

PD-L1 monoclonal antibody

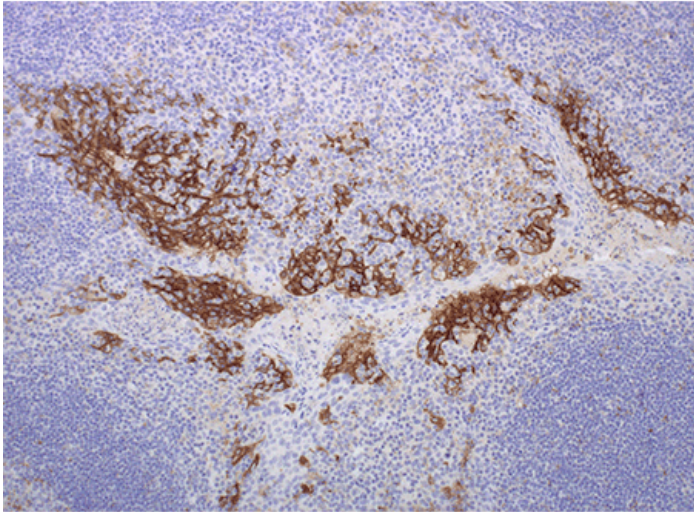
Monoclonal antibody to PD-L1 for use with IHC applications

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.

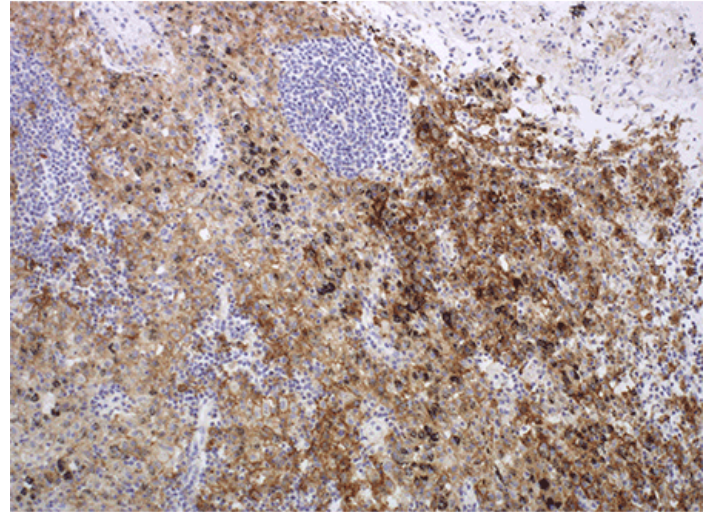
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

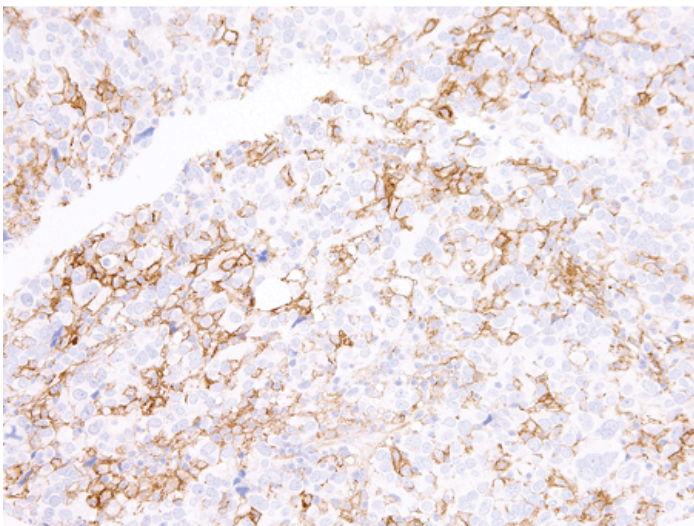
[View Online »](#)



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



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)