

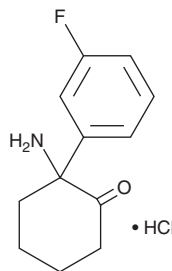
PRODUCT INFORMATION



3-fluoro Deschloronorketamine (hydrochloride)

Item No. 42932

Formal Name: 2-amino-2-(3-fluorophenyl)cyclohexan-1-one, monohydrochloride
Synonyms: 3-fluoro Deschloroketamine metabolite, Fluorexetamine metabolite
MF: $C_{12}H_{14}FNO \cdot HCl$
FW: 243.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Description

3-fluoro Deschloronorketamine (hydrochloride) (Item No. 42932) is an analytical reference standard categorized as an arylcyclohexylamine. 3-fluoro Deschloronorketamine is a potential metabolite of 3-fluoro deschloroketamine (Item No. 29756) and fluorexetamine (Item No. 35118) based on the published metabolism of 2-fluoro deschloroketamine (Item Nos. 41176 | 19786).¹ At the time 3-fluoro deschloronorketamine (hydrochloride) (Item No. 42932) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Mestria, S., Odoardi, S., Bioss, G., *et al.* Method development for the identification of methoxpropamine, 2-fluoro-deschloroketamine and deschloroketamine and their main metabolites in blood and hair and forensic application. *Forensic Sci. Int.* **323**, 110817 (2021).

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, 02/13/2025

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