



PRODUCT DATA SHEET

CD1a Recombinant monoclonal antibody (C1A/1506R)

ENZ-ABS344

Product Number/Sizes

ENZ-ABS344-0100 100 µg

Product Details

ALTERNATIVE NAME:	T-cell surface glycoprotein CD1a, T-cell surface antigen T6/Leu-6
CLONE:	C1A/1506R
HOST:	Rabbit
ISOTYPE:	IgGκ
IMMUNOGEN:	Recombinant full length human CD1a.
UNIPROT ID:	P06126
GENE/PROTEIN IDENTIFIER:	909 (Gene ID)
SPECIES REACTIVITY:	Human
APPLICATIONS:	Flow Cytometry, IF, IHC (PS)
RECOMMENDED 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.
PURITY DETAIL:	Protein A affinity purified.
FORMULATION:	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING:	Blue Ice
LONG TERM STORAGE:	+4°C
SCIENTIFIC BACKGROUND:	CD1a is a non-polymorphic MHC Class 1 related cell surface glycoprotein expressed in association with β-2 microglobulin.
REGULATORY STATUS:	RUO - Research Use Only

Revised 19-Dec-23

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
F +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com