

# MHC Class I monoclonal antibody (W6/32)

HLA-class I major histocompatibility (MHC) antigens are intrinsic membrane glycoproteins expressed on nucleated cells and noncovalently associated with an invariant beta2 microglobulin. They carry foreign determinants important for immune recognition by cytotoxic T cells, thus important for anti-viral and anti-tumour defence. Human HLA-class I antigens are represented by HLA-A, HLA-B and HLA-C molecules.

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

Citations: 19

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-805-711-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	HLA Class I
Application	ELISA, Flow Cytometry, FUNC, ICC, IHC (FS), IP, WB
Clone	W6/32
Crossreactivity	Does not cross-react with rabbit MHC class I.
Formulation	Liquid. In PBS, pH 7.4. 15mM sodium azide.
Host	Mouse
Immunogen	Membrane of human tonsil cells.
Isotype	IgG2a
Purity Detail	Purified by protein-A affinity chromatography.
Recommendation Dilutions/Conditions	Western Blot (non-reducing)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Bovine, Cat, Human, Monkey
Specificity	Recognizes virtually all nucleated human cells. Suitable as positive control for HLA tissue typing and crossmatching.
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

Last modified: February 25, 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-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)

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.