

PRODUCT INFORMATION



Melanostatin DM (acetate)

Item No. 45765

Formal Name: L-histidyl-D-arginyl-L-alanyl-L-tryptophyl-D-phenylalanyl-L-lysineamide, acetate

Synonym: His-D-Arg-Ala-Trp-D-Phe-Lys-NH₂

Peptide Sequence: HrAWfK-NH₂

MF: C₄₁H₅₈N₁₄O₆ • XC₂H₄O₂

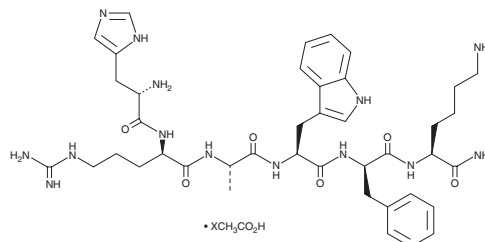
FW: 843.0

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

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

Description

Melanostatin DM is a peptide α -melanocyte-stimulating hormone (α -MSH) antagonist ($pA_2 = 7$ in *U. ornatus* melanocytes).¹ It inhibits skin darkening induced by α -MSH (Item No. 29923) or [Nle⁴,D-Phe⁷]- α -MSH (Item Nos. 41548 | 24475) in white background-adapted *U. ornatus*.

Reference

1. De L. Castrucci, A.M., Sherbrooke, W.C., Sawyer, T.K., *et al.* Discovery of an α -melanotropin antagonist effective in vivo. *Peptides* **15**(4), 627-632 (1994).

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, 07/29/2026

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