

SEW-2871

S1P receptor ligand

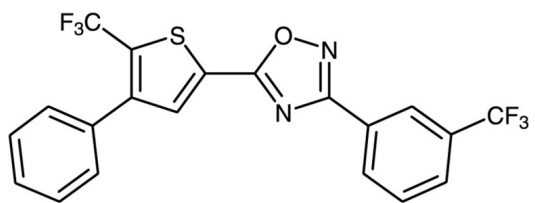
SEW-2871 is a novel, potent EDG receptor agonist structurally unrelated to sphingosine-1-phosphate and is highly selective for the human S1P1 receptor. It inhibits lymphocyte recirculation *in vivo* and produces immunosuppression without associated bradycardia which is induced by other less selective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

Citations: 5

[View Online »](#)

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

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