

PRODUCT INFORMATION



[pGlu⁴]-Myelin Basic Protein Peptide (trifluoroacetate salt)

Item No. 40149

Formal Name: 5-oxo-L-prolyl-L-lysyl-L-arginyl-L-prolyl-L-seryl-L-glutamyl-L-arginyl-L-seryl-L-lysyl-L-tyrosyl-L-leucine, trifluoroacetate salt

Synonym: [pGlu⁴]-MBP Peptide

Peptide Sequence: pEKRPQRSKYL-OH pGlu—Lys—Arg—Pro—Ser—Gln—Arg—Ser—Lys—Tyr—Leu—OH

MF: C₆₀H₁₀₀N₂₀O₁₇ • XCF₃COOH • XCF₃COOH

FW: 1,373.6

Purity: ≥98%

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

[pGlu⁴]-Myelin basic protein (MBP) peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the [pGlu⁴]-MBP peptide (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

[pGlu⁴]-MBP peptide is a substrate for PKC.¹ It is selectively phosphorylated by PKC over PKA and phosphorylase kinase.

Reference

1. Toomik, R. and Ek, P. A potent and highly selective peptide substrate for protein kinase C assay. *Biochem. J.* 322(Pt 2), 455-460 (1997).

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, 02/22/2024

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