

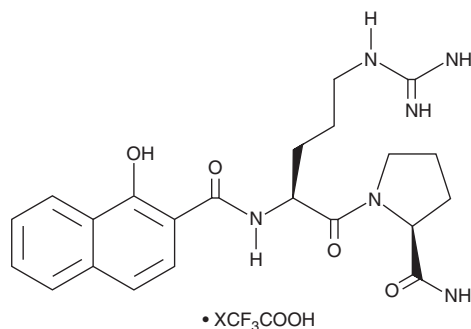
# PRODUCT INFORMATION



## APC 366 (trifluoroacetate salt)

Item No. 25990

**Formal Name:** N<sup>2</sup>-[(1-hydroxy-2-naphthalenyl) carbonyl]-L-arginyl-L-prolinamide, trifluoroacetate salt  
**MF:** C<sub>22</sub>H<sub>28</sub>N<sub>6</sub>O<sub>4</sub> • XCF<sub>3</sub>COOH  
**FW:** 440.5  
**Purity:** ≥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

APC 366 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the APC 366 (trifluoroacetate salt) in the solvent of choice. APC 366 (trifluoroacetate salt) is soluble in a 1:5 solution of ethanol:water at a concentration of 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

APC 366 is a mast cell tryptase inhibitor ( $K_i = 7.1 \mu\text{M}$  for human recombinant enzyme).<sup>1</sup> It reduces maximum contractile response to histamine in isolated bronchial tissue from ovalbumin-sensitized guinea pigs when used at a concentration of 1  $\mu\text{M}$ . Aerosolized APC 366 (5 mg in 1 ml of water) reduces airway resistance and increases dynamic lung compliance in *A. suum* antigen-challenged pigs.<sup>2</sup> It also reduces collagen content and expression of protease-activated receptor 2 (PAR-2) and  $\alpha$ -smooth muscle actin ( $\alpha$ -SMA) in rat liver in a model of bile duct ligation-induced liver fibrosis.<sup>3</sup>

### References

1. Barrios, V.E., Middleton, S.C., Kashem, M.A., *et al.* Tryptase mediates hyperresponsiveness in isolated guinea pig bronchi. *Life Sci.* **63(26)**, 2295-2303 (1998).
2. Sylvin, H., Dahlbäck, M., Van Der Ploeg, I., *et al.* The tryptase inhibitor APC-366 reduces the acute airway response to allergen in pigs sensitized to *Ascaris suum*. *Clin. Exp. Allergy* **32(6)**, 967-971 (2002).
3. Lu, J., Chen, B., Li, S., *et al.* Tryptase inhibitor APC 366 prevents hepatic fibrosis by inhibiting collagen synthesis induced by tryptase/protease-activated receptor 2 interactions in hepatic stellate cells. *Int. Immunopharmacol.* **20(2)**, 352-357 (2014).

#### 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, 11/29/2022

#### 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