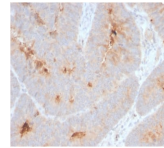


IgA Secretory Component Recombinant Monoclonal Rabbit Antibody (ECM1/2889R)



Product Description

This MAb reacts with a reduction-resistant epitope present in both free and SIgA bound Secretory Component. It does not react with the cell lines lacking secretory component. The antibody is useful for studying the distribution and level of both free and bound secretory component. Secretory component is differentially expressed in epithelium, and the antibody is a popular marker for identifying subpopulations of epithelial cells and epithelial differentiation. The Secretory component antibody is a useful research tool for studying mucosal immunity, inflammation, remodeling, differentiation and tumorigenesis, all processes associated with differential secretory component expression.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2889, Anti-Secretory Component / ECM1 (ECM1/2889R)

Product attributes

Antibody number	#2889
Antibody reactivity (target)	ECM1, Secretory Component
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	ECM1/2889R
Isotype	IgG
Molecular weight	~80 kDa
Synonyms	ECM1, Extracellular Matrix Protein 1, Secretory Component p85, URBWD
Human gene symbol	ECM1
Entrez gene ID	1893
SwissProt	Q16610
Unigene	81071
Immunogen	Recombinant full-length human ECM1 protein
Antibody target cellular	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human, Rat
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistology (formalin): 0.5-1 ug/mL 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, Optimal dilution for a specific application should be determined by user
Positive control	Stomach, Lung, or breast tumor
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Immunology
Cell/tissue expression	Epithelial cells

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
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
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, of LI-COR Bioscience.