PD-L1 monoclonal antibody

Monoclonal antibody to PD-L1 for use with IHC applications

PD-L1, also known as B7-H1 and CD274, is a 290 amino acid transmembrane glycoprotein in the B7 family of immune regulatory molecules. It has an extracellular domain that can interact with PD-1, which belongs to the CD28/CTLA4 family expressed on activated lymphoid cells. The PD1/PD-L1 interaction ensures that activation of the immune system occurs at the appropriate time. This will minimize the possibility of chronic autoimmune inflammation. PD-L1 is known to inhibit proliferation of T cells via apoptosis. It also plays an important role in arresting cell-cycle progression. This makes PD-L1 an attractive therapeutic target for cancer and immunologic diseases.

This antibody is covered by our Worry-Free Guarantee.

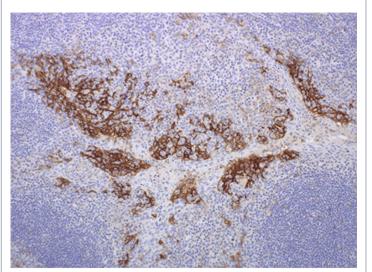
Ordering Information

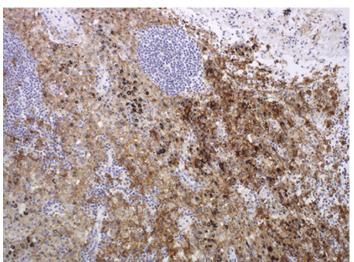
Order Online »

ENZ-ABS693-0100 100μl

Manuals, SDS & CofA

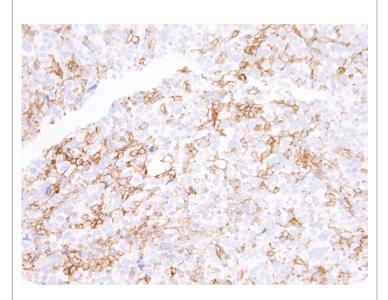
View Online »





Formalin-fixed paraffin-embedded tonsil tissue stained using Enzo's PD-L1 antibody (ENZ-ABS693). IHC staining was performed on an automated Leica BOND III.

Formalin-fixed paraffin-embedded lung tissue stained using Enzo's PD-L1 antibody (ENZ-ABS693). IHC staining was performed on an automated Leica BOND III.



IHC using anti-PD-L1 (Prod. No. ENZ-ABS693) on lung adenocarcinoma tissue.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage +4°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 IHC, WB

Application Notes For IHC applications: Dilute the concentrate as

appropriate in antibody diluent, and then apply 150 µL to

the slide, as indicated in the user's IHC protocol.

Formulation Liquid. In Tris buffer, pH 7.5, containing 1% BSA and

<0.1% sodium azide.

Host Rabbit

Immunogen Recombinant PD-L1.

Isotype IgG

Purity Detail Affinity-purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:50-1:100)Optimal conditions

must be determined individually for each application.

Species Reactivity Human

UniProt ID Q9NZQ7

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com