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.

Ordering Information

Order Online »

ENZ-ABS344-0100

100µg

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/106 cellsImmunofluorescence

(1-2 μ g/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 μ g/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



info-

eu@enzolifesciences.com