# CD38 Recombinant monoclonal antibody (HB7) (Mouse IgG1k)

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**

Isotype

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

lgG1κ

(Human Burkitt lymphoma B cell line), which was were injected intraperitoneally into 8-week-old BALB/c mice.

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

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France Phone: +33 472 440 655 fr@enzolifesciences.com Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900

uk@enzolifesciences.com