

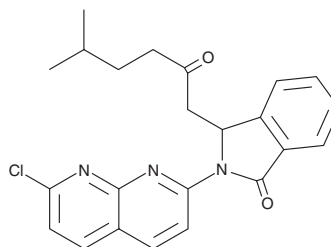
PRODUCT INFORMATION



Pagoclone

Item No. 27749

CAS Registry No.: 133737-48-1
Formal Name: 2-(7-chloro-1,8-naphthyridin-2-yl)-2,3-dihydro-3-(5-methyl-2-oxohexyl)-1H-isoindol-1-one
Synonym: RP 59037
MF: C₂₃H₂₂ClN₃O₂
FW: 407.9
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

Pagoclone (Item No. 27749) is an analytical reference standard categorized as a cyclopyrrolone.¹ Pagoclone has anxiolytic-like, anticonvulsant, and anti-aggressive activities in rodents. This product is intended for research and forensic applications.

Reference

1. Doble, A., Canton, T., Piot, D.O., *et al.* RP 59037 and RP 60503: Anxiolytic cyclopyrrolone derivatives with low sedative potential. Interaction with the γ -aminobutyric acid_A/benzodiazepine receptor complex and behavioral effects in the rodent. *J. Pharmacol. Exp. Ther.* **266**(3), 1213-1226 (1993).

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, 07/24/2023

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