

# PD-1 monoclonal antibody (4F12)

PDCD1 is an immunoreceptor that belongs to the immunoglobulin (Ig) superfamily, also known as Programmed Cell Death 1. It is expressed on activated T cell, B cells, myeloid cells and a subset of thymocytes. PDCD1 interacts with expressed PDL1 and PDL2 to inhibit the proliferation and activation of T cells and B cells. Both PDL1 and PDL2 are expressed in a wide variety of normal tissues including heart, lung, placenta, spleen and kidney.

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

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Programmed cell death 1, CD279, Programmed death receptor 1
<b>Application</b>	ELISA, Flow Cytometry, ICC, IF, WB
<b>Clone</b>	4F12
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human PDCD1 (aa 21-167) purified from E. coli.
<b>Isotype</b>	IgG2bk
<b>Purity Detail</b>	Protein G affinity purified
<b>Recommendation Dilutions/Conditions</b>	Western Blot (1:500).Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from ascites.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q15116
<b>Worry-free Guarantee</b>	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)