

# PRODUCT INFORMATION



## VTP-27999 (trifluoroacetate salt)

Item No. 27689

CAS Registry No.: 1013937-63-7

Formal Name: N-[2-[(R)-(3-chlorophenyl)[(3R)-1-[[[(2S)-2-(methylamino)-3-[(3R)-tetrahydro-2H-pyran-3-yl]propyl]amino]carbonyl]-3-piperidinyl]methoxy]ethyl]-carbamic acid, methyl ester, trifluoroacetate salt

MF:  $C_{26}H_{41}ClN_4O_5 \cdot CF_3COOH$

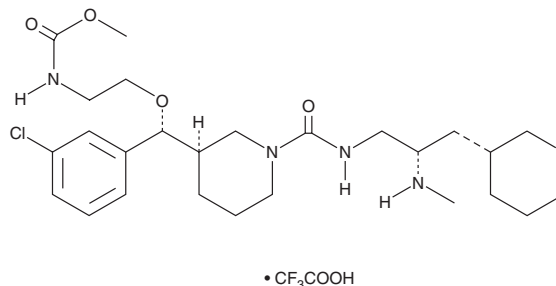
FW: 639.1

Purity:  $\geq 98\%$

Supplied as: A crystalline 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

VTP-27999 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the VTP-27999 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. VTP-27999 (trifluoroacetate salt) is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of VTP-27999 (trifluoroacetate salt) in these solvents is approximately 10 and 1 mg/ml, respectively.

VTP-27999 (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VTP-27999 (trifluoroacetate salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. VTP-27999 (trifluoroacetate salt) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

VTP-27999 is a renin inhibitor ( $IC_{50} = 0.47$  nM for the recombinant human enzyme).<sup>1</sup> It is selective for renin over the aspartyl proteases  $\beta$ -secretase, cathepsin D, and cathepsin E at 10  $\mu M$  and is greater than 1,000-fold selective for renin over a panel of more than 250 receptors, ion channels, and enzymes. VTP-27999 (10 mg/kg, p.o.) reduces mean arterial pressure in rats expressing human renin and angiotensinogen, a double-transgenic model for angiotensin II-dependent hypertension.

### Reference

1. Jia, L., Simpson, R.D., Yuan, J., *et al.* Discovery of VTP-27999, an alkyl amine renin inhibitor with potential for clinical utility. *ACS Med. Chem. Lett.* **2**(10), 747-751 (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, 09/29/2022

#### 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