



# PRODUCT DATA SHEET

## CD8 (α chain) Recombinant monoclonal antibody (C8/1779R)

ENZ-ABS323

### Product Number/Sizes

ENZ-ABS323-0100 100 µg

### Product Details

ALTERNATIVE NAME:	T-cell surface glycoprotein CD8 (α chain), T-lymphocytid differentiation antigen T8/Leu-2
CLONE:	C8/1779R
HOST:	Rabbit
ISOTYPE:	IgGκ
IMMUNOGEN:	Recombinant full length human CD8a.
UNIPROT ID:	P01732
SPECIES REACTIVITY:	Human
APPLICATIONS:	Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS:	Flow Cytometry (0.5-1 µg/10 <sup>6</sup> cells) Immunofluorescence (0.5-1 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 G 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:	CD8 is a cell surface receptor expressed either as a heterodimer with the CD8 β chain (CD8 α/β) or as a homodimer (CD8 α/α). A majority of thymocytes and a subpopulation of mature T cells and NK cells express CD8a. CD8 binds to MHC class 1 and through its association with protein tyrosine kinase p56lck plays a role in T cell development and activation of mature T cells. For mature T-cells, CD4 and CD8 are mutually exclusive, so anti-CD8, generally used in conjunction with anti-CD4. It is a useful marker for distinguishing helper/inducer T-lymphocytes, and most peripheral T-cell lymphomas are CD4+/CD8-. Anaplastic large cell lymphoma is usually CD4+ and CD8-, and in T-lymphoblastic lymphoma/leukemia, CD4 and CD8 are often co-expressed. CD8 is also found in littoral cell angioma of the spleen.
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](mailto:info-usa@enzolifesciences.com)  
[www.enzolifesciences.com](http://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](mailto:info-ch@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)