

PRODUCT INFORMATION



N-Nitroso Fluoxetine

Item No. 34879

CAS Registry No.: 150494-06-7

Formal Name: N-methyl-N-nitroso-γ-[4-(trifluoromethyl)phenoxy]-benzenepropanamine

MF: C₁₇H₁₇F₃N₂O₂

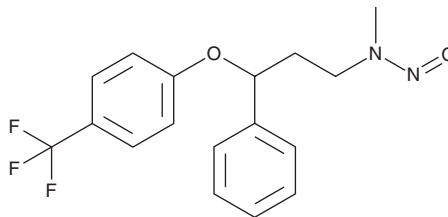
FW: 338.3

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

N-Nitroso fluoxetine is supplied as a solid. A stock solution may be made by dissolving the N-nitroso fluoxetine in the solvent of choice, which should be purged with an inert gas. N-Nitroso fluoxetine is slightly soluble in chloroform and dichloromethane.

Description

N-Nitroso fluoxetine is a derivative of fluoxetine.¹ It is genotoxic to *S. typhimurium* when used at concentrations ranging from 0.06 to 0.12 mg/ml.

Reference

1. Özhan, G. and Alpertunga, B. Genotoxic activities of drug-nitrite interaction products. *Drug Chem. Toxicol.* **26**(4), 295-308 (2003).

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, 11/28/2022

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