

# PRODUCT INFORMATION



## RTDLDSLRTYTL (trifluoroacetate salt)

Item No. 37455

**Formal Name:** L-arginyl-L-threonyl-L- $\alpha$ -aspartyl-L-leucyl-L- $\alpha$ -aspartyl-L-seryl-L-leucyl-L-arginyl-L-threonyl-L-tyrosyl-L-threonyl-L-leucine, trifluoroacetate salt

**Peptide Sequence:** RTDLDSLRTYTL-OH

**MF:** C<sub>62</sub>H<sub>104</sub>N<sub>18</sub>O<sub>22</sub> • XCF<sub>3</sub>COOH

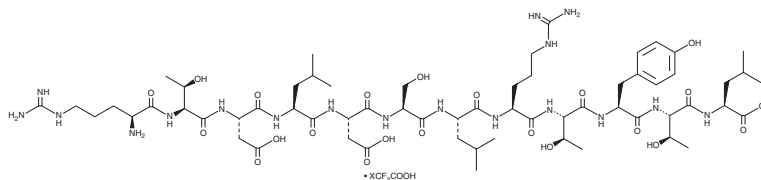
**FW:** 1,453.6

**Purity:**  $\geq$ 98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

RTDLDSLRTYTL (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the RTDLDSLRTYTL (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

RTDLDSLRTYTL is an  $\alpha$ V $\beta$ 6 integrin ligand.<sup>1</sup> It selectively inhibits the binding of fibronectin to  $\alpha$ V $\beta$ 6 integrin over vitronectin to  $\alpha$ V $\beta$ 3 or  $\alpha$ V $\beta$ 5 and fibrinogen to  $\alpha$ II $\beta$ 3 (IC<sub>50</sub>s = 0.02, >50, >50, and >50  $\mu$ M, respectively). RTDLDSLRTYTL also inhibits the binding of HT-29 colon cancer cells to fibronectin in a concentration-dependent manner.

### Reference

1. Kraft, S., Diefenbach, B., Mehta, R., *et al.* Definition of an unexpected ligand recognition motif for  $\alpha$ V $\beta$ 6 integrin. *J. Biol. Chem.* **274**(4), 1979-1985 (1999).

#### 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, 10/16/2023

#### 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