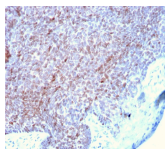


p73 Monoclonal Mouse Antibody (P73/2531)



Product Description

p73 shares considerable sequence similarity with the widely-studied tumor suppressor gene p53. The p73 gene maps to a region in chromosome 1 that is frequently deleted in neuroblastomas. However, unlike p53, p73 does not appear to be activated by DNA damaging agents.

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 2531, Anti-p73 (P73/2531)

Product attributes

Antibody number	#2531
Antibody reactivity (target)	p73
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	P73/2531
Isotype	IgG2b, kappa
Molecular weight	73 kDa
Synonyms	p53-like transcription factor; p53-related protein; p73; TP73; Tumor protein p73
Human gene symbol	TP73
Entrez gene ID	7161
SwissProt	O15350
Unigene	697294; 706990
Immunogen	Recombinant full-length human TP73 protein
Antibody target cellular localization	Golgi apparatus, Nucleus
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human
Positive control	HeLa, Jurkat, HEK293 cell lysates. Human kidney or bladder carcinoma.
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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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.

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 August 15, 2025 at 04:53:28 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/p73-monoclonal-mouse-antibody-p73-2531/>