



# CD100 Monoclonal Mouse Antibody (SEMA4D/933)

## Product Description

CD100 (also known as Semaphorin 4D), is a homodimeric protein comprised of 50 kDa subunits. It is expressed on resting T cells and weakly on other lymphocytes, and its expression is increased on B and T cells after stimulation. CD100 is a cell surface receptor that binds to CD72, Plexin B1 and Plexin B2. It promotes B cell aggregation, cell survival, and cell proliferation.

### Product attributes

Antibody number	#0933
Antibody reactivity (target)	CD100, Semaphorin-4D
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SEMA4D/933
Isotype	IgG1, kappa
Molecular weight	50 kDa (Monomer)
Synonyms	CD100; Semaphorin-4D; SEMA4D
Human gene symbol	SEMA4D
Entrez gene ID	10507
SwissProt	Q92854
Unigene	494406
Immunogen	Human recombinant SEMA4D protein
Antibody target cellular localization	Plasma membrane
Species reactivity	Human, Non-human primate
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	Daudi, Raji, HUT-78, Kg1a, U937, and human lymphocytes. Human tonsils and lymph nodes.
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	Hematopoietic cells
Antibody research areas	Immunology, Neuroscience, Signal transduction