

# Lactate Dehydrogenase A monoclonal antibody (AT1A4)

Lactate dehydrogenase A (LDHA), which is encoded by a target gene of c-Myc and hypoxia-inducible factor (HIF-1). LDHA is the predominant form in skeletal muscle and kinetically favors the conversion of pyruvate to lactate. Therefore, LDH-A is an attractive target for cancer therapy since its expression is largely relegated to skeletal muscle.

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** L-lactate dehydrogenase A chain, LDH1

**Application** ELISA, Flow Cytometry, ICC, IF, WB

**Clone** AT1A4

**Formulation** Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.

**Host** Mouse

**Immunogen** Recombinant human LDHA (aa 1-332) purified from E. coli.

**Isotype** IgG2bk

**Purity Detail** Protein G affinity purified

**Recommendation Dilutions/Conditions** Western Blot (1:1,000). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from ascites.

**Species Reactivity** Human

**UniProt ID** P00338

**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)