

PRODUCT INFORMATION



TET 830-modified T Helper Cell Epitope from Tetanus Toxin (trifluoroacetate salt)

Item No. 43815

Formal Name: L-alanyl-L-glutamyl-L-tyrosyl-L-isoleucyl-L-lysyl-L-alanyl-L-asparagyl-L-seryl-L-lysyl-L-phenylalanyl-L-isoleucylglycyl-L-isoleucyl-L-threonyl-L- α -glutamyl-L-leucine, trifluoroacetate salt

H—Ala—Gln—Tyr—Ile—Lys—Ala—Asn—Ser—Lys—Phe—
Ile—Gly—Ile—Thr—Glu—Leu—OH

Peptide Sequence: AQYIKANSKFIGITEL-OH

MF: $C_{83}H_{134}N_{20}O_{24} \cdot XCF_3COOH$

FW: 1,796.1

$\cdot XCF_3COOH$

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

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

Laboratory Procedures

TET 830-modified T helper cell epitope from tetanus toxin (trifluoroacetate salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

TET 830-modified T helper cell epitope from tetanus toxin is an immunogenic T helper (Th) cell epitope derived from the tetanus toxin helper peptide.¹ It increases serum levels of IgG, IgA, IL-4, and IFN- γ in combination with self-amplifying RNA lipid nanoparticles (saLNPs) in a mouse model of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) infection when administered as an adjuvant.

Reference

1. Keikha, R. and Jebali, A. The effect of a peptide adjuvant (YLQPRTFLL) on immune response against SARS-CoV-2 in Balb/C mice when vaccinated with saRNA LNPs. *Appl. Biosci.* (2025).

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, 11/05/2025

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