

PRODUCT INFORMATION



Netupitant N-oxide

Item No. 34005

CAS Registry No.: 910808-11-6

Formal Name: N, α , α -trimethyl-N-[6-(4-methyl-4-oxido-1-piperazinyl)-4-(2-methylphenyl)-3-pyridinyl]-3,5-bis(trifluoromethyl)-benzeneacetamide

MF: $C_{30}H_{32}F_6N_4O_2$

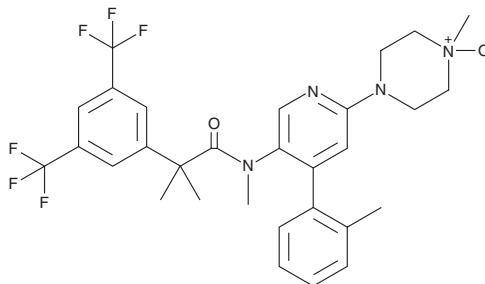
FW: 594.6

Purity: $\geq 95\%$

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

Netupitant N-oxide is supplied as a solid. A stock solution may be made by dissolving the netupitant N-oxide in the solvent of choice, which should be purged with an inert gas. Netupitant N-oxide is slightly soluble in methanol and DMSO (heated).

Description

Netupitant N-oxide is a metabolite of the neurokinin-1 (NK_1) receptor antagonist netupitant (Item No. 23809).¹ Netupitant N-oxide is formed from netupitant by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Lanzarotti, C. and Rossi, G. Effect of netupitant, a highly selective NK_1 receptor antagonist, on the pharmacokinetics of midazolam, erythromycin and dexamethasone. *Support Care Cancer* **21**(10), 2783-2791 (2013).

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, 11/19/2024

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