

# S1P1 (human) polyclonal antibody

Sphingosine-1-phosphate (S1P), a bioactive lipid, stimulates signaling pathways via G-protein-coupled receptors such as S1P1.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                |       |
|----------------|-------|
| BML-SA557-0100 | 100µg |
|----------------|-------|

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

|                      |                                                                        |
|----------------------|------------------------------------------------------------------------|
| Alternative Name     | Sphingosine 1-phosphate receptor 1, EDG-1 receptor                     |
| Application          | ICC, IP, WB                                                            |
| Host                 | Rabbit                                                                 |
| Immunogen            | Synthetic peptide corresponding to aa 359-372 of mouse S1P1.           |
| Species Reactivity   | Human                                                                  |
| UniProt ID           | O08530                                                                 |
| Worry-free Guarantee | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> . |

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)