

# SEW-2871

## S1P receptor ligand

SEW-2871 is a novel, potent EDG receptor agonist structurally unrelated to sphingosine-1-phosphate and is highly selective for the human S1P1 receptor. It inhibits lymphocyte recirculation *in vivo* and produces immunosuppression without associated bradycardia which is induced by other less selective S1P agonists. In mice SEW-2871 blocked TNF $\alpha$  induced monocyte adhesion to endothelium.

Citations: 5

[View Online »](#)

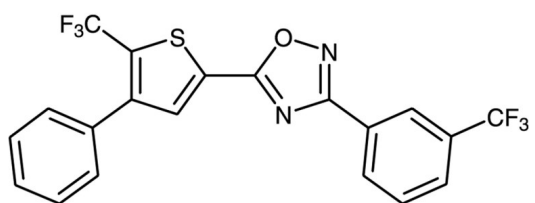
## Ordering Information

[Order Online »](#)

BML-LP106-0050	50mg
----------------	------

## 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** 5-[4-Phenyl-5-(trifluoromethyl)thiophen-2-yl]-3-[3-(trifluoromethyl)phenyl]-1,2,4-oxadiazole

**Appearance** White solid.

**CAS** 256414-75-2

**Couple Target** S1P receptor

**Couple Type** Ligand

**Formula**  $C_{20}H_{10}F_6N_2OS$

**MW** 440.4

**Purity**  $\geq 98\%$

**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-  
fr@enzolifesciences.com](mailto:info-fr@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)