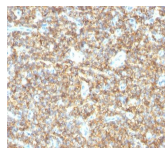


## CD99 Monoclonal Mouse Antibody (HO36-1.1)



### Product Description

Recognizes a sialoglycoprotein of 27-32 kDa, identified as CD99, or MIC2 gene product, or E2 antigen. MIC2 gene is located in the pseudo-autosomal region of the human X and Y chromosome. MIC2 gene encodes two distinct proteins, which are produced by alternative splicing of the CD99 gene transcript and are identified as bands of 30 and 32 kDa (p30/32). Although its function is not fully understood, CD99 is implicated in various cellular processes including homotypic aggregation of T cells, upregulation of T cell receptor and MHS molecules, apoptosis of immature thymocytes and leukocyte diapedesis. CD99 is expressed on the cell membrane of some lymphocytes, cortical thymocytes, and granulosa cells of the ovary. Most pancreatic islet cells, Sertoli cells of the testis, and some endothelial cells express this antigen. Mature granulocytes express very little or no CD99. MIC2 is strongly expressed on Ewing's sarcoma cells and primitive peripheral neuroectodermal tumors. This MAb shows a very similar reactivity to other CD99 MABs (e.g. O13, 12E7, or HBA-71) and is excellent for immunohistochemical staining of formalin-fixed, paraffin-embedded tissues. This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

### Product attributes

Antibody number	#0869
Antibody reactivity (target)	CD99
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HO36-1.1
Isotype	IgM, kappa
Molecular weight	27-32 kDa
Synonyms	12E7; E2 antigen; MIC 2X; MIC 2Y; MIC2; Protein MIC2; Surface antigen MIC2; T-cell surface glycoprotein E2
Human gene symbol	CD99
Entrez gene ID	4267
SwissProt	P14209
Unigene	653349
Immunogen	Purified E-rosette forming cells from human peripheral blood lymphocytes
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human, Rat
Antibody target cellular localization	Plasma membrane
Antibody application notes	Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 1-2 ug/mL, Flow Cytometry 5-10 uL/million cells in 0.1 mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody research areas	Cancer, Endocrinology, Immunology
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
Positive control	Pancreas or Ewing's sarcoma
Cell/tissue expression	Lymphocytes, Ovary, Pancreatic islet cells, T-cells, Testis
Tumor expression	Sarcoma