

HLA-DR polyclonal antibody

The anti-HLA-DR antibody detects human HLA-DR, a major histocompatibility complex (MHC) class II cell surface receptor expressed on antigen-presenting cells. This antibody is optimized for use in immunohistochemistry (IHC).

HLA-DR is a heterodimeric protein that belongs to the MHC class II family and is predominantly expressed on professional antigen-presenting cells, including dendritic cells, macrophages, and B cells. It plays a central role in adaptive immunity by presenting extracellular antigens to CD4⁺ T cells, initiating an immune response. HLA-DR is critical for immune system function and is widely studied in the context of autoimmune diseases, transplantation immunology, and infectious diseases. Dysregulation of HLA-DR expression is associated with conditions such as rheumatoid arthritis, multiple sclerosis, and chronic inflammatory disorders. Additionally, HLA-DR is often upregulated in activated microglia, making it a key marker for studying neuroinflammation in diseases like Alzheimer's and Parkinson's.

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

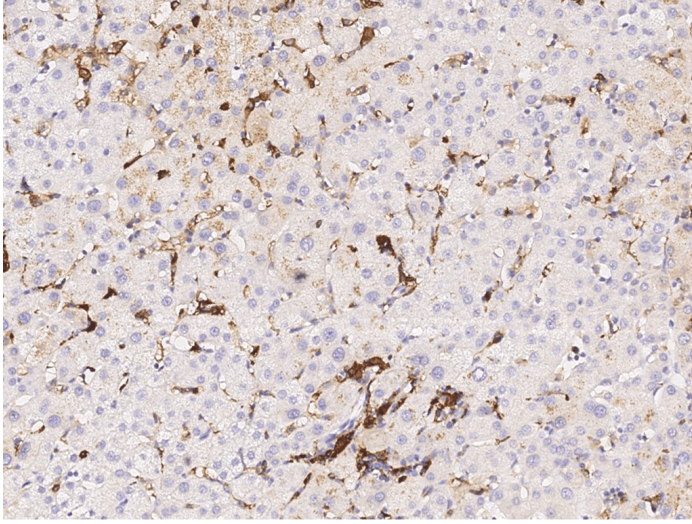
Ordering Information

[Order Online »](#)

ENZ-ABS1045-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human HLA-DR in human liver with rabbit polyclonal antibody (1:1000, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS)

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free.
Preservative-free.

Host Rabbit

Immunogen Synthetic peptide corresponding to the center region of the Human HLA-DR.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:500-1:2'000)

Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.

Species Reactivity Human

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

Last modified: January 9, 2025



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)