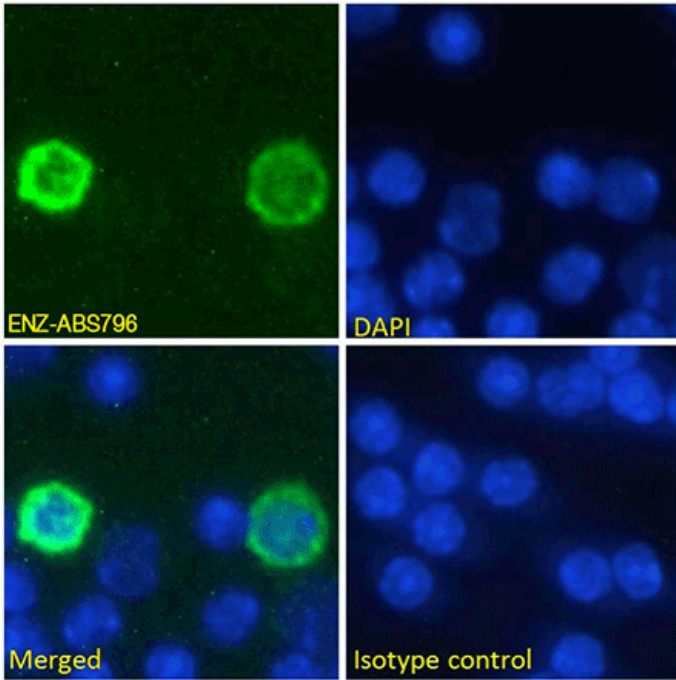


PD-1 Recombinant monoclonal antibody (RMP1-14)

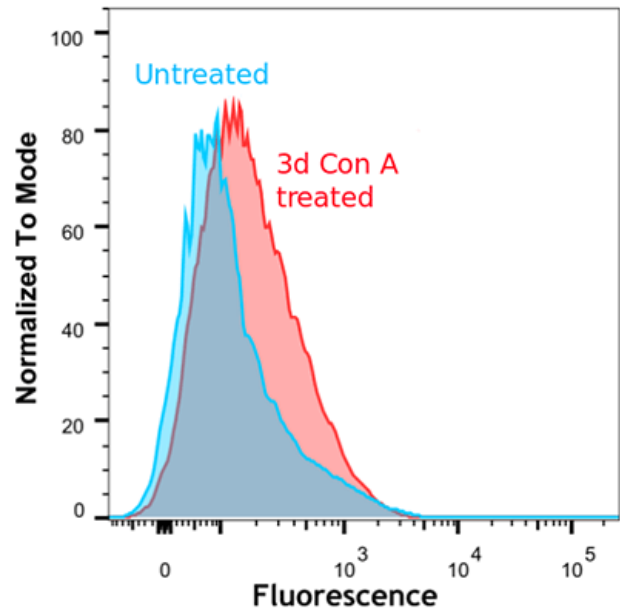
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of fixed mouse splenocytes with PD-1 (programmed cell death protein 1) antibody ENZ-ABS796. Immunofluorescence analysis of paraformaldehyde fixed mouse (*Mus musculus*) splenocytes. The nuclear stain is DAPI (blue). Panels show from left-right, top-bottom ENZ-ABS796, DAPI, merged channels and an isotype control.



Flow-cytometry using ENZ-ABS796. Staining of murine splenocytes (unstimulated and 3-day Con A stimulated) at a concentration of 1 $\mu\text{g/ml}$ for 30 mins at RT.

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Programmed death receptor 1, Programmend cell death-1, CD279
Application	Flow Cytometry, IF
Clone	RMP1-14
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	The original rat hybridoma was generated by immunizing Sprague Dawley rats with mouse PD-1-transfected BHK cells. A P3U1 myeloma was used as the fusion partner.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Source	This chimeric rabbit antibody was made using the variable domain sequences of the original Rat IgG2a format, for improved compatibility with existing reagents, assays and techniques.
Species Reactivity	Mouse
UniProt ID	Q02242
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

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)