

# PRODUCT INFORMATION



## RAD17-derived Peptide (trifluoroacetate salt)

Item No. 37512

**Formal Name:** L-alanyl-L-seryl-L- $\alpha$ -glutamyl-L-leucyl-L-prolyl-L-alanyl-L-seryl-L-glutamyl-L-prolyl-L-glutamyl-L-prolyl-L-phenylalanyl-L-seryl-L-alanyl-L-lysyl-L-lysyl-L-Lysine, trifluoroacetate salt  
H—Ala—Ser—Glu—Leu—Pro—Ala—Ser—Gln—Pro—Gln—Pro—Phe—Ser—Ala—Lys—Lys—Lys—OH

**Peptide Sequence:** ASELPASQPQPFSAKKK-OH

**MF:**  $C_{81}H_{132}N_{22}O_{25} \cdot XCF_3COOH$

**FW:** 1,814.1 •  $XCF_3COOH$

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

**Supplied as:** A solid

**Storage:**  $-20^\circ C$

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

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

### Laboratory Procedures

RAD17-derived peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the RAD17-derived peptide (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

RAD17-derived peptide is a peptide substrate for ataxia-telangiectasia and RAD3-related protein/kinase (ATR).<sup>1</sup> It has been used to identify inhibitors of ATR.

### Reference

1. Charrier, J.-D., Durrant, S.J., Golec, J.M.C., *et al.* Discovery of potent and selective inhibitors of ataxia telangiectasia mutated and Rad3 related (ATR) protein kinase as potential anticancer agents. *J. Med. Chem.* **54**(7), 2320-2330 (2011).

#### 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, 10/12/2023

#### 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