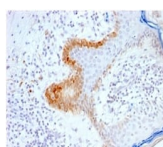


Cytokeratin 15 Recombinant Monoclonal Rabbit Antibody (KRT15/2103R)



Product Description

Keratin 15 is a type I keratin which is expressed only in basal keratinocytes in stratified epithelia and does not appear to have a natural type II expression partner. Keratin 15 is downregulated in activated keratinocytes. Cytokeratin 15 is a specific marker of stem cells of the hair-follicle bulge and may be a useful marker for diagnosis between basal cell carcinoma (BCC) and trichoepithelioma. Trichoblastoma are benign neoplasms of follicular differentiation frequently found in nevus sebaceous. Many morphologic features are shared with nodular basal cell carcinoma, sometimes rendering a diagnosis difficult. Trichoblastoma and BCC show variable expression of Cytokeratin 15 and Cytokeratin 19, and absence of hair keratins.

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 2103, Anti-Cytokeratin 15 (KRT15/2103R)

Product attributes

Antibody number	#2103
Antibody reactivity (target)	Cytokeratin 15
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	KRT15/2103R
Isotype	IgG
Molecular weight	52 kDa
Synonyms	CK15; Cytokeratin 15; K1CO; Ka15; Keratin 15 basic; Keratin 15 beta; Keratin complex 1 acidic gene 15; Keratin type I cytoskeletal 15; KRT15; KRTB; KRTL15; Type I cytoskeletal 15; Type I keratin Ka15
Human gene symbol	KRT15
Entrez gene ID	3866
SwissProt	P19012
Unigene	654570
Immunogen	Recombinant human KRT15 protein
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Cow, Human, Rat
Positive control	A431 cells. Skin or Colon
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Cancer, Cytoskeleton
Product origin	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.
Cell/tissue expression	Epidermal cells

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 August 1, 2025 at 07:48:48 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cytokeratin-15-recombinant-monoclonal-rabbit-antibody-krt15-2103r/>