

PRODUCT INFORMATION



WRW4-OH (trifluoroacetate salt)

Item No. 29913

Formal Name: L-tryptophyl-L-arginyl-L-tryptophyl-L-tryptophyl-L-tryptophyl-L-tryptophan, trifluoroacetate salt

Peptide Sequence: WRWWWW-OH

MF: $C_{61}H_{64}N_{14}O_7 \cdot XCF_3COOH$

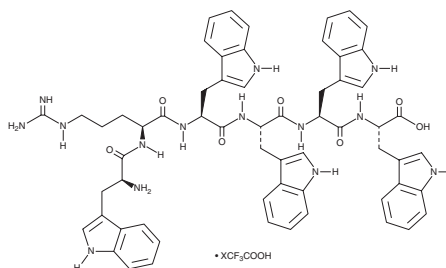
FW: 1,105.3

Purity: $\geq 95\%$

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

WRW4-OH (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the WRW4-OH (trifluoroacetate salt) in water. The solubility of WRW4-OH (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

WRW4-OH is a peptide and the free acid form of the formyl peptide receptor 2 (FPR2) antagonist WRW4-NH₂. It has been used as a model peptide in the study of peptide-carbon nanotube interactions.¹

Reference

1. Davis, T.A. and Holland, L.A. Peptide probe for multiwalled carbon nanotubes: Electrophoretic assessment of the binding interface and evaluation of surface functionalization. *ACS Appl. Mater. Interfaces* **10**(13), 11311-11318 (2018).

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, 08/07/2025

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