

PRODUCT INFORMATION



TB-500 (acetate)

Item No. 45081

Formal Name: N-acetyl-L-leucyl-L-lysyl-L-lysyl-L-threonyl-L- α -glutamyl-L-threonyl-L-glutamine, acetate

Synonym: Ac-Leu-Lys-Lys-Thr-Glu-Thr-Gln-OH

Peptide Sequence: Ac-LKKTETQ-OH

MF: C₃₈H₆₈N₁₀O₁₄ • XC₂H₄O₂

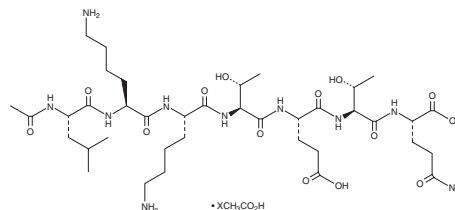
FW: 889.0

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

TB-500 (acetate) is supplied as a solid. Aqueous solutions of TB-500 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. TB-500 (acetate) is soluble (\geq 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

TB-500 (acetate) is a peptide fragment corresponding to amino acid residues 17-23 of thymosin β 4 (T β 4; Item No. 28605), an actin-sequestering peptide.¹ TB-500 inhibits proliferation and migration of PDGF-BB-stimulated hepatic stellate cells at a concentration of 10 ng/ml. Topical application of TB-500 (15 μ g/30 μ l) reduces wound closure and healing times of skin wounds in mice.²

References

1. Shah, R., Reyes-Gordillo, K., and Rojkind, M. Thymosin β 4 inhibits PDGF-BB induced activation, proliferation, and migration of human hepatic stellate cells via its actin-binding domain. *Expert Opin. Biol. Ther.* **18(sup1)**, 177-184 (2018).
2. Huang, C.M., Wang, C.C., Barnes, S., et al. In vivo detection of secreted proteins from wounded skin using capillary ultrafiltration probes and mass spectrometric proteomics. *Proteomics* **6(21)**, 5805-5814 (2006).

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, 03/10/2026

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