

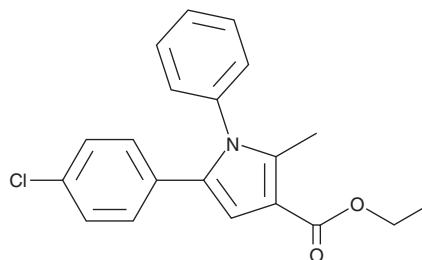
PRODUCT INFORMATION



NAP-1

Item No. 28719

CAS Registry No.: 131721-28-3
Formal Name: 5-(4-chlorophenyl)-2-methyl-1-phenyl-1H-pyrrole-3-carboxylic acid, ethyl ester
MF: C₂₀H₁₈ClNO₂
FW: 339.8
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 281 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

NAP-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the NAP-1 in the solvent of choice. NAP-1 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of NAP-1 in these solvents is approximately 10 and 30 mg/ml, respectively.

NAP-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NAP-1 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NAP-1 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NAP-1 is a compound with anesthetic activity.¹ It increases paired-pulse inhibition in the CA1 region of rat hippocampal brain slices when used at a concentration of 100 μM. NAP-1 also induces loss of righting reflex in tadpoles (EC₅₀ = 0.53 μM).

Reference

1. Bertaccini, E.J. and Davies, M.F. Novel methods, compounds, and compositions for anesthesia. *Stanford Univ. and US Govrnt. Dept. Vet. Affairs PCT/US2015/056067* (2016).

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, 03/19/2024

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