

CD38 Recombinant monoclonal antibody (HB7) (Rabbit IgGκ)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	ADP-ribosyl cyclase 1, Cyclic ADP-ribose hydrolase, Cyclic ADP-ribose hydrolase 1, cADPr hydrolase 1, T10, p45, NAD(+) nucleosidase
Application	Flow Cytometry, WB
Clone	HB7
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	Viable cells (1 x 10E7) of the human cell line, BJAB (Human Burkitt lymphoma B cell line), which were injected intraperitoneally into 8-week-old BALB/c mice.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Human

Specificity

Recognizes an epitope between aa 273-285 of human CD38, an approximately 45 kDa type II transmembrane protein, expressed on essentially all pre-B lymphocytes, plasma cells, and thymocytes. Also present on activated T lymphocytes, natural killer (NK) lymphocytes, myeloblasts, and erythroblasts. Bimodally expressed during B cell development, modulating from high in immature cells to low in intermediate ones and back to high on mature B cells. This antibody competes with clone AT13/5

Technical Info / Product Notes

This chimeric rabbit antibody was made using the variable domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID

P28907

Worry-free Guarantee

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

Last modified: May 29, 2024



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
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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