

PRODUCT INFORMATION

MOG (92-106) (mouse) (trifluoroacetate salt)

Item No. 45096

Synonym:	Myelin Oligodendrocyte Glycoprotein (92-106)	
Peptide Sequence:	H-DEGGYTCFFRDHSYQ-NH ₂	H-Asp-Glu-Gly-Gly-Tyr-Thr-Cys-Phe-Phe-Arg-
MF:	C ₈₀ H ₁₀₅ N ₂₁ O ₂₇ S • XCF ₃ COOH	
FW:	1,824.9	Asp-His-Ser-Tyr-Gln-NH ₂
Purity:	≥95%	• XCF ₃ COOH
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) (92-106) (mouse) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the MOG (92-106) (mouse) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. MOG (92-106) (mouse) (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Description

(MOG) (92-106) is an encephalitogenic peptide and a peptide fragment of MOG, a protein involved in demyelinating CNS disorders.¹ MOG (92-106) induces relapsing-remitting experimental autoimmune encephalomyelitis (RR-EAE) in SJL/J mice that lack NK1.1⁺ T cells and express a Th2-low phenotype. It induces a primary progressive (PP) form of EAE when used alone but induces the secondary progressive (SP) form when supplemented with *B. pertussis* in A.SW mice that have a Th2-responsive phenotype.

Reference

1. Tsunoda, I., Kuang, L.Q., Theil, D.J., *et al.* Antibody association with a novel model for primary progressive multiple sclerosis: induction of relapsing-remitting and progressive forms of EAE in H2s mouse strains. *Brain Pathol.* **10**(3), 402-418 (2000).

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, 06/04/2026

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