

CD38 Monoclonal Mouse Antibody (FS02)



Product Description

This antibody reacts with a type II membrane glycopeptide with a transmembrane sequence near the NH₂-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+ cells. The CD34+CD38- population of hematopoietic stems cells defines the most pluripotent cells (e.g. blast colony forming cells).

Product attributes

Antibody number	#0171
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
Cellular localization	Cytoplasmic, Membrane/cell surface, Nucleus
Species reactivity	Human
Applications	Immunofluorescence, Flow cytometry
Application notes	Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user
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)
Supplied As	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, BSA-free: 1 mg/mL in PBS without azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% 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