

T0901317

LXR activator

Potent agonist ($EC_{50}=50\text{nM}$) of the orphan nuclear hormone receptor (liver X receptor; LXR). Inhibitor of human sex hormone-dependent cancer cell lines.

Citations: 9

[View Online »](#)

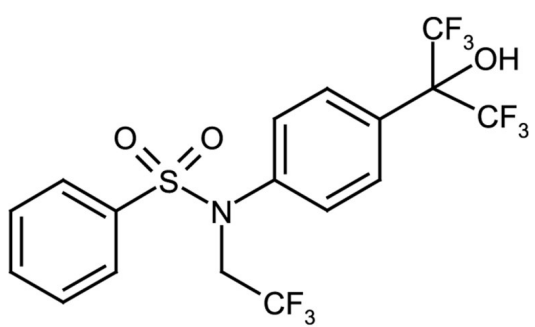
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-270-309-M010 | 10mg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|---|
| Use/Stability | As indicated on product label or CoA when stored as recommended. We do not recommend storing the aqueous solutions for more than one day. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------|--|
| Alternative Name | N-(2,2,2-Trifluoroethyl)-N-[4-[2,2,2-trifluoro-1-hydroxy-1(trifluoromethyl)ethyl]phenyl]-benzenesulfonamide |
| Appearance | White to off-white crystalline solid. |
| CAS | 293754-55-9 |
| Couple Target | LXR |
| Couple Type | Activator, Ligand |
| Formula | $C_{17}H_{12}NF_9O_3S$ |
| Identity | Determined by MS. |
| MW | 481.3 |
| Purity | ≥98% (HPLC) |
| Solubility | 50mg/ml soluble in 100% ethanol, DMSO, dimethyl formamide or any organic solvent; also soluble in a 1:1 solution of DMSO:PBS, pH 7.2 (500µg/ml). |

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-be@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