

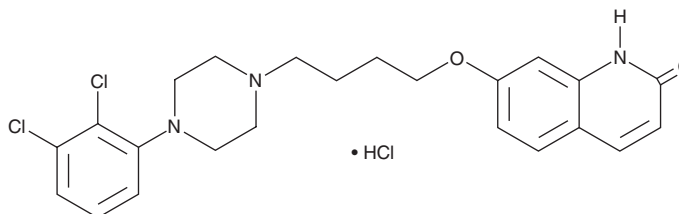
PRODUCT INFORMATION



Dehydro Aripiprazole (hydrochloride)

Item No. 19539

CAS Registry No.: 1008531-60-9
Formal Name: 7-[4-[4-(2,3-dichlorophenyl)-1-piperazinyl]butoxy]-2(1H)-quinolinone, monohydrochloride
MF: C₂₃H₂₅Cl₂N₃O₂ • HCl
FW: 482.8
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 258, 326, 341 nm
Supplied as: A crystalline 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

Dehydro aripiprazole (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the dehydro aripiprazole (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Dehydro aripiprazole (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of dehydro aripiprazole (hydrochloride) in these solvents is approximately 20 and 25, respectively.

Dehydro aripiprazole (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dehydro aripiprazole (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Dehydro aripiprazole (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dehydro Aripiprazole is a metabolite of the atypical antipsychotic aripiprazole (Item No. 19989).¹ It is formed from aripiprazole by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2D6.

Reference

1. Spina, E. and de Leon, J. Metabolic drug interactions with newer antipsychotics: A comparative review. *Basic Clin. Pharmacol. Toxicol.* **100**(1), 4-22 (2007).

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, 09/08/2023

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