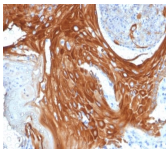




Glowing products for science

# Cytokeratin 6A Monoclonal Mouse Antibody (KRT6A/2368)



## Product Description

This MAb recognizes a protein of 56 kDa, identified as cytokeratin 6A (KRT6A). In humans, multiple isoforms of Cytokeratin 6 (6A-6F), encoded by several highly homologous genes, have distinct tissue expression patterns. Cytokeratin 6A is the dominant form in epithelial tissue. Cytokeratin 6 and 16 are expressed in keratinocytes, which are undergoing rapid turnover in the suprabasal region (also known as hyper-proliferation-related keratins). Cytokeratin 6 is found in hair follicles, suprabasal cells of a variety of internal stratified epithelia, in epidermis, in both normal and hyper-proliferative situations. Epidermal injury results in activation of keratinocytes, which express KRT6 and KRT16. KRT6 is strongly expressed in about 75% of head and neck squamous cell carcinomas..

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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2368, Anti-Cytokeratin 6A (KRT6A/2368)**

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

## Product attributes

Antibody number	#2368
Antibody reactivity (target)	Cytokeratin 6A
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	KRT6A/2368
Isotype	IgG2a, kappa
Molecular weight	56 kDa
Synonyms	58kDa Cytokeratin; CK5; Cytokeratin-5; DDD1; Epidermolysis Bullosa Simplex 2 (EBS2); Keratin 5; Keratin, Type II Cytoskeletal 5; Keratin-5; KRT5; Type-II Cytoskeletal 5; Type-II keratin Kb5
Human gene symbol	KRT6A
Entrez gene ID	3853
SwissProt	P02538
Unigene	700779
Immunogen	Recombinant full-length human Cytokeratin 6A (KRT6A) protein
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
Positive control	Tonsil or Basal Cell Carcinoma (BCC)
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
Validated in protein array	Monospecific
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
Tumor expression	Squamous cell carcinoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 12, 2025 at 12:34:56 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cytokeratin-6a-monoclonal-mouse-antibody-kr6a-2368/>