

Tamoxifen . citrate

PKC inhibitor

Antiestrogen with anti-cancer activity. Selective protein kinase C (PKC) inhibitor and induces PKC epsilon translocation. Induces apoptosis in estrogen receptor-negative human cancer cell lines.

Citations: 25

[View Online »](#)

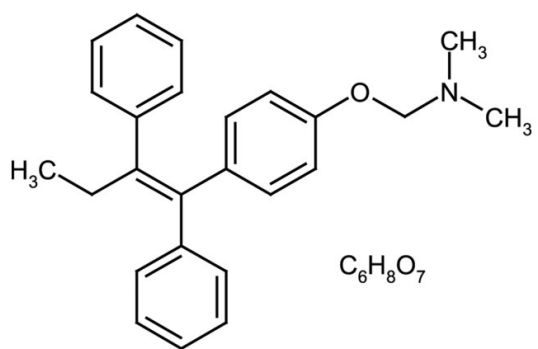
Ordering Information

[Order Online »](#)

ALX-550-095-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (Z)-2-[4-(1,2-Diphenyl-1-butenyl)phenoxy]-N,N-dimethyl-ethanamine . citrate

Appearance White to off-white powder.

CAS 54965-24-1

Couple Target PKC

Couple Type Inhibitor

Formula $C_{26}H_{29}NO \cdot C_6H_8O_7$

Identity Determined by IR.

MI 14: 9048

MW 371.5 . 192.1

Purity ≥99% (Assay)

RTECS KH2387000

Solubility Soluble in methanol or ethanol; very slightly soluble in water or acetone.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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