

# CD1a Recombinant monoclonal antibody (C1A/1506R)

CD1a is a non-polymorphic MHC Class 1 related cell surface glycoprotein expressed in association with  $\beta$ -2 microglobulin.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS344-0100	100 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** T-cell surface glycoprotein CD1a, T-cell surface antigen T6/Leu-6

**Application** Flow Cytometry, IF, IHC (PS)

**Clone** C1A/1506R

**Formulation** Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.

**Gene/Protein Identifier** 909 (Gene ID)

**Host** Rabbit

**Immunogen** Recombinant full length human CD1a.

**Isotype** IgGk

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (0.5-1 µg/10<sup>6</sup> cells) Immunofluorescence (1-2 µg/mL) Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL for 30 min at RT) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**UniProt ID** P06126

**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-  
fr@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)