

# PRODUCT INFORMATION



## P9R (trifluoroacetate salt)

Item No. 37463

**Formal Name:** L-asparaginylglycyl-L-alanyl-L-isoleucyl-L-cysteinyl-L-tryptophylglycyl-L-prolyl-L-cysteinyl-L-prolyl-L-threonyl-L-alanyl-L-phenylalanyl-L-arginyl-L-glutamyl-L-isoleucylglycyl-L-asparaginy-L-cysteinylglycyl-L-arginyl-L-phenylalanyl-L-arginyl-L-valyl-L-arginyl-L-cysteinyl-L-cysteinyl-L-arginyl-L-isoleucyl-L-arginine, trifluoroacetate

**Peptide Sequence:** NGAICWGPCPTAFRQIGNCGRFRVRCRIR-OH

**MF:**  $C_{144}H_{232}N_{52}O_{35}S_5 \cdot XCF_3COOH$

**FW:** 3,412.1

**Purity:**  $\geq 95\%$

**Supplied as:** A solid

**Storage:**  $-20^\circ\text{C}$

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

H—Asn—Gly—Ala—Ile—Cys—Trp—Gly—Pro—Cys—Pro—

Thr—Ala—Phe—Arg—Gln—Ile—Gly—Asn—Cys—Gly—

Arg—Phe—Arg—Val—Arg—Cys—Cys—Arg—Ile—Arg—OH

•  $XCF_3COOH$

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

### Laboratory Procedures

P9R (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the P9R (trifluoroacetate salt) in the solvent of choice. P9R (trifluoroacetate salt) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

P9R is a peptide with antiviral activity.<sup>1</sup> It inhibits viral replication of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2), Middle East respiratory syndrome coronavirus (MERS-CoV), SARS-CoV, influenza A subtypes H1N1 and H7N9, and rhinovirus in plaque reduction assays ( $IC_{50}$ s = 0.9, 2.2, 4.2, 0.6, 0.9, and 5.7  $\mu\text{g/ml}$ , respectively). *In vivo*, P9R (50  $\mu\text{g/animal}$  three times over two days) increases survival and reduces H1N1 particle-forming units (PFUs) in the lungs to a similar extent as the influenza neuraminidase inhibitor zanamivir (Item No. 15123) in H1N1-infected mice.

### Reference

1. Zhao, H., To, K.K.W., Sze, K.-H., *et al.* A broad-spectrum virus- and host-targeting peptide against respiratory viruses including influenza virus and SARS-CoV-2. *Nat. Commun.* **11**(1), 4252 (2020).

#### 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, 08/01/2024

#### 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](http://WWW.CAYMANCHEM.COM)