

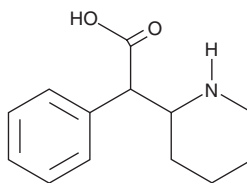
PRODUCT INFORMATION



Ritalinic Acid

Item No. 15915

CAS Registry No.: 19395-41-6
Formal Name: α -phenyl-2-piperidineacetic acid
MF: C₁₃H₁₇NO₂
FW: 219.3
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Description

Ritalinic acid is an inactive, major metabolite of methylphenidate (Item No. 11639), a schedule II drug in the United States commonly used as a psychostimulant for the treatment of attention-deficit hyperactivity disorder, also known as ADHD.¹ Ritalinic acid is reported to reach plasma concentrations of 1-2 orders of magnitude greater than that of the parent drug and has been monitored in urine to ensure compliance with treatment programs.² This product is intended for forensic and biological research purposes.

References

1. Markowitz, J.S., Devane, C.L., Boulton, D.W., *et al.* Ethylphenidate formation in human subjects after the administration of a single dose of methylphenidate and ethanol. *Drug Metab. Dispos.* **28(6)**, 620-624 (2000).
2. del Mar Ramirez Fernandez, M., Laloup, M., Wood, M., *et al.* Liquid chromatography-tandem mass spectrometry method for the simultaneous analysis of multiple hallucinogens, chlorpheniramine, ketamine, ritalinic acid, and metabolites, in urine. *J. Anal. Toxicol.* **31(8)**, 497-504 (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, 11/21/2022

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