

# PRODUCT INFORMATION



## F9170 (trifluoroacetate salt)

Item No. 36962

**Formal Name:** l-cyl-L-tryptophyl-L- $\alpha$ -glutamyl-L-alanyl-L-leucyl-L-lysyl-L-tyrosyl-L-leucyl-L-tryptophyl-L-asparaginy-L-leucyl-L-leucyl-L-glutamyl-L-tyrosyl-L-tryptophan, trifluoroacetate salt

**Peptide Sequence:** GWEAKLKYLWNLLQYW-NH<sub>2</sub>

**MF:** C<sub>100</sub>H<sub>135</sub>N<sub>21</sub>O<sub>22</sub> • XCF<sub>3</sub>COOH

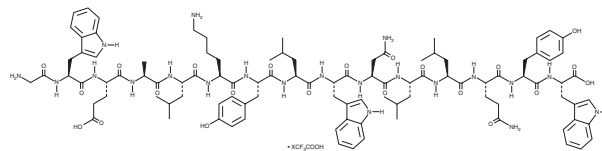
**FW:** 1,983.3

**Purity:** ≥90%

**Supplied as:** A 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

F9170 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the F9170 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. F9170 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of F9170 (trifluoroacetate salt) in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml and DMSO.

### Description

F9170 is an antiviral peptide corresponding to amino acids 789-803 of the HIV-1 envelope glycoprotein.<sup>1</sup> It binds to the LLP1 domain within the HIV-1 envelope protein cytoplasmic tail and inhibits HIV-1 IIIB infection of MT-2 cells ( $EC_{50} = 0.19 \mu\text{M}$ ). F9170 (3 mg/kg) reduces plasma viral load in a rhesus macaque model of simian HIV (SHIV) infection.

### Reference

1. Wang, Q., Su, S., Xue, J., *et al.* An amphipathic peptide targeting the gp41 cytoplasmic tail kills HIV-1 virions and infected cells. *Sci. Transl. Med.* **12(546)**, eaaz2254 (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, 12/18/2023

#### 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