

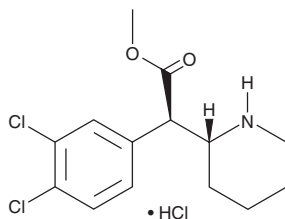
PRODUCT INFORMATION



(±)-*threo*-3,4-Dichloromethylphenidate (hydrochloride)

Item No. 18588

CAS Registry No.: 214149-42-5
Formal Name: (αR,2R)-*rel*-α-(3,4-dichlorophenyl)-2-piperidineacetic acid, methyl ester, monohydrochloride
MF: C₁₄H₁₇Cl₂NO₂ • HCl
FW: 338.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

(±)-*threo*-3,4-Dichloromethylphenidate (hydrochloride) (Item No. 18588) is an analytical reference material that is structurally categorized as an amphetamine. Its *threo*-diastereomer, 3,4-CTMP, can inhibit dopamine uptake through the dopamine transporter with an IC₅₀ value of 65 nM in rat striatal tissue.¹ This product is intended for research and forensic studies.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

Reference

1. Wayment, H.K., Deutsch, H., Schveri, M.M., *et al.* Effects of methylphenidate analogues on phenethylamine substrates for the striatal dopamine transporter: Potential as amphetamine antagonists? *J. Neurochem.* **72**, 1266-1274 (1999).

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, 08/14/2019

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