

PRODUCT INFORMATION



Befetupitant

Item No. 43004

CAS Registry No.: 290296-68-3

Formal Name: N, α , α -trimethyl-N-[4-(2-methylphenyl)-6-(4-morpholinyl)-3-pyridinyl]-3,5-bis(trifluoromethyl)-benzeneacetamide

Synonym: Ro 67-5930

MF: $C_{29}H_{29}F_6N_3O_2$

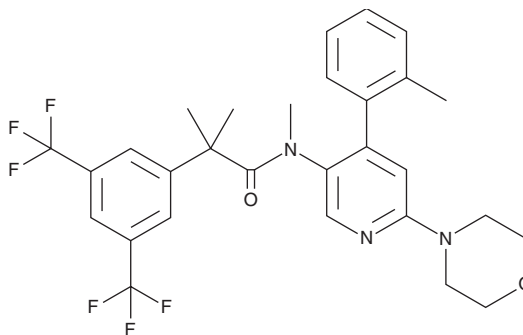
FW: 565.6

Purity: $\geq 98\%$

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

Befetupitant is supplied as a solid. A stock solution may be made by dissolving the befetupitant in the solvent of choice, which should be purged with an inert gas. Befetupitant is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Befetupitant is a neurokinin-1 (NK_1) receptor antagonist ($K_i = 1$ nM).¹ It inhibits GR73632-induced foot tapping in gerbils ($ID_{50} = 0.2$ mg/kg). Topical administration of befetupitant (0.4 or 1.6 mg/ml for four days) decreases hemangiogenesis and lymphangiogenesis but induces corneal opacity, stromal leukocyte infiltration, and corneal toxicity in a mouse model of corneal alkali burn.²

References

- Hoffmann, T., Börs, M., Stadler, H., *et al.* Design and synthesis of a novel, achiral class of highly potent and selective, orally active neurokinin-1 receptor antagonists. *Bioorg. Med. Chem. Lett.* **16**(5), 1362-1365 (2006).
- Bignami, F., Giacomini, C., Lorusso, A., *et al.* NK1 receptor antagonists as a new treatment for corneal neovascularization. *Invest. Ophthalmol. Vis. Sci.* **55**(10), 6783-6794 (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, 03/20/2025

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