

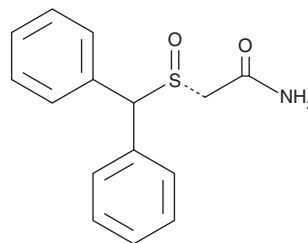
PRODUCT INFORMATION



(R)-Modafinil

Item No. 29910

CAS Registry No.: 112111-43-0
Formal Name: 2-[(R)-(diphenylmethyl)sulfinyl]-acetamide
Synonyms: CEP-10952, CRL 40982, (-)-Modafinil
MF: C₁₅H₁₅NO₂S
FW: 273.4
Purity: ≥98%
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

(R)-Modafinil (Item No. 29910) is an analytical reference standard categorized as a nootropic and stimulant.¹⁻³ (R)-Modafinil substitutes for cocaine (Item Nos. ISO60176 | 16186 | 22165) in a two-lever drug discrimination test. Formulations containing (R)-modafinil have been used in the treatment of narcolepsy, obstructive sleep apnea, and shift work disorder. (R)-Modafinil is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Loland, C.J., Mereu, M., Okunola, O.M., *et al.* R-Modafinil (armodafinil): A unique dopamine uptake inhibitor and potential medication for psychostimulant abuse. *Biol. Psychiatry* **72**(5), 405-413 (2012).
2. Schmitt, K.C. and Reith, M.E.A. The atypical stimulant and nootropic modafinil interacts with the dopamine transporter in a different manner than classical cocaine-like inhibitors. *PLoS One* **6**(10), e25790 (2011).
3. Czeisler, C.A., Walsh, J.K., Wesnes, K.A., *et al.* Armodafinil for treatment of excessive sleepiness associated with shift work disorder: A randomized controlled study. *Mayo Clin. Proc.* **84**(11), 958-972 (2009).

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, 07/06/2023

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