

CD1a Recombinant monoclonal antibody (C1A/1506R)

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

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

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	IgGκ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1 µg/10 ⁶ 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

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)