

BADGE

PPAR γ inhibitor

Antagonist of PPAR γ . Binds to PPAR γ with an apparent K_d of 100 μ M.

Interferes with the ability of 3T3-L1 and 3T3-F442A cells to undergo hormone-mediated cell differentiation.

Citations: 5

[View Online »](#)

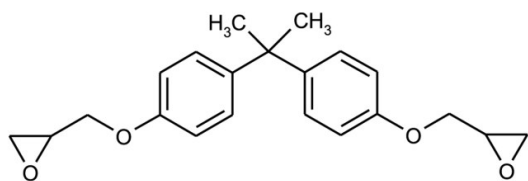
Ordering Information

[Order Online »](#)

ALX-270-353-G025	25g
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in 100% ethanol, DMSO or dimethyl formamide are stable for at least 1 year if stored at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Bisphenol A diglycidyl ether, 2,2'-[(1-Methylethylidene)bis(4,1-phenyleneoxymethylene)]bis-oxirane
Appearance	Colorless viscous liquid.
CAS	1675-54-3
Couple Target	PPAR
Couple Type	Inhibitor
Formula	$C_{21}H_{24}O_4$
Identity	Determined by IR.
MW	340.4
MeltingPoint	40-44°C
Purity	≥85%
Solubility	Miscible in organic solvents as 100% ethanol, DMSO or dimethyl formamide; insoluble in PBS, pH 7.2.