PD-L1 Recombinant monoclonal antibody (YDC 127.1.1)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

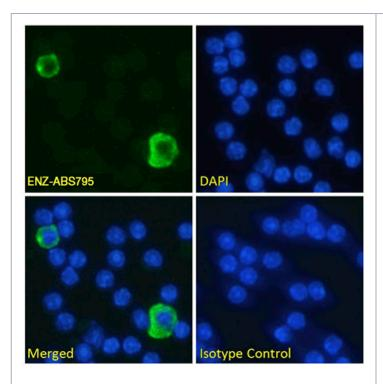
Order Online »

ENZ-ABS795-0200

200µg

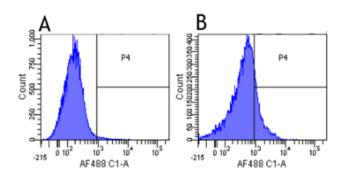
Manuals, SDS & CofA

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 spleenocytes were stained at a concentration of 1 µg/ml.



Flow-cytometry using ENZ-ABS795. Mouse lymphocytes were stained at a concentration of 1 µg/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

ImmunogenThis 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com