

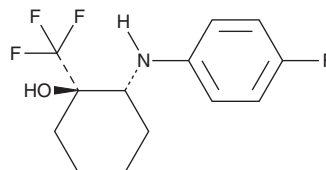
PRODUCT INFORMATION



BAY-390

Item No. 41313

CAS Registry No.: 2741956-55-6
Formal Name: 2R-[(4-fluorophenyl)amino]-1R-(trifluoromethyl)-cyclohexanol
MF: $C_{13}H_{15}F_4NO$
FW: 277.3
Purity: $\geq 95\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

BAY-390 is supplied as a solution in ethanol. A stock solution may be made by dissolving the BAY-390 in the solvent of choice, which should be purged with an inert gas. BAY-390 is slightly soluble (0.1-1 mg/ml) in acetonitrile and ethanol.

Description

BAY-390 is an inhibitor of transient receptor potential ankyrin 1 (TRPA1; $IC_{50}s = 16$ and 63 nM for the human and rat receptor, respectively).¹ It is selective for TRPA1 over a variety of receptors, ion channels, transporters, and enzymes at $10 \mu M$ but weakly inhibits estrogen receptor α (ER α), progesterone receptor B (PR-B), and the dopamine transporter (DAT; $IC_{50}s = 5,000$, $5,000$, and $1,200$ nM, respectively). BAY-390 (3-10 mg/kg) reduces the number of flinches and the duration of paw licking in a rat model of nociceptive behavior induced by the TRPA1 agonist cinnamaldehyde and reduces inflammatory pain induced by complete Freund's adjuvant (CFA) in rats. It also reverses mechanical allodynia in a rat model of neuropathic pain induced by spinal nerve ligation when administered at a dose of 90 mg/kg twice per day.

Reference

1. Mesch, S., Walter, D., Laux-Biehlmann, A., et al. Discovery of BAY-390, a selective CNS penetrant chemical probe as transient receptor potential ankyrin 1 (TRPA1) antagonist. *J. Med. Chem.* **66**(2), 1583-1600 (2023).

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, 05/11/2026

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