TAG-72 / CA72.4 Recombinant Monoclonal Mouse Antibody (rB72.3)



Product Description

Recognizes an oncofetal antigen of 220 kDa, identified as a tumor-associated glycoprotein (TAG-72) with properties of a mucin. This MAb defines the mucin-carried sialylated-Tn epitope. TAG-72 is usually expressed by adenocarcinomas, but is negative in mesotheliomas. Studies have reported that this antibody has 80% sensitivity and 93% specificity for pulmonary adenocarcinoma. Therefore, TAG-72 is a useful marker to distinguish between mesothelioma and adenocarcinoma. However, false positive reactions can occur so results must be interpreted with the utmost caution. This antibody may be useful in the differentiation of non-small cell carcinomas from small cell carcinomas of the lung. The combined use of anti-TAG-72 and anti-GCDFP-15 is valuable in the diagnosis of apocrine carcinoma.

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 3177, Anti-CA72.4|TAG-72 (rB72.3)

Product attributes

| Product attributes | | | | |
|---------------------------------------|---|--|--|--|
| Antibody number | #3177 | | | |
| Antibody reactivity (target) | CA72.4, TAG-72 | | | |
| Antibody type | Primary | | | |
| Host species | Mouse | | | |
| Clonality | Recombinant Monoclonal | | | |
| Clone | rB72.3 | | | |
| Isotype | IgG1, kappa | | | |
| Molecular weight | 220 kDa | | | |
| Synonyms | CA72.4; Tumor associated glycoprotein 72 | | | |
| Human gene symbol | Not Known | | | |
| Entrez gene ID | 182875 | | | |
| SwissProt | Not Known | | | |
| Unigene | Not Known | | | |
| Immunogen | Membrane-enriched fraction of a human breast carcinoma liver metastasis | | | |
| Verified antibody applications | IHC (FFPE) (verified) | | | |
| Antibody target cellular localization | Plasma membrane | | | |
| Species reactivity | Cow, Dog, Hamster, Human, Rat | | | |
| Positive control | Jurkat cells, Breast or lung carcinoma. | | | |
| Shipping condition | Room temperature | | | |
| Storage Conditions | Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ C, Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light | | | |
| 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 | | | |
| 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. | | | |
| Cell/tissue expression | Epithelial cells | | | |
| Tumor expression | Adenocarcinoma, Lung cancer | | | |

Call us: 800-304-5357 Email: btinfo@biotium.com

| 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.

This datasheet was generated on November 4, 2025 at 02:09:09 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tag-72-ca72-4-recombinant-monoclonal-mouse-antibody-rb72-3/