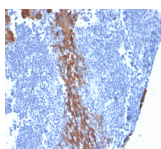


Cytokeratin 13 Monoclonal Mouse Antibody (KRT13/2213)



Product Description

Cytokeratin 13 (KRT13) is the major acidic keratin, which together with KRT4, its basic partner, is expressed in the suprabasal layers of non-cornified stratified epithelia including tongue mucosa, esophagus, anal canal epithelium, tracheal epithelium, uterine cervix, and urothelium. Defects in the KRT13 gene are a cause of white sponge nevus of cannon (WSN), a rare autosomal dominant disorder, which predominantly affects non-cornified stratified squamous epithelia and is characterized by the presence of soft, white and spongy plaques in the oral mucosa. KRT13 has been used as a marker for non-keratinized squamous epithelium. It is also expressed in various squamous metaplasia, but it is down regulated in squamous dysplasia and squamous carcinoma.

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 2659, Anti-Cytokeratin 13 (KRT13/2213)

Product attributes

Antibody number	#2659
Antibody reactivity (target)	Cytokeratin 13
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KRT13/2213
Isotype	IgG
Molecular weight	52 kDa
Synonyms	CK13; Cytokeratin-13; Keratin Type I Cytoskeletal 13; Keratin-13; KRT13; Type I Cytoskeletal 13; WSN2
Human gene symbol	KRT13
Entrez gene ID	3860
SwissProt	P13646
Unigene	654550
Immunogen	Recombinant full-length human KRT13 protein
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human
Positive control	HeLa or A431 cells. Skin or Squamous cell 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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Epidermal cells
Antibody research areas	Cancer, Cytoskeleton

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 27, 2025 at 12:30:34 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cytokeratin-13-monoclonal-mouse-antibody-krt13-2213/>