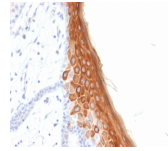


# Cytokeratin 10 Recombinant Monoclonal Mouse Antibody (rKRT10/844)



## Product Description

This MAb recognizes a protein of 56.5 kDa, identified as cytokeratin 10 (CK10). CK10 is expressed in all suprabasal layers of the epidermis. In the epidermis, expression of CK10 strictly parallels the extent of differentiation; it is absent in the basal layer, appears in the first suprabasal layers and increases in concentration towards the granular layer. However, CK10 is rarely detected in early stages of vulvar squamous carcinomas (tumors less than 2 cm, clinical stage I) regardless of the tumor grade. In larger and more advanced tumors (greater than 2 cm, clinical stages II and III), CK10 is detected very frequently. Expression of CK10 is related to maturation of malignant keratinocytes, being preferentially detected in more-differentiated parts. 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 1989, Anti-Cytokeratin 10 (rKRT10/844)**

## Product attributes

Antibody number	#1989
Antibody reactivity (target)	Cytokeratin 10, KRT10
Antibody type	Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rKRT10/844
Isotype	IgG1, kappa
Molecular weight	56.5 kDa
Synonyms	BCIE; BIE; EHK; Keratin Type I Cytoskeletal 10; KRT10
Human gene symbol	KRT10
Entrez gene ID	3858
SwissProt	P13645
Unigene	99936
Immunogen	Recombinant full-length human KRT10 protein.
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoskeleton
Species reactivity	Human, Mouse
Positive control	A431, HeLa, MCF7 cells. Esophagus or tonsil.
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
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>
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, LI-COR Bioscience.