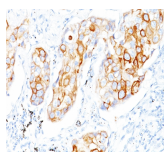


# Cytokeratin 7/17 Monoclonal Mouse Antibody (C-46)



## Product Description

Recognizes intermediate filament proteins (IFP's) of 55 kDa and 46 kDa, which are identified as cytokeratin 7 and Cytokeratin 17 respectively. On formalin-fixed, paraffin embedded sections, C-46 reacts with keratin 7 only. Cytokeratins comprise a diverse group of intermediate filament proteins (IFPs) that are expressed as pairs in both keratinized and non-keratinized epithelial tissue. Cytokeratins play a critical role in differentiation and tissue specialization and function to maintain the overall structural integrity of epithelial cells. Cytokeratins have been found to be useful markers of tissue differentiation, which is directly applicable to the characterization of malignant tumors. Cytokeratin 7 is a basic cytokeratin, which is found in most glandular and transitional epithelia but not in the stratified squamous epithelia. Cytokeratin 7 is expressed in the epithelial cells of ovary, lung, and breast but not of colon, prostate, or gastrointestinal tract. Cytokeratin 17 is expressed in epithelial cells of various origins, such as bronchial epithelial cells and skin appendages. Cytokeratin 17 antibody marks basal cell 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0949, Anti-Cytokeratin 7/17 (C-46)**

## Product attributes

Antibody number	#0949
Antibody reactivity (target)	Cytokeratin 7/17
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	C-46
Isotype	IgG1, kappa
Molecular weight	55 kDa (CK7); 46 kDa (CK17)
Synonyms	CK-7; K2C7; Keratin 55K Type II Cytoskeletal; Keratin Simple Epithelial Type 1 K7; Keratin Type II Cytoskeletal 7; Krt2-7; KRT7; Sarcolectin; SCL; Type II Mesothelial Keratin K7; Type-II Keratin Kb7; CK17; K17; Keratin Type I Cytoskeletal 17; Keratin-17; KRT17; PCHC1
Human gene symbol	KRT7 & KRT17
Entrez gene ID	3855; 3872
SwissProt	P08729 ; Q04695
Unigene	411501 ; 2785
Immunogen	Cytoskeletal preparation from HeLa cells (human cervical carcinoma cells)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Cow, Human
Positive control	HeLa or ME-180 cells. Skin or hepatocellular 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
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 ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Epithelial cells
Antibody research areas	Cancer, Cytoskeleton
Tumor expression	Ovarian 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>
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 November 25, 2024 at 04:33:30 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cytokeratin-7-17-monoclonal-mouse-antibody-c-46/>