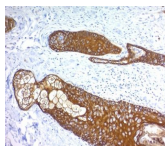


Cytokeratin, HMW Monoclonal Mouse Antibody (KRTH/1076)



Product Description

This MAbs recognizes basic (Type II or HMW) cytokeratins, which include 67 kDa (CK1); 64 kDa (CK3); 59 kDa (CK4); 58 kDa (CK5); 56 kDa (CK6); 52 kDa (CK8). Twenty human keratins are resolved with two-dimensional gel electrophoresis into acidic (pI 6.0) subfamilies. The acidic keratins have molecular weights (MW) of 56.5, 55, 51, 50, 50, 48, 46, 45, and 40 kDa. Many studies have shown the usefulness of keratins as markers in cancer research and tumor diagnosis.

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 1076, Anti-Cytokeratin, HMW (KRTH/1076)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#1076
Antibody reactivity (target)	Cytokeratin, HMW
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KRTH/1076
Isotype	IgG1, kappa
Molecular weight	67 kDa (CK1); 64 kDa (CK3); 59 kDa (CK4); 58 kDa (CK5); 56 kDa (CK6); 52 kDa (CK8)
Synonyms	KRT2B; KRT2P; HUMCYT2A; Keratin, Type II Cytoskeletal 2 Oral; K2P; K76; CK-2P; Keratin 2p; Keratin-76; Cytokeratin 2; Cytokeratin-2P; Type-II Keratin Kb9
Human gene symbol	KRT76
Entrez gene ID	51350
SwissProt	Q01546
Unigene	654392
Immunogen	Recombinant human KRT76 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human, Rat
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, Western blotting 0.5-1 ug/mL, ELISA For coating, order Ab without BSA, Predicted to show broad species reactivity
Positive control	Skin, Adenocarcinomas
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, Epithelial 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 December 11, 2025 at 11:40:40 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cytokeratin-hmw-krth1076/>