

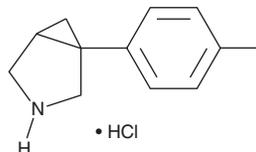
PRODUCT INFORMATION



Bicifadine (hydrochloride)

Item No. 43813

CAS Registry No.: 66504-75-4
Formal Name: 1-(4-methylphenyl)-3-azabicyclo[3.1.0]hexane,
monohydrochloride
MF: C₁₂H₁₅N • HCl
FW: 209.7
Purity: ≥90%
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

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

Description

Bicifadine is an inhibitor of the serotonin transporter (SERT), dopamine transporter (DAT), and norepinephrine transporter (NET).¹ It binds these transporters (K_s = 2.4, 5.2, and 5.4 μM, respectively) and inhibits serotonin, dopamine, and norepinephrine reuptake in intact cells expressing the human transporters (IC₅₀s = 117, 910, and 55 nM, respectively). Bicifadine increases the mechanical withdrawal threshold in a rat model of acute inflammatory pain induced by subplantar injection of brewer's yeast (ED₅₀ = 9.2 mg/kg) and decreases writhing in mouse model of persistent abdominal pain induced by phenyl-*p*-quinone (PPQ; ED₅₀ = 13.2 mg/kg). It also reduces paw licking in the first and second phases of the formalin test in mice and rats when administered 30 minutes prior to formalin.

Reference

1. Basile, A.S., Janowsky, A., Golembiowska, K., *et al.* Characterization of the antinociceptive actions of bicifadine in models of acute, persistent, and chronic pain. *J. Pharmacol. Exp. Ther.* **321(3)**, 1208-1225 (2007).

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, 02/06/2026

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