

# PRODUCT INFORMATION



## PTD2 (trifluoroacetate salt)

Item No. 38842

**Formal Name:** (S)-2-(2-((5S,8S)-8-(2-acetamidoacetamido)-9-methyl-6,7-dioxodecan-5-yl)hydrazineyl)-N-((2S,3S)-1-amino-3-methyl-1-oxopentan-2-yl)-5-guanidinopentanamide, trifluoroacetate salt  
**Peptide Sequence:** Ac-GV[Nle]RI-NH<sub>2</sub> Ac-Gly-Val-Nle-Arg-Ile-NH<sub>2</sub>  
**MF:** C<sub>27</sub>H<sub>51</sub>N<sub>9</sub>O<sub>6</sub> • XCF<sub>3</sub>COOH • XCF<sub>3</sub>COOH  
**FW:** 597.8  
**Purity:** ≥98%  
**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

PTD2 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PTD2 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. PTD2 (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PTD2 (trifluoroacetate salt) in these solvents is approximately 10 and 2 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of PTD2 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of PTD2 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PTD2 is a peptide inhibitor of the histone methyltransferase Wolf-Hirschhorn syndrome candidate 1 (WHSC1; IC<sub>50</sub> = 22 μM) that was derived from the histone 4 sequence.<sup>1</sup>

### Reference

1. Morrison, M.J., Boriack-Sjodin, P.A., Swinger, K.K., *et al.* Identification of a peptide inhibitor for the histone methyltransferase WHSC1. *PLoS One* **13**(5), e0197082 (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, 08/23/2023

#### 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