

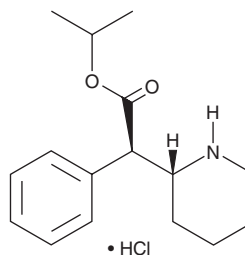
PRODUCT INFORMATION



(±)-*threo*-Isopropylphenidate (hydrochloride)

Item No. 19264

CAS Registry No.: 1262795-94-7
Formal Name: (αR,2R)-*rel*-α-phenyl-2-piperidineacetic acid, 1-methylethyl ester, monohydrochloride
Synonyms: Dimethylphenidate, Ritalinic Acid isopropyl ester
MF: C₁₆H₂₃NO₂ • HCl
FW: 297.8
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*-Isopropylphenidate (hydrochloride) (Item No. 19264) is an analytical reference standard that is structurally categorized as a phenethylamine. It is closely related to methylphenidate (Item No. 11639) and has been speculated to act as both a dopamine reuptake inhibitor and norepinephrine reuptake inhibitor.¹ This product is intended for forensic and research uses.

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

Reference

1. Markowitz, J.S., Zhu, H.-J., and Patrick, K.S. Isopropylphenidate: An ester homolog of methylphenidate with sustained and selective dopaminergic activity and reduced drug interaction liability. *J. Child Adolesc. Psychopharmacol.* **23**(10), 648-654 (2013).

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, 08/26/2019

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