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

CD8 is a cell surface receptor expressed either as a heterodimer with the CD8  $\beta$  chain (CD8  $\alpha/\beta$ ) or as a homodimer (CD8  $\alpha/\alpha$ ). 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.

This antibody is covered by our Worry-Free Guarantee.

# **Ordering Information**

**Order Online** »

ENZ-ABS323-0100	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 CD8 (α chain), T-lymphocyted

differentiation antigen T8/Leu-2

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

**Clone** C8/1779R

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Recombinant full length human CD8a.

Isotype IgGk

Purity Detail Protein G affinity purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (0.5-1 μg/106 cellsImmunofluorescence

(0.5-1  $\mu$ g/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 P01732

### **Worry-free Guarantee**

ENZO LIFE SCIENCES, INC.
Phone: 800.942.0430 info-usa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024