

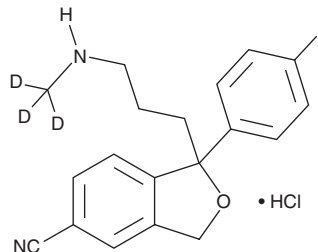
PRODUCT INFORMATION



Desmethylcitalopram-d₃ (hydrochloride)

Item No. 15840

CAS Registry No.: 1189650-18-7
Formal Name: 1-(4-fluorophenyl)-1,3-dihydro-1-[3-((methyl-d₃)amino)propyl]-5-isobenzofurancarboxitrile, monohydrochloride
MF: C₁₉H₁₆D₃FN₂O • HCl
FW: 349.9
Chemical Purity: ≥98% (Desmethylcitalopram)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
UV/Vis.: λ_{max}: 238, 284 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Desmethylcitalopram-d₃ (hydrochloride) (Item No. 15840) is intended for use as an internal standard for the quantification of desmethylcitalopram (Item No. 15905) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Desmethylcitalopram is an active metabolite of citalopram (Item Nos. 23252 | 14572).¹ This product is intended for research and forensic applications.

Reference

1. Weisskopf, E., Panchaud, A., Nguyen, K.A., *et al.* Stereoselective determination of citalopram and desmethylcitalopram in human plasma and breast milk by liquid chromatography tandem mass spectrometry. *J. Pharm. Biomed. Anal.* **131**, 233-245 (2016).

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, 06/23/2023

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