PRODUCT INFORMATION



VUF 11207 (trifluoroacetate salt)

Item No. 19812

Formal Name:	N-[(2E)-3-(2-fluorophenyl)-2-methyl-2-propen-1-yl]-		
	3,4,5-trimethoxy-N-[2-(1-methyl-2-pyrrolidinyl)ethyl]-	o o	
	benzamide 2,2,2-trifluoroacetate		
MF:	$C_{27}H_{35}FN_2O_4 \bullet CF_3COOH$		
FW:	584.6		
Purity:	≥95%		/
UV/Vis.:	λ _{max} : 209 nm	\bigvee \bigvee \bigvee \bigvee \bigvee \bigcirc \bigcirc	
Supplied as:	A crystalline solid	O	
Storage:	-20°C	• CF ₃ COOH	
Stability:	≥4 years	5	

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

Laboratory Procedures

VUF 11207 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the VUF 11207 (trifluoroacetate salt) in the solvent of choice. VUF 11207 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of VUF 11207 (trifluoroacetate salt) in these solvents is approximately 20 mg/ml.

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 VUF 11207 (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of VUF 11207 (trifluoroacetate salt) in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

VUF 11207 is a potent agonist of the G protein-coupled chemokine receptor 7 (CXCR7; EC_{50} = 1.6 nM for β -arrestin recruitment to CXCR7).¹ It also reduces CXCR7 cell surface expression (EC₅₀ = 14.1 nM) in vitro.

Reference

1. Wijtmans, M., Maussang, D., Sirci, F., et al. Synthesis, modeling and functional activity of substituted styrene-amides as small-molecule CXCR7 agonists. Eur. J. Med. Chem. 51, 184-192 (2012).

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.

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, 11/02/2022

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