

PRODUCT INFORMATION



HTLV-1 Tax (11-19) (trifluoroacetate salt)

Item No. 44035

Formal Name: L-leucyl-L-leucyl-L-phenylalanylglycyl-L-tyrosyl-L-prolyl-L-valyl-L-tyrosyl-L-valine, trifluoroacetate salt

Synonyms: HTLV-1 Tax-1 (11-19), Human T Cell Leukemia Virus Type 1 Tax (11-19), Human T Lymphotropic Virus Type 1 Tax (11-19), Leu-Leu-Phe-Gly-Tyr-Pro-Val-Tyr-Val

Peptide Sequence: LLFGYPVYV-OH

MF: C₅₆H₇₉N₉O₁₂ • XCF₃COOH

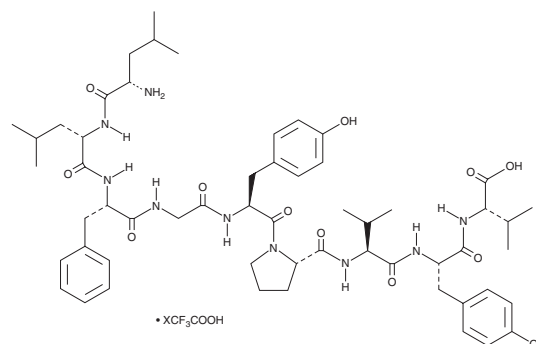
FW: 1,070.3

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

HTLV-1 Tax (11-19) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the HTLV-1 Tax (11-19) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. HTLV-1 Tax (11-19) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Human T cell leukemia virus type 1 (HTLV-1) Tax (11-19) is an immunogenic T cell epitope derived from the HTLV-1 oncoprotein Tax.¹⁻³ It binds to HLA-A*0201 in a cell-free assay when used at a concentration of 1 nM.¹ Co-culture of HTLV-1 Tax (11-19)-primed isolated human HLA-A2-positive mature dendritic cells with naïve autologous peripheral blood lymphocytes induces IFN-γ production in the CD8⁺ T cell population.² HTLV-1 Tax (11-19) has been used to evaluate the role of dendritic cells in the cytotoxic CD8⁺ T cell response in transgenic mice.³

References

1. Pereira-Santos, T.A., da Rocha, A.S., Lopes-Ribeiro, Á., *et al.* Diversity of HLA-A2-restricted and immunodominant epitope repertoire of human T-lymphotropic virus type 1 (HTLV-1) Tax protein: Novel insights among N-terminal, central and C-terminal regions. *Biomolecules* **13**(3), 545 (2023).
2. Manuel, S.L., Schell, T.D., Acheampong, E., *et al.* Presentation of human T cell leukemia virus type 1 (HTLV-1) Tax protein by dendritic cells: The underlying mechanism of HTLV-1-associated neuroinflammatory disease. *J. Leukoc. Biol.* **86**(5), 1205-1216 (2009).
3. Sagar, D., Masih, S., Schell, T., *et al.* In vivo immunogenicity of Tax(11-19) epitope in HLA-A2/DTR transgenic mice: Implication for dendritic cell-based anti-HTLV-1 vaccine. *Vaccine* **32**(26), 3274-3284 (2014).

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, 10/14/2025

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