Cytokeratin 6 Monoclonal Mouse Antibody (KRT6/1702)

Product Description

This MAb recognizes a protein of 56 kDa, identified as cytokeratin 6 (CK6). In humans, multiple isoforms of Cytokeratin 6 (6A-6F), encoded by several highly homologous genes, have distinct tissue expression patterns, and Cytokeratin 6A is the dominant form in epithelial tissue. The gene encoding human Cytokeratin 6A maps to chromosome 12q13, and mutations in this gene are linked to several inheritable hair and skin pathologies. Keratins 6 and 16 are expressed in keratinocytes, which are undergoing rapid turnover in the suprabasal region (also known as hyper-proliferation-related keratins). Keratin 6 is found in hair follicles, suprabasal cells of a variety of internal stratified epithelia, in epidermis, in both normal and hyper-proliferative situations. Epidermal injury results in activation of keratinocytes, which express CK6 and CK16. CK6 is strongly expressed in about 75% of head and neck squamous cell carcinomas. Expression of CK6 is particularly associated with differentiation.

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 1702, Anti-Cytokeratin 6 (KRT6/1702)

Call us: 800-304-5357

| Product attributes | | | | |
|---------------------------------------|---|--|--|--|
| Antibody number | #1702 | | | |
| Antibody reactivity (target) | Cytokeratin 6 | | | |
| Antibody type | Primary | | | |
| Host species | Mouse | | | |
| Clonality | Monoclonal | | | |
| Clone | KRT6/1702 | | | |
| Isotype | IgG2a, kappa | | | |
| Molecular weight | 56 kDa | | | |
| Synonyms | CK6A, CK6B, CK6C, CK6D, CK6E, Keratin Type II Cytoskeleta 6A, Keratin Type II Cytoskeletal 6B, Keratin Type II Cytoskeleta 6C, Keratin Type II Cytoskeletal 6D, Keratin Type II Cytoskeleta 6E, KRT6, KRT6A, KRT6B, KRT6C, KRT6D, KRT6E | | | |
| Human gene symbol | KRT6A; KRT6B; KRT6C | | | |
| Entrez gene ID | 286887-KRT6C, 3853-KRT6A, 3854-KRT6B | | | |
| SwissProt | P02538-KRT6A, P04259-KRT6B, P48668-KRT6C | | | |
| Unigene | 700779 | | | |
| Immunogen | Recombinant human full-length keratin 6 protein | | | |
| Antibody target cellular localization | Cytoskeleton | | | |
| Verified antibody applications | IHC (FFPE) (verified) | | | |
| Species reactivity | Human, Mouse | | | |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology (formalin) 0.1-0.2 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user | | | |
| Positive control | Tonsil or basal cell carcinoma | | | |
| 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 | Conjugator: 0.1 mg/ml in PRS/0.1% RSA/0.05% azido HRP | | | |

| | 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 | Cancer, Cytoskeleton | | | |
| 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 | Epidermal cells | | | |
| Tumor expression | Squamous cell carcinoma | | | |

| 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 December 14, 2025 at 04:23:13 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mouse-monoclonal-anti-cytokeratin-6-krt6170