

# GW501516

PPAR $\delta$  agonist

GW501516 is a specific agonist for PPAR $\delta$  ( $EC_{50}$ =1.1nM) with a 1'000-fold selectivity over the other human subtypes.

Citations: 44

[View Online »](#)

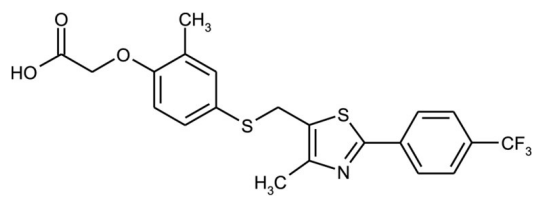
Ordering Information

[Order Online »](#)

ALX-420-032-M001	1mg
ALX-420-032-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

Appearance	White to off-white solid.
CAS	317318-70-0
Couple Target	PPAR
Couple Type	Activator, Ligand
Formula	$C_{21}H_{18}F_3NO_3S_2$
Identity	Identity determined by 1H-NMR and 13C-NMR.
MW	453.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (50mg/ml), dimethyl formamide (25mg/ml), 100% ethanol (12mg/ml) or methanol (50mg/ml); slightly soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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