

PD-L1 recombinant monoclonal antibody (485)

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

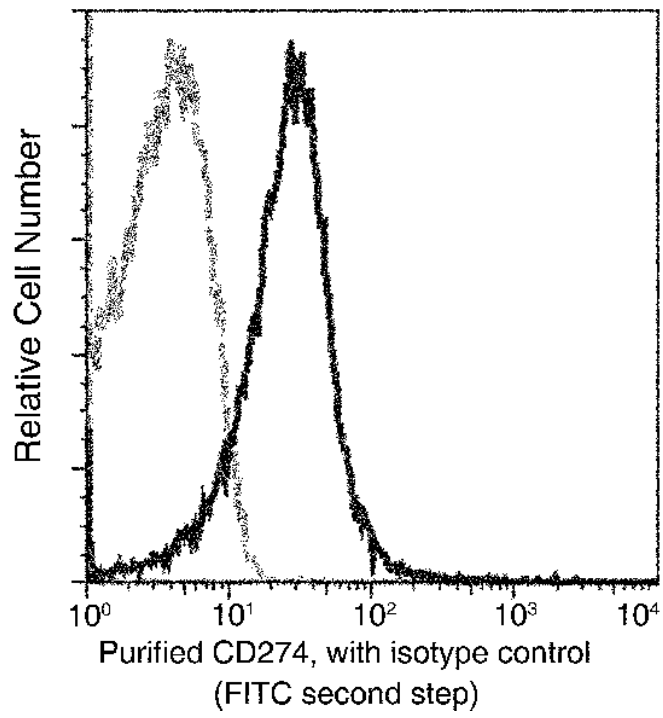
Ordering Information

[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



Last modified: September 17, 2024

ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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