

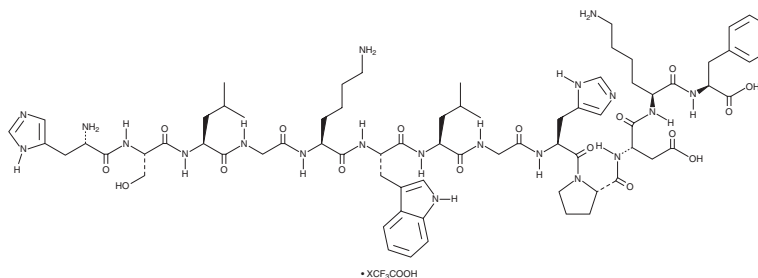
PRODUCT INFORMATION



PLP (139-151) (trifluoroacetate salt)

Item No. 33159

Formal Name: L-histidyl-L-seryl-L-leucylglycyl-L-lysyl-L-tryptophyl-L-leucylglycyl-L-histidyl-L-prolyl-L- α -aspartyl-L-lysyl-L-phenylalanine, trifluoroacetate salt
Synonyms: linPLP, Proteolipid Peptide (139-151)
MF: C₇₂H₁₀₄N₂₀O₁₇ • XCF₃COOH
FW: 1,521.7
Purity: ≥95%
Supplied as: A crystalline 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

PLP (139-151) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the PLP (139-151) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

PLP (139-151) is an encephalitogenic peptide fragment of myelin proteolipid protein (PLP).¹ *In vivo*, PLP (139-151) (150 μ g/animal) induces the formation of inflammatory foci, axonal injury, and demyelination in the spinal cord, as well as motor deficits in mice. It is commonly used to induce experimental autoimmune encephalomyelitis (EAE), a model of multiple sclerosis (MS), in mice.¹⁻³

References

1. Loubopoulos, A., Matsoukas, M.-T., Katsara, M., *et al.* Cyclization of PLP139-151 peptide reduces its encephalitogenic potential in experimental autoimmune encephalomyelitis. *Bioorg. Med. Chem.* **26(9)**, 2221-2228 (2018).
2. Reddy, J., Bettelli, E., Nicholson, L., *et al.* Detection of autoreactive myelin proteolipid protein 139-151-specific T cells by using MHC II (IA^b) tetramers. *J. Immunol.* **170(2)**, 870-877 (2003).
3. Mohindru, M., Kang, B., and Kim, B.S. Functional maturation of proteolipid protein₁₃₉₋₁₅₁-specific Th1 cells in the central nervous system in experimental autoimmune encephalomyelitis. *J. Neuroimmunol.* **155(1-2)**, 127-135 (2004).

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, 11/14/2022

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