## PD-1 Recombinant monoclonal antibody (RMPD1-1)

PDCD1 (Programmed cell death-1 protein), also called PD1 and CD279, is a type I transmembrane receptor and a member of the immunoglobin gene superfamily. It is expressed on activated T-cells, B-cells, and myeloid cells. Anti-PDCD1 is a marker of angioimmunoblastic lymphoma and suggests a unique cell of origin for this neoplasm. Unlike CD10 and BCL6, PDCD1 is expressed by few B-cells, so may be a more specific and useful diagnostic marker in angioimmunoblastic lymphoma.

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

**ENZ-ABS337-0100** 100μg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Programmend cell death-1, CD279, Programmed death

receptor 1

**Application** IF, IHC (PS)

Clone RMPD1-1

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Recombinant full length human PD-1.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunofluorescence (0.5-1 µg/mL)Immunohistochemistry

(formalin-fixed paraffin sections) (0.5-1  $\mu$ g/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

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 infoLast modified: Musy 29 1/2 2024 com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com