

PRODUCT INFORMATION



MOG Peptide (human, bovine) (acetate)

Item No. 36906

Formal Name:	L-methionyl-L- α -glutamyl-L-valylglycyl-L-tryptophyl-L-tyrosyl-L-arginyl-L-prolyl-L-prolyl-L-phenylalanyl-L-seryl-L-arginyl-L-valyl-L-valyl-L-histidyl-L-leucyl-L-tyrosyl-L-arginyl-L-asparaginylglycyl-L-lysine, acetate	H—Met—Glu—Val—Gly—Trp—Tyr—Arg—Pro—Pro—Phe— Ser—Arg—Val—Val—His—Leu—Tyr—Arg—Asn—Gly— Lys—OH • XCH ₃ CO ₂ H
Synonym:	Myelin Oligodendrocyte Glycoprotein Peptide	
Peptide Sequence:	MEVGWYRPPFSRVVHLYRNGK-OH	
MF:	C ₁₂₀ H ₁₇₉ N ₃₅ O ₂₈ S • XC ₂ H ₄ O ₂	
FW:	2,592.0	
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

MOG Peptide (human, bovine) (acetate) is supplied as a solid. Aqueous solutions of MOG peptide (human, bovine) (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of MOG peptide (human, bovine) (acetate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Myelin oligodendrocyte glycoprotein (MOG) peptide is an endogenous peptide truncation of MOG, which is found on the extracellular surface of myelin sheaths.¹ When coupled to the antigen serotype HLA-DR2, MOG peptide reduces cortical and striatal infarct volumes, decreases sensorimotor impairments in tube- and corner turn tests, increases time spent investigating novel odors, and increases the duration of cage mate-directed calls in female, but not male, mice in a model of ischemic stroke induced by middle cerebral artery occlusion (MCAO).

Reference

1. Pan, J., Palmateer, J., Schallert, T., *et al.* Novel humanized recombinant T cell receptor ligands protect the female brain after experimental stroke. *Transl. Stroke Res.* **5**(5), 577-585 (2014).

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, 08/15/2023

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