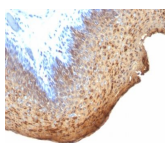


# Cytokeratin 10/13 Monoclonal Mouse Antibody (DE-K13)



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

## 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](mailto: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

<b>Antibody number</b>	#0696
<b>Antibody reactivity (target)</b>	Cytokeratin 10/13
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	DE-K13
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight of antigen</b>	56.5 kDa (CK10) & 53 kDa (CK13)
<b>Synonyms</b>	BCIE, BIE, EHK, Keratin Type I Cytoskeletal 10, KRT10, CK13, K13, keratin type I cytoskeletal 13, Keratin-13, Krt1-13, KRT1
<b>Human gene symbol</b>	KRT10
<b>Entrez gene ID</b>	3858
<b>SwissProt</b>	P13645
<b>Unigene</b>	99936
<b>Immunogen</b>	Cytoskeletal preparation extracted from human ectocervical epithelium
<b>Antibody target cellular localization</b>	Cytoskeleton
<b>Verified antibody applications</b>	IF (verified), IHC (FFPE) (verified)
<b>Species reactivity</b>	Cat, Human
<b>Antibody application notes</b>	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
<b>Positive control</b>	A431, HeLa, MCF7 cells. Tonsil, esophageal or bladder carcinoma.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer, Cytoskeleton
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	Epidermal cells
<b>Tumor expression</b>	Squamous cell carcinoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 June 16, 2026 at 01:10:10 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-cytokeratin-1013-de-k13/>