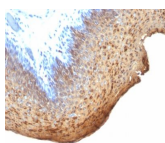


Cytokeratin 10/13 Monoclonal Mouse Antibody (DE-K13)



Product Description

This antibody recognizes cytoke­ratin 10 (56.5 kDa) and cytoke­ratin 13 (53 kDa) in Western blotting. It recognizes only cytoke­ratin 13 in formalin-fixed, paraffin-embedded tissue sections. It does not react with cytoke­ratin 10 positive, cytoke­ratin 13 negative epithelia such as epidermis. However, on frozen sections this MA­b serves as differentiation-related marker of all stratified epithelia; it stains all suprabasal cells in both cornifying and non-cornifying stratified epithelia and more differentiated cells of squamous carcinomas.

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 0696, Anti-Cytokeratin 10 (DE-K13)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #0696 |
| Antibody reactivity (target) | Cytokeratin 10/13 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | DE-K13 |
| Isotype | IgG2a, kappa |
| Molecular weight | 56.5 kDa (CK10) & 53 kDa (CK13) |
| Synonyms | BCIE, BIE, EHK, Keratin Type I Cytoskeletal 10, KRT10, CK13, K13, keratin type I cytoskeletal 13, Keratin-13, Krt1-13, KRT1 |
| Human gene symbol | KRT10 |
| Entrez gene ID | 3858 |
| SwissProt | P13645 |
| Unigene | 99936 |
| Immunogen | Cytoskeletal preparation extracted from human ectocervical epithelium |
| Antibody target cellular localization | Cytoskeleton |
| Verified antibody applications | IF (verified), IHC (FFPE) (verified) |
| Species reactivity | Cat, Human |
| 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, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Western blotting 0.25-0.5 ug/mL |
| Positive control | A431, HeLa, MCF7 cells. Tonsil, esophageal or bladder 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 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 November 24, 2024 at 06:40:14 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-cytokeratin-1013-de-k13/>