

# PD-L1 recombinant monoclonal antibody (485)

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

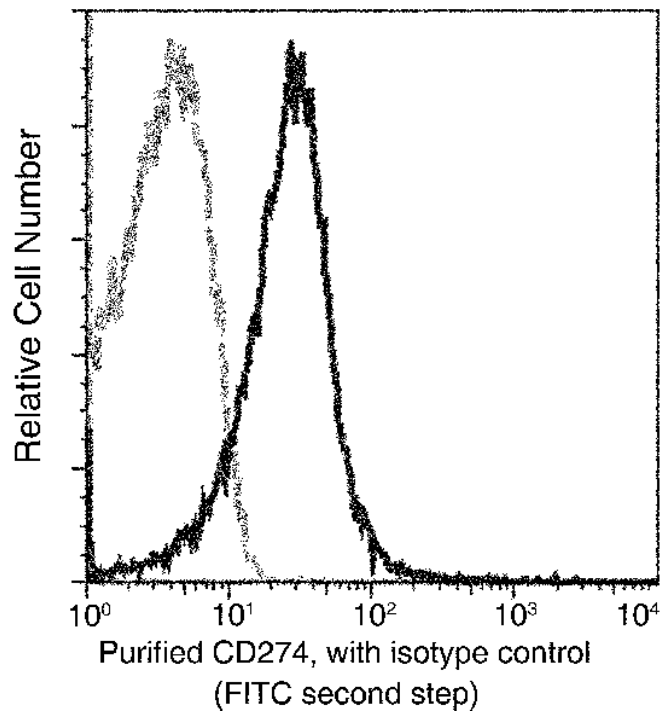
## Ordering Information

[Order Online »](#)

ENZ-ABS1003-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse CD274 expression on BABL/c splenocytes. Cells were stained with purified anti-Mouse CD274, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

## Handling & Storage

**Handling** Protected from light. Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Programmed cell death 1 ligand 1, CD274, B7-H1, Programmed death ligand 1

**Application** Flow Cytometry

**Clone** 485

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free.

**Host** Rabbit

**Immunogen** Recombinant mouse PD-L1 / B7-H1 / CD274 protein.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (1:25-1:100) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Mouse

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



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)