## CD99 / MIC2 Monoclonal Mouse Antibody (12E7 + MIC2/877)

## **Product Description**

This antibody 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. Catalog number key for antibody number 0878, Anti-CD99|MIC2 (12E7 MIC2/877)

## Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#0878		
Antibody reactivity (target)	CD99, MIC2		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	12E7 + MIC2/877		
Isotype	IgG1, 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	Human acute lymphocytic leukemia T-cells (12E7); Recombinant human MIC2 protein (MIC2/877)		
Antibody target cellular localization	Plasma membrane		
Verified antibody applications	IHC (FFPE) (verified)		
Species reactivity	Human		
Antibody application notes	Flow cytometry: 5-10 uL/million cells, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-lixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user		
Antibody application notes  Positive control	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be		
	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunchistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user		
Positive control	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.		
Positive control Shipping condition	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Positive control Shipping condition Storage Conditions	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT., Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 ° C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 ° C  For research use only (RUO)  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/Purified; 0.5 mg/mL in PBS/0		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  For research use only (RUO)  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/ree: 1 mg/mL in PBS without azide  Guaranteed for at least 24 months from date of receipt when		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life	be required for direct detection using primary antibody. Initial projugates than for indirect detection with secondary antibody, Inmunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Inmunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  For research use only (RUO)  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 without azide  Guaranteed for at least 24 months from date of receipt when stored as recommended		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life Cell/tissue expression	be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistochemistry (formalin-fixed): 1:100-1:200 for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Immunofluorescence 1:100-1:200, Optimal dilution for a specific application should be determined by user  MOLT-4 cells. Pancreas or Ewing's sarcoma.  Room temperature  Store at 2 to 8 ° C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 ° C  For research use only (RUO)  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 surface as recommended  Guaranteed for at least 24 months from date of receipt when stored as recommended  Lymphocytes, Ovary, Pancreatic islet cells, T-cells, Testis  Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in		

Email: techsupport@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
Antibody # prefix	Conjugation	EX/EIII (IIIII)	Laser IIIIe	Detection chainlei	Dye realules
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on October 29, 2025 at 09:37:27 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cd99-mic2-monoclonal-mouse-antibody-12e7-mic2-877/">https://biotium.com/product/cd99-mic2-monoclonal-mouse-antibody-12e7-mic2-877/</a>