

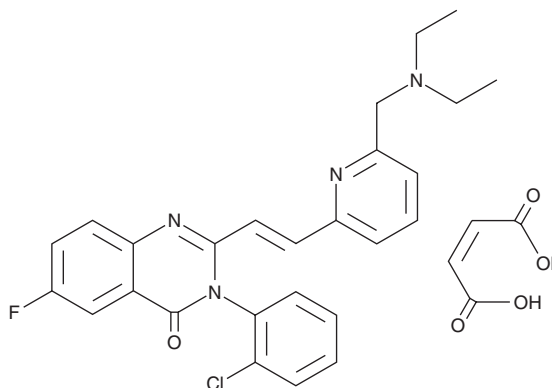
PRODUCT INFORMATION



CP-392,110 (maleate)

Item No. 20473

CAS Registry No.: 199656-46-7
Formal Name: 3-(2-chlorophenyl)-2-[2-[6-[(diethylamino)methyl]-2-pyridinyl]ethenyl]-6-fluoro-4(3H)-quinazolinone, (2Z)-2-butenedioate
MF: C₂₆H₂₄ClFN₄O • C₄H₄O₄
FW: 579.0
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 334 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CP-392,110 (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the CP-392,110 (maleate) in the solvent of choice, which should be purged with an inert gas. CP-392,110 (maleate) is soluble in the organic solvent DMSO.

Description

CP-392,110 is an AMPA receptor antagonist.¹ It inhibits AMPA receptor-mediated calcium uptake in rat cerebellar neurons (IC₅₀ = 74 nM). CP-392,110 inhibits seizures induced by pentylenetetrazol (Item No. 18682) or AMPA (Item No. 14571) in mice with ID₅₀ values of 3.8 and 3.6 mg/kg, respectively.

Reference

1. Welch, W.M., Ewing, F.E., Huang, J., *et al.* Atropisomeric quinazolin-4-one derivatives are potent noncompetitive α-amino-3-hydroxy-5-methyl-4-isoxazolepropionic acid (AMPA) receptor antagonists. *Bioorg. Med. Chem.* **11(2)**, 177-181 (2001).

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, 11/04/2019

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