

Item No. 24578 * Batch No. 0664990

Purity Specification: $\geq 95\%$ Molecular formula: C₂₁₆H₃₇₁N₇₅O₅₉S₆ • XCF₃COOH

CAS number: Formula weight: 5155.12

Expiration date: 04NOV2029

Overview

Tests	Results
Mass Spec	[M+6H] ⁶⁺ : 860.2
HPLC	Purity: 95.0 %
Peptide content	72.50

Reviewed and approved by: Justine Albertson

S. Subburay

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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.

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, 10/12/2018

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