

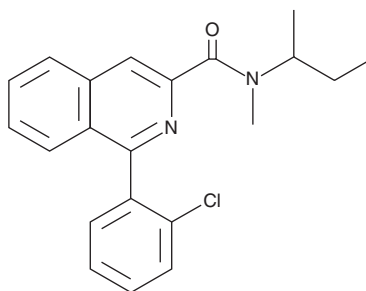
PRODUCT INFORMATION



PK 11195

Item No. 10525

CAS Registry No.: 85532-75-8
Formal Name: 1-(2-chlorophenyl)-N-methyl-N-(1-methylpropyl)-3-isoquinolinecarboxamide
Synonym: RP 52028
MF: C₂₁H₂₁ClN₂O
FW: 352.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

PK 11195 is supplied as a solid. A stock solution may be made by dissolving the PK 11195 in the solvent of choice. PK 11195 is soluble in organic solvents such as methanol, chloroform, and acetonitrile, which should be purged with an inert gas. The solubility of PK 11195 in these solvents is approximately 16, 33, and 10 mg/ml, respectively.

Description

Benzodiazepines bind two types of receptors, the central γ -aminobutyric acid A receptor and the peripheral translocator protein (TSPO). PK 11195 binds the peripheral benzodiazepine receptor, TSPO, with selectivity and high affinity ($K_i = 3.1$ nM in cerebellum, 4.1 nM in spinal cord).^{1,2} PK 11195, which lacks the 7-member heterocycle of diazepines, blocks binding of typical benzodiazepines to TSPO.¹ In addition, the binding of labeled PK 11195 to tissues is used to detect the presence of TSPO by various methods.^{2,3} PK 11195 has been used to study the ligand binding site of the *B. cereus* TSPO.⁴

References

1. Watanabe, Y., Shibuya, T., Khatami, S., *et al.* Comparison of typical and atypical benzodiazepines on the central and peripheral benzodiazepine receptors. *Japan J. Pharmacol.* **42**, 189-197 (1986).
2. Pike, V.W., Halldin, C., Crouzel, C., *et al.* Radioligands for PET studies of central benzodiazepine receptors and PK (peripheral benzodiazepine) binding sites-current status. *Nucl. Med. Biol.* **20**(4), 503-525 (1993).
3. Olson, J.M.M., Junck, L., Young, A.B., *et al.* Isoquinoline and peripheral-type benzodiazepine binding in gliomas: Implications for diagnostic imaging. *Cancer Res.* **48**, 5837-5841 (1988).
4. Guo, Y., Kalathur, R.C., Liu, Q., *et al.* Protein structure. Structure and activity of tryptophan-rich TSPO proteins. *Science* **347**(6221), 551-555 (2015).

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, 03/22/2018

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