

# Trifluoperazine . dihydrochloride

Antipsychotic

Facilitates agonist induced elevation of cytosolic calcium at low concentrations (10 $\mu$ M). Antagonizes calmodulin at higher concentrations. Antipsychotic and sedative dopamine antagonist. Potent and irreversible inhibitor of cAMP-gated cationic channels. Inhibits DNA repair. Cell permeable. Potent neurotoxin.

Citations: 10

[View Online »](#)

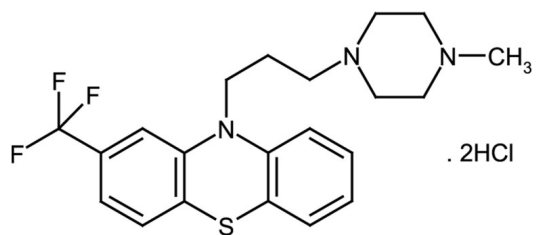
Ordering Information

[Order Online »](#)

ALX-550-310-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White to off-white powder.
CAS	440-17-5
Couple Target	Calmodulin, CNG channel, Dopamine
Couple Type	Inhibitor
Formula	$C_{21}H_{24}F_3N_3S \cdot 2HCl$
Identity	Identity determined by IR.
MI	14: 9680
MW	407.5 . 73.0
Purity	≥98% (UHPLC)
RTECS	SP1750000
Solubility	Soluble in water (50mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)