

Mucin 5AC / MUC5AC Monoclonal Mouse Antibody (1-13M1)

Product Description

This mucin is present in primary ovarian mucinous cancer but usually absent in colorectal adenocarcinoma, thus showing an expression pattern opposite to MUC2. Together with a panel of antibodies, anti-MUC5AC may be useful for differential identification of primary mucinous ovarian tumors from colon adenocarcinoma metastatic to the ovary. MUC5AC antibodies may also be useful for identification of intestinal metaplasia as well as in the identification of pancreatic carcinoma and pre-cancerous changes vs. normal pancreas.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Am J Respir Crit Care Med (2001) 163: 517-523. (IHC, FFPE)
2. Oncotarget (2015) 6(6): 4274-4285. (WB)
3. Int J Cancer (1998) 75:767-773. (RIA)
4. Exp Eye Res (2007) 84(5): 939-950. (WB)
5. Am J Gastroenterol (2017) 112(1): 172-183. (ELISA)

Product attributes

Antibody number	#0432
Antibody reactivity (target)	MUC5AC
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	1-13M1
Isotype	IgG1, kappa
Molecular weight	>1,000 kDa
Synonyms	Apomucin Major Airway Glycoprotein, Mucin 5 subtype AC tracheobronchial, Mucin 5 Subtypes A And C, Mucin 5AC oligomeric mucus/gel forming, Tracheobronchial Mucin (TBM)
Human gene symbol	MUC5AC
Entrez gene ID	4586
SwissProt	P98088
Unigene	534332
Immunogen	M1 mucin preparation from the fluid of an ovarian mucinous cyst belonging to an O Le(a-b) patient
Antibody target cellular localization	Secreted (extracellular)
Expected antibody applications	RIA (published for clone), IHC (FFPE) (published for clone), WB (published for clone)
Species reactivity	Cat, Chicken, Cow, Human, Monkey, Mouse, Rabbit, Rat
Antibody application notes	Does not react with Pig or Hedgehog, ELISA: For coating, purchase antibody without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution for a specific application should be determined by user
Positive control	Human colon or stomach (IHC).
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	Epithelial cells
Tumor expression	Ovarian cancer
Antibody research areas	Cancer, Mucins