

# PRODUCT INFORMATION



## FITC-Stearyl-R8

Item No. 43816

**Formal Name:** N-((6S,9S,12S,15S,18S,21S,24S,27S,33S)-1-amino-33-carbamoyl-39-((3',6'-dihydroxy-3-oxo-3H-spiro[isobenzofuran-1,9'-xanthen]-5-yl)amino)-9,12,15,18,21,24,27-heptakis(3-guanidinopropyl)-1-imino-7,10,13,16,19,22,25,28,31-nonaoxo-39-thioxo-2,8,11,14,17,20,23,26,29,32,38-undecaazanotriacontan-6-yl)stearamide

**Synonyms:** FITC-Octadecyl-R8, FITC-Stearyl-octaarginine,

**Peptide Sequence:** CH<sub>3</sub>(CH<sub>2</sub>)<sub>16</sub>-CONH-RRRRRRRRGK(FITC)-NH<sub>2</sub>

**MF:** C<sub>95</sub>H<sub>159</sub>N<sub>37</sub>O<sub>16</sub>S

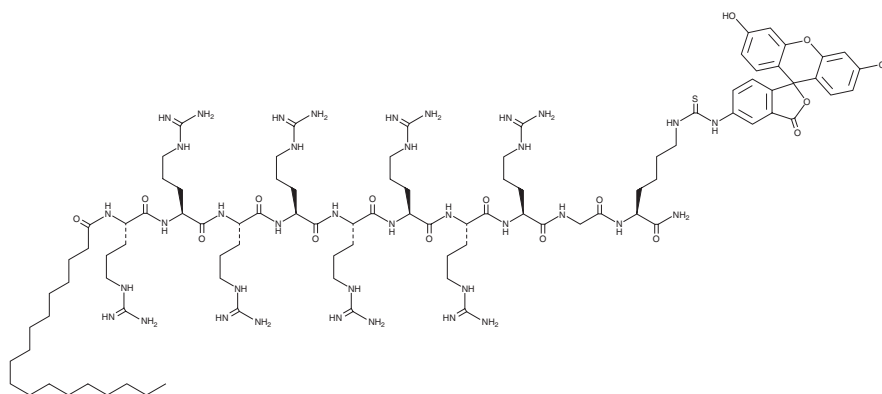
**FW:** 2,107.6

**Purity:** ≥90%

**Supplied as:** A semi-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

FITC-Stearyl-R8 is supplied as a semi-solid. A stock solution may be made by dissolving the FITC-stearyl-R8 in the solvent of choice, which should be purged with an inert gas. FITC-Stearyl-R8 is soluble in a 2:1 solution acetonitrile:water.

### Description

FITC-Stearyl-R8 is a fluorescently labeled form of the stearic acid-conjugated cell-penetrating peptide stearyl-R8. FITC displays excitation/emission maxima of 485/535 nm, respectively. FITC-Stearyl-R8 has been used as a positive control to monitor cellular uptake of cell-penetrating peptidomimetics.<sup>1</sup>

### Reference

1. Madhu, C., Voshavar, C., Rajasekhar, K., et al. Cyclic dipeptide based cell-penetrating peptidomimetics for effective DNA delivery. *Org. Biomol. Chem.* **15**(15), 3170-3174 (2017).

#### 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, 12/05/2025

#### 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