

CD38 Recombinant monoclonal antibody (HB7) (Mouse IgG1κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS438-0200	200µg
-----------------	-------

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	Mouse
Immunogen	Viable cells (1 x 10E7) of the human cell line, BJAB (Human Burkitt lymphoma B cell line), which was were injected intraperitoneally into 8-week-old BALB/c mice.
Isotype	IgG1κ
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
UniProt ID	P28907

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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