 µL PBS, mixed with 100 µL CFA containing 500 µg of *Mycobacterium tuberculosis* (Difco, Detroit, MI) – [Sharp, A. et al. J Neuroinflamm 5, 33 \(2008\).](#)

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

subcutaneously in the flanks of the hind legs (100µl/deposit) – [Frausto, R. et al. J Neuroimmuno 192, 124 \(2007\).](#)

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