

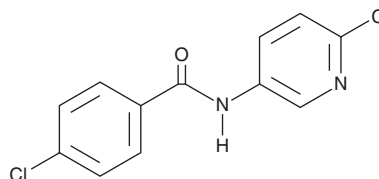
PRODUCT INFORMATION



ICA 110381

Item No. 37294

CAS Registry No.: 325457-99-6
Formal Name: 4-chloro-N-(6-chloro-3-pyridinyl)-benzamide
MF: C₁₂H₈Cl₂N₂O
FW: 267.1
Purity: ≥98%
UV/Vis.: λ_{max}: 268 nm
Supplied as: A 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

ICA 110381 is supplied as a solid. A stock solution may be made by dissolving the ICA 110381 in the solvent of choice, which should be purged with an inert gas. ICA 110381 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ICA 110381 in these solvents is approximately 1, 25, and 10 mg/ml, respectively.

Description

ICA 110381 is an activator of voltage-gated potassium channel K_v7.2.¹ It induces rubidium efflux in an agonist-induced rubidium efflux assay using CHO cells (EC₅₀ = 0.38 μM for the human channel). ICA 110381 also inhibits rubidium efflux from CHO cells expressing K_v7.1 (EC₅₀ = 15 μM for the human channel). It reduces seizure number and increases the current threshold for seizure induction by amygdala-kindling in rats when administered at a dose of 30 mg/kg.²

References

1. Amato, G., Roeloffs, R., Rigdon, G.C., *et al.* N-Pyridyl and pyrimidine benzamides as KCNQ2/Q3 potassium channel openers for the treatment of epilepsy. *ACS Med. Chem. Lett.* **2(6)**, 481-484 (2011).
2. Boehlen, A., Schwake, M., Dost, R., *et al.* The new KCNQ2 activator 4-Chlor-N-(6-chlor-pyridin-3-yl)-benzamid displays anticonvulsant potential. *Br. J. Pharmacol.* **168(5)**, 1182-1200 (2012).

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, 10/17/2023

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