

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](#).

Ordering Information

[Order Online »](#)

ENZ-ABS737-0100	100µl
-----------------	-------

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

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