

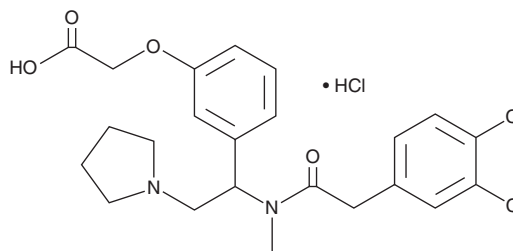
PRODUCT INFORMATION



ICI 204448 (hydrochloride)

Item No. 45206

CAS Registry No.: 121264-04-8
Formal Name: 2-[3-[1-[2-(3,4-dichlorophenyl)acetyl]methylamino]-2-(1-pyrrolidinyl)ethyl]phenoxy]-acetic acid, monohydrochloride
MF: C₂₃H₂₆Cl₂N₂O₄ • HCl
FW: 501.8
Purity: ≥98%
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

ICI 204448 is supplied as a solid. A stock solution may be made by dissolving the ICI 204448 in the solvent of choice, which should be purged with an inert gas. ICI 204448 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

ICI 204448 is a κ -opioid receptor agonist.¹ It binds to the κ -opioid receptor ($K_i = 2.69$ nM) and is 162- and 69-fold selective for the κ -opioid receptor over the μ - and δ -opioid receptors in membranes prepared from cells expressing the human κ -, μ - and δ -opioid receptors, respectively. ICI 204448 reduces acetic acid-induced writhing and chloroquine-induced scratching in mice.^{1,2} It also reduces infarct size in rat hearts when administered ten minutes prior to the onset of ischemia.³

References

1. Kumar, V., Guo, D., Cassel, J.A., *et al.* Synthesis and evaluation of novel peripherally restricted κ -opioid receptor agonists. *Bioorg. Med. Chem. Lett.* **15**(4), 1091-1095 (2005).
2. Inan, S. and Cowan, A. Kappa opioid agonists suppress chloroquine-induced scratching in mice. *Eur. J. Pharmacol.* **502**(3), 233-237 (2004).
3. Peart, J.N., Gross, E.R., and Gross, G.J. Effect of exogenous kappa-opioid receptor activation in rat model of myocardial infarction. *J. Cardiovasc. Pharmacol.* **43**(3), 410-415 (2004).

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, 04/10/2026

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