

PRODUCT INFORMATION



[Ala¹⁰⁷]-Myelin Basic Protein (104-118) (acetate)

Item No. 43009

Formal Name: (S)-2-((5S,11S,14S,17S,20S,23S,26S,29S,32S,38S)-32-((1H-indol-3-yl)methyl)-26-benzyl-23-(3-((diaminomethylene)amino)propyl)-38-(hydroxyamino)-14,20,29-tris(hydroxymethyl)-11,17-diisobutyl-5-methyl-4,7,10,13,16,19,24,25,28,31,34,37-dodecaoxo-3,6,9,12,15,18,21,22,27,30,33,36-dodecaazanona triacontanamido)-6-amino-N-(2-oxoethyl)hexanamide, acetate
Synonyms: [Ala¹⁰⁷]-MBP (104-118)

Peptide Sequence: GKGAGLSLSRFSWGA-OH

MF: C₆₇H₁₀₄N₂₀O₁₉ • XC₂H₄O₂

FW: 1,493.7

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H — Gly — Lys — Gly — Ala — Gly — Leu — Ser — Leu — Ser — Arg —

Phe — Ser — Trp — Gly — Ala — OH

• XCH₃CO₂H

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

Laboratory Procedures

[Ala¹⁰⁷]-Myelin basic protein (MBP) (104-118) (acetate) is supplied as a solid. A stock solution may be made by dissolving the [Ala¹⁰⁷]-MBP (104-118) (acetate) in the solvent of choice, which should be purged with an inert gas. [Ala¹⁰⁷]-MBP (104-118) (acetate) is soluble (≥10 mg/ml) in DMSO.

Description

[Ala¹⁰⁷]-Myelin basic protein (MBP) (104-118) is a peptide inhibitor of PKC.¹ It is selective for PKC over myosin light-chain kinase (MLCK) and PKA at 100 μM. [Ala¹⁰⁷]-MBP (104-118) inhibits the phosphorylation of histone H1, MBP, [Ala¹⁰⁵]-MBP (104-118), MBP (104-118), and MBP (104-123) by PKC with IC₅₀ values of 126, 145, 46, 72, and 63 μM, respectively.

Reference

1. Su, H.D., Kemp, B.E., Turner, R.S., et al. Synthetic myelin basic protein peptide analogs are specific inhibitors of phospholipid/calcium-dependent protein kinase (protein kinase C). *Biochem. Biophys. Res. Commun.* **134**(1), 78-84 (1986).

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, 04/02/2025

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