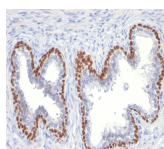


# p63 Monoclonal Mouse Antibody (TP63/11)



## Product Description

p63 is a homolog of the tumor suppressor p53. It is identified in basal cells in the epithelial layers of a variety of tissues, including epidermis, cervix, urothelium, breast and prostate. p63 was detected in nuclei of the basal epithelium in normal prostate glands; however, it is not expressed in malignant tumors of the prostate. As a result, p63 has been reported as a useful marker for differentiating benign from malignant lesions in the prostate, particularly when used in combination with markers of high molecular weight cytokeratins and the prostate-specific marker AMACR (P504S). p63 has also been shown to be a sensitive marker for lung squamous cell carcinomas (SqCC), with a sensitivity of ~90%. Specificity for lung SqCC, vs. lung adenocarcinoma (LADC), is approximately 80%. In breast tissue, p63 has been identified in myoepithelial cells of normal ducts.

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 0011, Anti-p63 (TP63/11)**

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

## Product attributes

<b>Isotype</b>	IgG2a, kappa
<b>Antibody number</b>	#0011
<b>Antibody reactivity (target)</b>	p63
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	TP63/11
<b>Molecular weight of antigen</b>	63 kDa
<b>Synonyms</b>	Amplified in squamous cell carcinoma (AIS); Chronic ulcerative stomatitis protein (CUSP); EEC3; Keratinocyte transcription factor KET; LMS; NBP; p40; P51/P63; p53 like transcription factor; p53-related protein p63; RHS; SHFM4; TAp63alpha; TP53CP; TP53L; TP63; TP73; TP73L; Transformation-related protein 63; Trp53rp1; Trp6; 3; Tumor protein 63; Tumor protein p53-like; tumor protein p73-like
<b>Human gene symbol</b>	TP63
<b>Entrez gene ID</b>	8626
<b>SwissProt</b>	Q9H3D4
<b>Unigene</b>	137569
<b>Immunogen</b>	Recombinant full-length human p63 protein
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Antibody target cellular localization</b>	Nucleus
<b>Species reactivity</b>	Human, Mouse
<b>Positive control</b>	HEK293 cells, Prostate Carcinoma or Lung or bladder squamous cell carcinoma.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<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>Tumor expression</b>	Lung cancer, Prostate cancer, 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 June 3, 2026 at 04:53:41 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/p63-monoclonal-mouse-antibody-tp63-11/>