

# CD99 Recombinant monoclonal antibody (MIC2/1495R)

CD99 is a cell surface glycoprotein involved in leukocyte migration, T-cell adhesion, ganglioside GM1 and transmembrane protein transport, and T-cell death by a caspase-independent pathway. In addition, it may have the ability to rearrange the actin cytoskeleton and may also act as an oncosuppressor in osteosarcoma. Cyclophilin A binds to CD99 and may act as a signaling regulator of the protein.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS357-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	Flow Cytometry, IF, IHC (PS)
Clone	MIC2/1495R
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 CD99.
Isotype	IgGκ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (2-4 µg/106 cells in 0.1 mL)Immunofluorescence (2-4 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (5-10 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	P14209
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>