

SR-12813

PXR activator

Potent and selective PXR agonist. (EC_{50} =200 nM for human PXR). A weak activator of rat and mouse PXR. Inhibits cholesterol biosynthesis in a number of species.

Citations: 9

[View Online »](#)

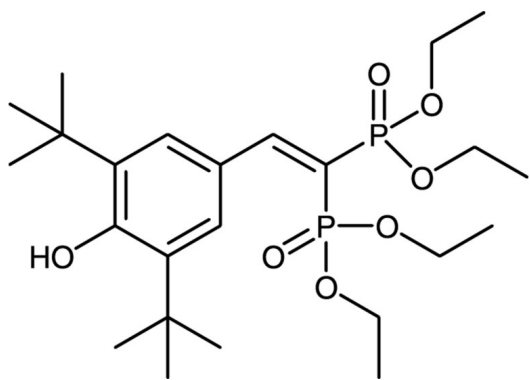
Ordering Information

[Order Online »](#)

BML-GR242-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tetraethyl 2-(3,5-Di-t-butyl-4-hydroxyphenyl)ethenyl-1,1-bisphosphonate

Appearance White solid.

CAS 126411-39-0

Couple Target PXR

Couple Type Activator, Ligand

Formula $C_{24}H_{42}O_7P_2$

MW 504.5

Purity $\geq 98\%$ (TLC)

Solubility Soluble in DMSO (25mg/ml).

Source Synthetic.

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