

## CD38 Monoclonal Mouse Antibody (FS02)

### Product Description

This antibody reacts with a type II membrane glycopeptide with a transmembrane sequence near the NH<sub>2</sub>-terminus. CD38 is a type II transmembrane glycoprotein that is present on early B- and T-cell lineages and activated B- and T-cells but is absent from most mature resting peripheral lymphocytes. CD38 is also found on thymocytes, pre-B cells, germinal center B-cells, mitogen-activated T-cells, monocytes and Ig-secreting plasma cells. CD38 is expressed on CD34<sup>+</sup> cells. The CD34<sup>+</sup>CD38<sup>-</sup> population of hematopoietic stems cells defines the most pluripotent cells (e.g. blast colony forming cells).

### Product attributes

Antibody number	#0171
Antibody reactivity (target)	CD38
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FS02
Isotype	IgG1, kappa
Molecular weight	~45 kDa (Glycoprotein); 35 kDa (protein core)
Synonyms	Acute Lymphoblastic Leukemia Cells Antigen; ADP Ribosyl Cyclase 1; cADP-ribose Hydrolase 1; CD38H; NAD(+) Nucleosidase; NIM-R5 Antigen; p45; T10
Human gene symbol	CD38
Entrez gene ID	952
SwissProt	P28907
Unigene	479214
Immunogen	Human CD38
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	CCRF-CEM cells, Tonsil, Spleen or Skin
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	B-cells, T-cells
Antibody research areas	Immunology