

MHC Class I monoclonal antibody (W6/32) (FITC conjugate)

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](#).

Ordering Information

[Order Online »](#)

ALX-805-711F-C100	FITC - 100µg
-------------------	--------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Do not freeze. Protect from light.
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	Flow Cytometry
Clone	W6/32
Formulation	Liquid. In PBS, containing 15mM sodium azide.
Host	Mouse
Immunogen	Membrane of human tonsil cells.
Isotype	IgG2a
Purity Detail	Protein G-affinity purified. The purified antibody is conjugated with fluorescein isothiocyanate (FITC) under optimum conditions. The conjugate is purified by size-exclusion chromatography.
Species Reactivity	Bovine, Cat, Human, Monkey
Specificity	The antibody W6/32 recognises an extracellular epitope of MHC Class I molecules (MHC Class Ia) that are expressed on the surface of all human nucleated cell types. This antibody W6/32 is also suitable as a positive control for HLA tissue typing and crossmatching.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)