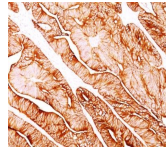


# Cytokeratin 5/8 Monoclonal Mouse Antibody (C-50)



## Product Description

This antibody reacts with keratin 5 (58 kDa) and keratin 8 (52.5 kDa). Simple epithelia express cytokeratin 8 in combination with 18. On the other hand, basal cells of stratified epithelia express cytokeratin 5 paired with 14. This antibody therefore, reacts with a wide range of epithelia and their 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 0948, Anti-Cytokeratin 5 (C-50)**

## Product attributes

<b>Antibody number</b>	#0948
<b>Antibody reactivity (target)</b>	Cytokeratin 5/8
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	C-50
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	58 kDa (CK5); 52.5 kDa (CK8)
<b>Synonyms</b>	CK5; DDD; EBS2; Epidermolysis bullosa simplex 2 Dowling-Meara/Kobner/Weber-Cockayne Types; K5; Keratin complex 2, basic, Gene 5; Keratin Type II Cytoskeletal 5; Keratin-5; KRT5; KRT5A; Type II Cytoskeletal 5; Type-II Keratin Kb5. CK8; CARD2; CYK8; CYKER; Cytokeratin Endo A; DreK8; EndoA; K2C8; K8; Keratin 8; Krt 2.8; KRT8; Type-II Keratin Kb8
<b>Human gene symbol</b>	KRT5
<b>Entrez gene ID</b>	3852 (CK5); 3856 (CK8)
<b>SwissProt</b>	P13647 (CK5); P05787 (CK8)
<b>Unigene</b>	433845 (CK5); 533782 & 708445 (CK8)
<b>Immunogen</b>	Cytoskeletal preparation from HeLa cells
<b>Antibody target cellular</b>	Cytoskeleton
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Species reactivity</b>	Cow, Dog, Ferret, Hamster, Human, Mouse, Pig, Rat, Sheep
<b>Antibody application notes</b>	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistology formalin-fixed 0.5-1.0 ug/mL. For staining of formalin-fixed tissues, digest sections with trypsin at 1 mg/mL PBS, 15 min at RT, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 0.5-1.0 ug/mL, Western blotting 0.5-1.0 ug/mL. Does not react with chicken, or Xenopus, Optimal dilution for a specific application should be determined by user
<b>Positive control</b>	HepG2 or A431 cells. Skin or Squamous cell 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>Cell/tissue expression</b>	Epithelial cells
<b>Tumor expression</b>	Breast cancer, Colorectal cancer, Lung cancer, Ovarian cancer, Thyroid cancer

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>
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, LI-COR Bioscience.

This datasheet was generated on July 22, 2024 at 12:29:13 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/monoclonal-anti-cytokeratin-58-c-50/>

