

PRODUCT INFORMATION



M2 Peptide

Item No. 40263

Overview and Properties

Contents: This vial contains 1 mg of a lyophilized peptide trifluoroacetate salt.
Synonyms: M2pep, TAM-targeting Peptide
Peptide Sequence: YWWKVGWPDQEYC-NH₂
Molecular Weight: 1,757.7
Form: Solid
Storage: -20°C (as supplied)
Stability: ≥1 year

Description

M2 peptide is an M2 tumor-associated macrophage-targeting peptide.^{1,2} It has been used in the generation of anticancer polymeric micelles encapsulating the STAT6 inhibitor AS1517499 (Item No. 29071) for delivery to M2 tumor-associated macrophages (TAMs) *in vitro* and *in vivo*. Cayman's M2 Peptide is intended for use in peptide conjugation to lipid-based drug delivery platforms containing maleimide-activated lipids such as DSPE-PEG(2000) maleimide for TAM-targeting applications.

References

1. Xiao, H., Guo, Y., Li, B., *et al.* M2-like tumor-associated macrophage-targeted codelivery of STAT6 inhibitor and IKK β siRNA induces M2-to-M1 repolarization for cancer immunotherapy with low immune side effects. *ACS Cent. Sci.* **6(7)**, 1208-1222 (2020).
2. Huang, T., Zhang, Q., Yi, J., *et al.* PEG-sheddable nanodrug remodels tumor microenvironment to promote effector T cell infiltration and revise their exhaustion for breast cancer immunotherapy. *Small* **19(38)**, e2301749 (2023).

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, 01/15/2025

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