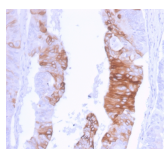


Cytokeratin 20 Recombinant Monoclonal Rabbit Antibody (KRT20/3129R)



Product Description

This MAbs recognizes an intermediate filament protein of 46 kDa, identified as cytokeratin 20 (KRT20). KRT20 is abundantly expressed in goblet cells and enterocytes of the gastrointestinal tract. It is a useful marker of pancreatic and colorectal cancer. KRT20 is expressed under normal, hyperplastic and neoplastic conditions. It has been detected in adenocarcinomas of the colon, stomach and biliary tract. Breast carcinomas are generally non-reactive.

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 3129, Anti-Cytokeratin 20 (KRT20/3129R)

Product attributes

Antibody number	#3129
Antibody reactivity (target)	Cytokeratin 20
Antibody type	Primary
Host species	Rabbit
Clonality	Monoclonal, Recombinant Monoclonal
Clone	KRT20/3129R
Isotype	IgG, kappa
Molecular weight of antigen	46 kDa
Synonyms	CK20; Cytokeratin-20; K20; KA20; Keratin 20; keratin 20; type I; Keratin type I cytoskeletal 20; Keratin-20; KRT20
Human gene symbol	KRT20
Entrez gene ID	54474
SwissProt	P35900
Unigene	84905
Immunogen	Recombinant fragment of human KRT20 protein (around aa 196-323) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human
Positive control	HT29 cell lysate (WB). Human colon carcinoma (IHC).
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	Goblet cells, Intestine
Antibody research areas	Cancer, Cytoskeleton
Tumor expression	Gastrointestinal cancer, Pancreatic cancer

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 June 20, 2026 at 02:27:27 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cytokeratin-20-recombinant-monoclonal-rabbit-antibody-krt20-3129r/>