

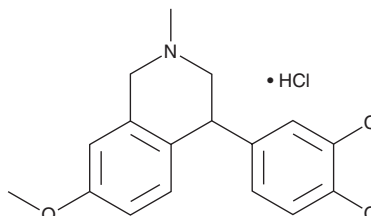
PRODUCT INFORMATION



Diclofensine (hydrochloride)

Item No. 15355

CAS Registry No.: 34041-84-4
Formal Name: 4-(3,4-dichlorophenyl)-1,2,3,4-tetrahydro-7-methoxy-2-methyl-isoquinoline, monohydrochloride
Synonym: Ro 8-4650
MF: C₁₇H₁₇Cl₂NO • HCl
FW: 358.7
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

Diclofensine is a potent inhibitor of monoamine reuptake, blocking the uptake of dopamine, noradrenaline, and serotonin by rat brain synaptosomes with IC₅₀ values of 0.74, 2.3, and 3.7 nM, respectively.¹ While it has applications as an antidepressant, diclofensine increases locomotor activity, decreases food intake, and has abuse potential.²⁻⁵ This product is intended for research and forensic applications.

References

1. Hyttel, J. and Larsen, J.J. Neurochemical profile of Lu 19-005, a potent inhibitor of uptake of dopamine, noradrenaline, and serotonin. *J. Neurochem.* **44(5)**, 1615-1622 (1985).
2. Jackson, S.H., Turner, P., and Ehsanullah, R.S. A comparison of diclofensine and desmethylinipramine using tyramine pressor tests in normal subjects. *Br. J. Clin. Pharmacol.* **16(4)**, 427-429 (1983).
3. Cilug, J., Ehsanullah, R.S., Hallett, C., et al. A clinical pharmacological comparison of diclofensine (Ro 8-4650) with nomifensine and amitriptyline in normal human volunteers. *Br. J. Clin. Pharmacol.* **15(5)**, 537-543 (1983).
4. Lamb, R.J. and Griffiths, R.R. Self-administration in baboons and the discriminative stimulus effects in rats of bupropion, nomifensine, diclofensine and imipramine. *Psychopharmacology (Berl)* **102(2)**, 183-190 (1990).
5. Nakachi, N., Kiuchi, Y., Inagaki, M., et al. Effects of various dopamine uptake inhibitors on striatal extracellular dopamine levels and behaviours in rats. *Eur. J. Pharmacol.* **281(2)**, 195-203 (1995).

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, 06/27/2023

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