Cytokeratin 8 Monoclonal Mouse Antibody (B22.1)



Product Description

Cytokeratin 8 (CK8) belongs to the type II (or B or basic) subfamily of high molecular weight cytokeratins and exists in combination with cytokeratin 18 (CK18). CK8 is primarily found in the non-squamous epithelia and is present in majority of adenocarcinomas and ductal carcinomas. It is absent in squamous cell carcinomas. Hepatocellular carcinomas are defined by the use of antibodies that recognize only cytokeratin 8 and 18. CK8 exists on several types of normal and neoplastic epithelia, including many ductal and glandular epithelia such as colon, stomach, small intestine, trachea, and esophagus as well as in transitional epithelium. Anti-CK8 does not react with skeletal muscle or nerve cells. Epithelioid sarcoma, chordoma, and adamantinoma show strong positivity corresponding to that of simple epithelia (with antibodies against CK8, CK18 and CK19). Reportedly, anti-CK8 is useful for the differentiation of lobular (ring-like, perinuclear) from ductal (peripheral-predominant) carcinoma of the breast.

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 1219, Anti-Cytokeratin 8 (B22.1)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#1219		
Antibody reactivity (target)	Cytokeratin 8		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	B22.1		
Isotype	IgG2a, kappa		
Molecular weight	52.5 kDa		
Synonyms	CARD2; CK8; CYK8; CYKER; Cytokeratin Endo A; DreK8; EndoA; K2C8; K8; Keratin 8; Krt 2.8; KRT8; Type-II Keratin Kb8		
Human gene symbol	KRT8		
Entrez gene ID	3856		
SwissProt	P05787		
Unigene	533782 & 708445		
Immunogen	Cytoskeletal preparation from HeLa cells		
Antibody target cellular localization	Cytoskeleton		
Verified antibody applications	IHC (FFPE) (verified)		
	Cow, Human, Pig, Rabbit, Sheep		
Species reactivity	Cow, Human, Pig, Rabbit, Sheep		
Species reactivity Antibody application notes	Cow, Human, Pig, Rabbit, Sheep Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/ml., Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL		
	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/ml., Does not react with mouse, rat, chicken or Xenopus; others not known,		
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL		
Antibody application notes Positive control	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma		
Antibody application notes Positive control Shipping condition	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Antibody application notes Positive control Shipping condition Storage Conditions	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Unffled: 0.2 mg/mL in PBS/0.05% BSA, Unffled: 0.2 mg/mL in PBS/0.05% BSA, Unffled: 9.2 mg/mL in PB		
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS without azide		
Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Antibody research areas	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse, rat, chicken or Xenopus; others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL MCF-7 or A431 cells. Skin, Colon, lung or breast carcinoma Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS without azide Cancer, Cytoskeleton Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in		

Email: techsupport@biotium.com

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 October 30, 2025 at 08:51:11 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cytokeratin-8-b22-1/