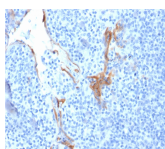


# Cytokeratin 16 Monoclonal Mouse Antibody (LL025)



## Product Description

Cytokeratins are a family of intermediate filament proteins that assemble into filaments through forming heterodimers of one type I Cytokeratins (Cytokeratins 9 to 23) and one type II Cytokeratins (keratins 1 to 8). The cytokeratin proteins play a critical role in differentiation, as well as tissue specialization and function, to maintain the overall structural integrity of epithelial cells. Cytokeratins are also useful markers in identifying the origin of metastatic tumors. Cytokeratin 16 is expressed in benign stratified squamous epithelium and squamous cell carcinoma of the head and neck, as well as luminal cells of mammary gland and sweat ducts. It is absent in non-invasive breast carcinomas and normal breast tissue.

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 1692, Anti-Cytokeratin 16 (LL025)**

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

## Product attributes

<b>Antibody number</b>	#1692
<b>Antibody reactivity (target)</b>	Cytokeratin 16
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	LL025
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight of antigen</b>	48 kDa
<b>Synonyms</b>	CK16; Cytokeratin-16; Focal non epidermolytic palmoplantar keratoderma (FNEPPK / NEPPK); K1CP; Keratin Type I Cytoskeletal 16; Keratin-16; KRT16A
<b>Human gene symbol</b>	KRT16
<b>Entrez gene ID</b>	3868
<b>SwissProt</b>	P08779
<b>Unigene</b>	655160
<b>Immunogen</b>	A synthetic peptide from the C-terminal of human Cytokeratin 16
<b>Verified antibody applications</b>	IHC (FPPE) (verified)
<b>Antibody target cellular localization</b>	Cytoskeleton
<b>Species reactivity</b>	Human
<b>Positive control</b>	HeLa cells, Tonsil, Skin, Bladder, or Cervix.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35°C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	Epidermal cells
<b>Antibody research areas</b>	Cancer, Cytoskeleton

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 June 4, 2026 at 11:12:45 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cytokeratin-16-monoclonal-mouse-antibody-ll025/>