

# PD-L1 Recombinant monoclonal antibody (YDC 127.1.1)

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

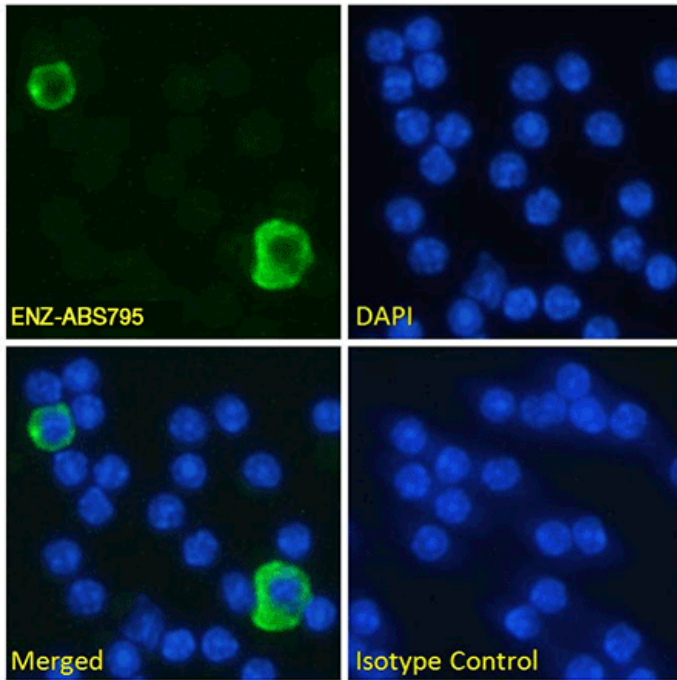
## Ordering Information

[Order Online »](#)

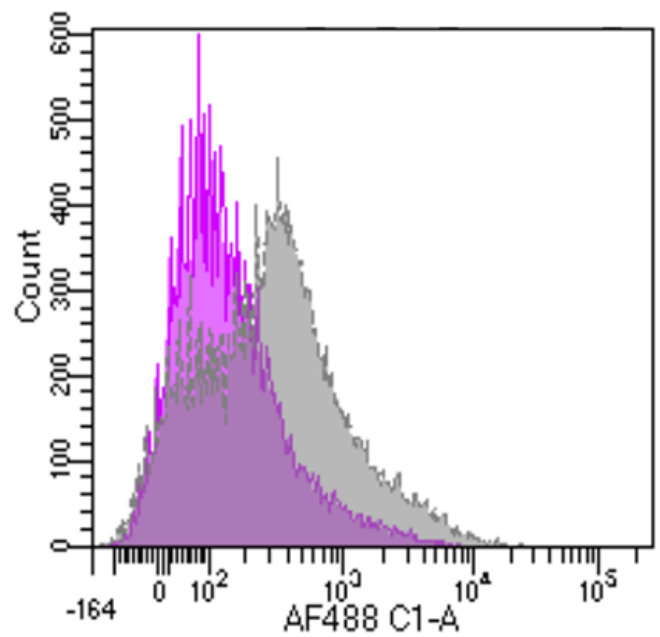
ENZ-ABS795-0200	200µg
-----------------	-------

## Manuals, SDS & CofA

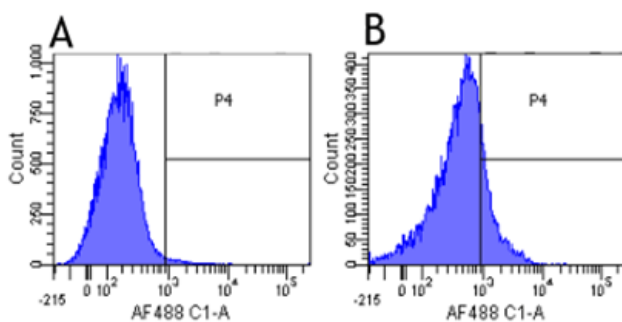
[View Online »](#)



Immunofluorescence staining of fixed mouse splenocytes with ENZ-ABS795. Immunofluorescence analysis of paraformaldehyde fixed mouse (*Mus musculus*) splenocytes. The isotype control was stained with an anti-Fluorescein antibody followed by Alexa Fluor® 488 secondary antibody.



Flow-cytometry using ENZ-ABS795. Mouse splenocytes were stained at a concentration of 1  $\mu\text{g/ml}$ .



Flow-cytometry using ENZ-ABS795. Mouse lymphocytes were stained at a concentration of 1  $\mu\text{g/ml}$ .

## 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 cell death 1 ligand 1, CD274, B7-H1, Programmed death ligand 1
Application	Flow Cytometry, IF
Clone	YDC 127.1.1
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	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.
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	Q9EP73



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 1 72 440 1335  
[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)

This antibody is covered by our **Worry-Free Guarantee**