

PRODUCT INFORMATION



PRAME (425-433) (human) (trifluoroacetate salt)

Item No. 44403

Formal Name: (S)-2-((2S,5S,8S,14S)-2-((1H-imidazol-4-yl)methyl)-8-((S)-sec-butyl)-14-(hydroxyamino)-5-isobutyl-16-methyl-4,7,10,13-tetraoxo-3,6,9,12-tetraazaheptadecanamido)-N1-((S)-1-(((S)-1-(((S)-1-hydroxy-3-oxopropan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)pentanediamide, trifluoroacetate salt H-Ser-Leu-Leu-Gln-His-Leu-Ile-Gly-Leu-OH

Synonyms: Cancer Testis Antigen 130 (425-433), CT130 (425-433), MAPE (425-433), Melanoma Antigen Preferentially Expressed in Tumors (425-433), OIP4 (425-433), Opa-interacting Protein 4 (425-433), Preferentially Expressed Antigen in Melanoma (425-433), Ser-Leu-Leu-Gln-His-Leu-Ile-Gly-Leu • XCF₃COOH

Peptide Sequence: SLLQHLIGL-OH

MF: C₄₆H₈₀N₁₂O₁₂ • XCF₃COOH

FW: 993.2

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

PRAME (425-433) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PRAME (425-433) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. PRAME (425-433) (human) (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO. PRAME (425-433) (human) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Preferentially expressed antigen in melanoma (PRAME) (425-433) is an immunogenic T cell epitope.¹ It binds to HLA-A*0201-positive JY B lymphoblastoid cells with an IC₅₀ value of 3.7 μM. Co-culture of PRAME (425-433)-primed isolated human HLA-A*0201-positive antigen-presenting cells with naïve autologous CD8⁺ T cells to generate PRAME (425-433)-specific cytotoxic T lymphocytes induces lysis of MCF-7 breast cancer cells.

Reference

1. Kessler, J.H., Beekman, N.J., Bres-Vloemans, S.A., *et al.* Efficient identification of novel HLA-A(*)0201-presented cytotoxic T lymphocyte epitopes in the widely expressed tumor antigen PRAME by proteasome-mediated digestion analysis. *J. Exp. Med.* **193**(1), 73-88 (2001).

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, 10/21/2025

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