

# PRODUCT INFORMATION



## Vapreotide (trifluoroacetate salt)

Item No. 24558

**Formal Name:** D-phenylalanyl-L-cysteinyl-L-tyrosyl-D-tryptophyl-L-lysyl-L-valyl-L-cysteinyl-L-tryptophanamide, cyclic (2→7)-disulfide, trifluoroacetate salt

**Synonyms:** BMY 41606, Octastatin, RC 160

**MF:** C<sub>57</sub>H<sub>70</sub>N<sub>12</sub>O<sub>9</sub>S<sub>2</sub> • XCF<sub>3</sub>COOH

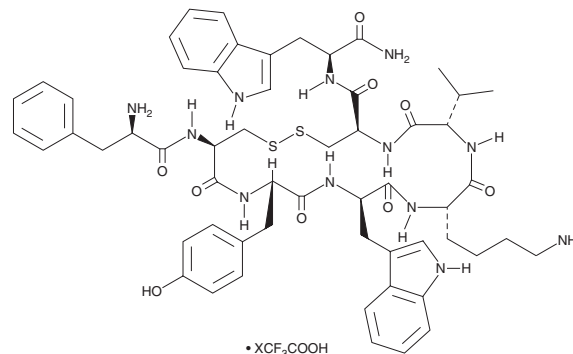
**FW:** 1,131.4

**Purity:** ≥95%

**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

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

### Description

Vapreotide is a peptide neurokinin-1 receptor (NK<sub>1</sub>) antagonist and analog of somatostatin (IC<sub>50</sub> = 330 nM in a radioligand binding assay).<sup>1</sup> It inhibits increases in vascular permeability stimulated by substance P (Item No. 24035) in a concentration-dependent manner in isolated guinea pig trachea and main bronchi. Vapreotide (8 nmol per animal) reduces substance P-induced biting and scratching in mice. It increases reaction time to heat stimuli in the hot plate and tail flick tests when administered at a dose of 512 µg/kg, indicating antinociceptive activity in mice.<sup>2</sup> Vapreotide reduces plasma growth hormone (GH) increases induced by phenobarbital (Item Nos. 9001494 | 20987), morphine (Item No. ISO60147), and chlorpromazine (Item No. 16129), hypoglycemia-induced glucagon elevation, and glucose-induced insulin secretion in rats.<sup>3</sup> It also decreases tumor weight and volume in a hamster model of pancreatic ductal adenocarcinoma and suppresses tumor growth and reduces GH secretion in SK-ES-1 and MNNG/HOS osteosarcoma mouse xenograft models.<sup>4,5</sup>

### References

1. Bétoin, F., Advenier, C., Fardin, V., *et al.* *In vitro* and *in vivo* evidence for a tachykinin NK<sub>1</sub> receptor antagonist effect of vapreotide, an analgesic cyclic analog of somatostatin. *Eur. J. Pharmacol.* **279(2-3)**, 241-249 (1995).
2. Eschalier, A., Aumaître, O., Ardid, D., *et al.* Long-lasting antinociceptive effect of RC-160, a somatostatin analog, in mice and rats. *Eur. J. Pharmacol.* **199(1)**, 119-121 (1991).
3. Karashima, T., Cai, R.Z., and Schally, A.V. Effects of highly potent octapeptide analogs of somatostatin on growth hormone, insulin and glucagon release. *Life Sci.* **41(8)**, 1011-1019 (1987).
4. Paz-Bouza, J.I., Redding, T.W., and Schally, A.V. Treatment of nitrosamine-induced pancreatic tumors in hamsters with analogs of somatostatin and luteinizing hormone-releasing hormone. *Proc. Natl. Acad. Sci. U.S.A.* **84(4)**, 1112-1116 (1987).
5. Pinski, J., Schally, A.V., Halmos, G., *et al.* Somatostatin analog RC-160 inhibits the growth of human osteosarcomas in nude mice. *Int. J. Cancer.* **65(6)**, 870-874 (1996).

#### 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, 11/30/2022

#### 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