

PRODUCT INFORMATION



VUF 11207 (fumarate)

Item No. 44309

Formal Name: N-(3-(2-fluorophenyl)-2-methylallyl)-3,4,5-trimethoxy-N-(2-(1-methylpyrrolidin-2-yl)ethyl)benzamide, fumarate

MF: C₂₇H₃₅FN₂O₄ • C₄H₄O₄

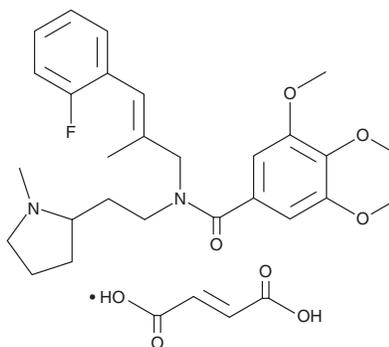
FW: 586.7

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

VUF 11207 (fumarate) is supplied as a solid. A stock solution may be made by dissolving the VUF 11207 (fumarate) in the solvent of choice, which should be purged with an inert gas. VUF 11207 (fumarate) is soluble (≥10 mg/ml) in ethanol and DMSO.

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 (fumarate) can be prepared by directly dissolving the solid in aqueous buffers. VUF 11207 (fumarate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

VUF 11207 is an agonist of chemokine (C-X-C motif) receptor 7 (CXCR7; EC₅₀ = 1.6 nM in a β-arrestin recruitment assay).¹ It also reduces CXCR7 cell surface expression in HEK293 cells expressing the human receptor (EC₅₀ = 14.1 nM). VUF 11207 (5 nM) induces the migration and proliferation of primary human pericytes isolated from patients with pulmonary arterial hypertension.² *In vivo*, VUF 11207 (5 mg/kg) increases thoracic adventitia thickness and aorta collagen deposition in a mouse model of hypertension induced by angiotensin II.³

References

1. Wijtmans, M., Maussang, D., Sirici, 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).
2. Bordenave, J., Tu, L., Berrebeh, N., *et al.* Lineage tracing reveals the dynamic contribution of pericytes to the blood vessel remodeling in pulmonary hypertension. *Arterioscler. Thromb. Vasc. Biol.* **40**(3), 766-782 (2020).
3. Song, B., Chen, D., Liu, Z., *et al.* Stromal cell-derived factor-1 exerts opposing roles through CXCR4 and CXCR7 in angiotensin II-induced adventitial remodeling. *Biochem. Biophys. Res. Commun.* **594**, 38-45 (2022).

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, 12/17/2025

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