

MHC Class I monoclonal antibody (W6/32) (R-PE 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](#).

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-711R0

R-PE - 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

Crossreactivity Does not cross-react with rabbit MHC class I.

Formulation Liquid. In PBS, pH 7.4, 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 R-Phycoerythrin (PE). The conjugate is purified by size-exclusion chromatography.

Recommendation Dilutions/Conditions Flow Cytometry (1:50) 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.

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)