

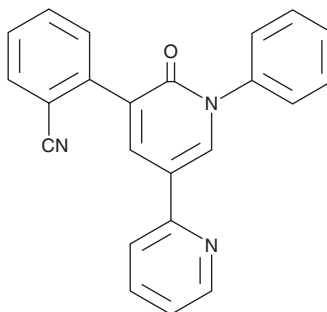
PRODUCT INFORMATION



Perampanel

Item No. 23003

CAS Registry No.: 380917-97-5
Formal Name: 2-(1',6'-dihydro-6'-oxo-1'-phenyl[2,3'-bipyridin]-5'-yl)-benzonitrile
MF: C₂₃H₁₅N₃O
FW: 349.4
Purity: ≥98%
UV/Vis.: λ_{max}: 295, 347 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

Perampanel (Item No. 23003) is an analytical reference standard categorized as an anticonvulsant.¹ Formulations containing perampanel induce sedation and euphoria at high doses.² Perampanel is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

References

1. Hanada, T., Hashizume, Y., Tokuhara, N., *et al.* Perampanel: A novel, orally active, noncompetitive AMPA-receptor antagonist that reduces seizure activity in rodent models of epilepsy. *Epilepsia* **52**(7), 1331-1340 (2011).
2. Hawkins, K.L. and Gidal, B.E. When adverse effects are seen as desirable: Abuse potential of the newer generation antiepileptic drugs. *Epilepsy Behav.* **77**, 62-72 (2017).

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