

PRODUCT INFORMATION



MOG (35-55) (mouse, rat) (trifluoroacetate salt)

Item No. 35644

Formal Name:	L-methionyl-L- α -glutamyl-L-valylglycyl-L-tryptophyl-L-tyrosyl-L-arginyl-L-seryl-L-prolyl-L-phenylalanyl-L-seryl-L-arginyl-L-valyl-L-valyl-L-histidyl-L-leucyl-L-tyrosyl-L-arginyl-L-asparaginylglycyl-L-lysine, trifluoroacetate salt	H—Met—Glu—Val—Gly—Trp—Tyr—Arg—Ser—Pro—Phe—Ser—
Synonyms:	MEVGWYRSPFSRVVHLYRNGK, Myelin Oligodendrocyte Glycoprotein (35-55)	Arg—Val—Val—His—Leu—Tyr—Arg—Asn—Gly—Lys—OH
MF:	C ₁₁₈ H ₁₇₇ N ₃₅ O ₂₉ S • XCF ₃ COOH	
FW:	2,582.0	• XCF ₃ COOH
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Myelin oligodendrocyte glycoprotein (MOG) (35-55) (mouse, rat) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the MOG (35-55) (mouse, rat) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

MOG (35-55) is an antigen peptide of MOG, a protein involved in demyelinating CNS disorders.^{1,2} MOG (35-55) in combination with heat-killed *M. tuberculosis* and complete Freund's adjuvant (CFA) induces demyelination of the spinal cord, degeneration of the optic nerve, and visual impairments in mice.¹ Immunization with MOG (35-55) in combination with adjuvants has been used to induce experimental autoimmune encephalomyelitis (EAE) in animal models of multiple sclerosis.^{1,2}

References

1. Guo, X., Harada, C., Namekata, K., *et al.* Delayed onset of experimental autoimmune encephalomyelitis in Olig1 deficient mice. *PLoS One* **5(9)**, e13083 (2010).
2. Zhang, F., Wei, W., Chai, H., *et al.* Aurintricarboxylic acid ameliorates experimental autoimmune encephalomyelitis by blocking chemokine-mediated pathogenic cell migration and infiltration. *J. Immunol.* **190(3)**, 1017-1025 (2013).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/07/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM