

# PRODUCT INFORMATION

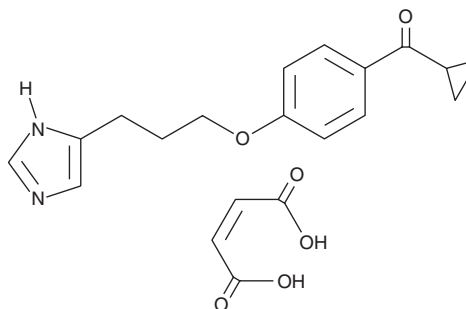


## Ciproxifan (maleate)

Item No. 29513

**CAS Registry No.:** 184025-19-2  
**Formal Name:** cyclopropyl[4-[3-(1H-imidazol-4-yl)propoxy]phenyl]-methanone (2Z)-2-butenedioate

**Synonym:** FUB-359  
**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>O<sub>2</sub> • C<sub>4</sub>H<sub>4</sub>O<sub>4</sub>  
**FW:** 386.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 272 nm  
**Supplied as:** A crystalline 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

Ciproxifan (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the ciproxifan (maleate) in the solvent of choice, which should be purged with an inert gas. Ciproxifan (maleate) is soluble in DMSO.

### Description

Ciproxifan is a histamine H<sub>3</sub> receptor antagonist (K<sub>i</sub> = 0.5 nM for inhibition of histamine release in rat synaptosomal preparations).<sup>1</sup> It is selective for histamine H<sub>3</sub> over histamine H<sub>1</sub> and H<sub>2</sub>, M<sub>3</sub> muscarinic, and α<sub>1D</sub>- and β<sub>1</sub>-adrenergic receptors (K<sub>i</sub>s = 3.15-25 μM), as well as the serotonin (5-HT) receptor subtypes 5-HT<sub>1B</sub>, 5-HT<sub>2A</sub>, 5-HT<sub>3</sub>, and 5-HT<sub>4</sub> (K<sub>i</sub>s = 1.58-15 μM) in radioligand binding assays. Ciproxifan (3-300 nM) inhibits relaxation in precontracted isolated guinea pig ileal longitudinal muscle induced by R(-)-α-methylhistamine (Item No. 25601). *In vivo*, ciproxifan reduces R(-)-α-methylhistamine-induced water consumption in rats (ED<sub>50</sub> = 0.09 mg/kg). It increases wake episode duration and latency to fall asleep in rats when administered at a dose of 2 mg/kg.

### Reference

1. Ligneau, X., Lin, J., Vanni-Mercier, G., *et al.* Neurochemical and behavioral effects of ciproxifan, a potent histamine H<sub>3</sub>-receptor antagonist. *J. Pharmacol. Exp. Ther.* **287**(2), 658-666 (1998).

**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, 12/13/2022

### 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