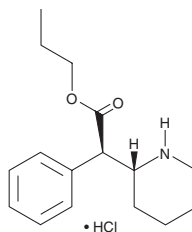


PRODUCT INFORMATION

(±)-*threo*-Propylphenidate (hydrochloride)

Item No. 19268

CAS Registry No.: 2514950-54-8
Formal Name: (R,R)-*rel*-α-phenyl-2-piperidineacetic acid, propyl ester, monohydrochloride
MF: C₁₆H₂₃NO₂ • HCl
FW: 297.8
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

(±)-*threo*-Propylphenidate (hydrochloride) (Item No. 19268) is an analytical reference standard that is structurally categorized as a phenethylamine. It is closely related to methylphenidate (Item No. 11639) but with the methyl ester replaced by a propyl ester. This product is intended for forensic and research uses.

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, 06/19/2023

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