

PRODUCT INFORMATION



Myr5A Peptide

Item No. 40257

Overview and Properties

Contents:	This vial contains 1 mg of lyophilized peptide.
Synonym:	Myristic Acid 5A Peptide
Peptide Sequence:	Myr-DWLKAFYDKVAEKLKEAFPDWAKAAYDKAAEKAKEAA (Myr=myristic acid)
Molecular Weight:	4,428.0
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥1 year

Description

Myr5A peptide is an acylated peptide composed of the apolipoprotein A1 (ApoA1) mimetic peptide 5A peptide conjugated to the saturated fatty acid myristic acid (Item No. 13351).¹ It self-assembles into lipid nanostructures and has been used to encapsulate the anthracyclines doxorubicin (Item No. 15007) and valrubicin (Item No. 22964) for *in vitro* compound-release studies.

Reference

1. Raut, S., Garud, A., Nagarajan, B., *et al.* Probing the assembly of HDL mimetic, drug carrying nanoparticles using intrinsic fluorescence. *J. Pharmacol. Exp. Ther.* **373**(1), 113-121 (2020).

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, 12/21/2023

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