



Glowing products for science

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. **Catalog number key for antibody number 0432, Anti-MUC5AC (1-13M1)**

Call us : [800-304-5357](tel:800-304-5357)

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 of antigen | >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 |
| Product origin | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Tumor expression | Ovarian cancer |
| Antibody research areas | Cancer, Mucins |

| 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 |
| 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.

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)

This datasheet was generated on June 4, 2026 at 08:25:50 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mucin-5ac-muc5ac-monoclonal-mouse-antibody-1-13m1/>

