

(Z)-4-Hydroxytamoxifen

Tamoxifen metabolite

Metabolite of tamoxifen (Prod. No. ALX-550-095). Exhibits a more potent estrogen agonist/antagonist activity than the parent drug. Has antioxidant properties. Intramembranous inhibitor of lipid peroxidation.

Citations: 11

[View Online »](#)

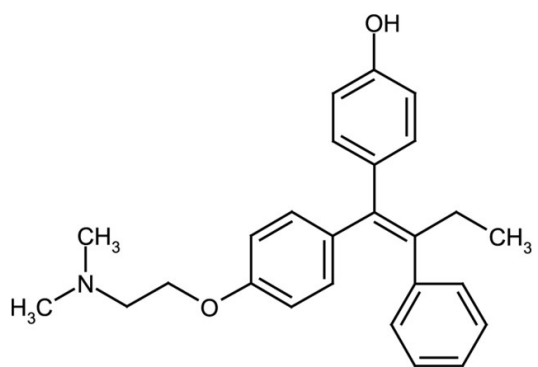
Ordering Information

[Order Online »](#)

ALX-550-361-M001	1mg
ALX-550-361-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(Z)-4-[1-(p-(Dimethylaminoethoxy)phenyl)-2-phenyl-1-butenyl]phenol
Appearance	White to off-white powder.
CAS	68047-06-3, 65213-48-1
Formula	$C_{26}H_{29}NO_2$
Identity	Determined by MS.
MW	387.5
Purity	≥98% (UHPLC)
Solubility	Soluble in 100% ethanol (20mg/ml) or methanol (10mg/ml); insoluble water.

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