

PRODUCT INFORMATION



Pep2 Peptide

Item No. 40258

Contents: This vial contains 1 mg of lyophilized peptide.

Peptide Sequence: RRGTIAFDNWVDTGTRVYDC

Molecular Weight: 2.3456 kDa

Storage: -20°C

Stability: ≥1 year

Description

Pep2 peptide is a peptide ligand for the transmembrane glycoprotein CD36, also known as scavenger receptor B2.¹ It selectively binds to CD36 over human serum albumin (HSA), IgG, and CD44 in a cell-free assay when used at a concentration of 1 mM. Pep2 peptide-functionalized micelles encapsulating the anticancer agent doxorubicin (Item No. 15007) increase cellular delivery of doxorubicin and enhance cellular toxicity in CD36-overexpressing HepG2 hepatocellular carcinoma cells compared with micelles lacking Pep2 peptide.

Reference

1. Zheng, Y., Fang, X., Yang, Y., *et al.* Peptide-directed delivery of drug-loaded nanocarriers targeting CD36 overexpressing cells. *Colloid. Surface. A* **610**, 12590 (2021).

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, 12/13/2023

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