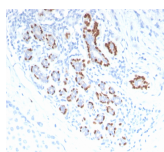


p40 Polyclonal Rabbit Antibody



Product Description

p40 (p63 delta) is a marker recently determined to be highly specific for squamous basal cells in the immunohistochemistry (IHC) application. The current more routinely recommended marker, p63, appears to have less specificity compared to p40, especially on squamous cell tumors. The ability to differentiate between lung adenocarcinoma vs. squamous cell carcinoma is difficult and has bearing on the different therapeutic avenues for each subtype treatment. p63 antibody's ability to distinguish between the tumor types appears to be inferior when compared to p40. The ability to utilize an antibody probe for p40 as a squamous cell marker bolsters its use for future sub-classification of lung cancers, especially by immunohistochemical techniques.

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 2314, Anti-p40

Product attributes

Antibody number	#2314
Antibody reactivity (target)	p40
Antibody type	Primary
Host species	Rabbit
Clonality	Polyclonal
Isotype	IgG, kappa
Molecular weight of antigen	40 kDa
Synonyms	deltaNp63; TP63
Human gene symbol	TP63
Entrez gene ID	8626
SwissProt	Q9H3D4
Unigene	137569
Immunogen	Recombinant human p40 protein fragment
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Positive control	Human prostate or lung tissue (IHC).
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.
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	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 June 20, 2026 at 07:40:08 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/p40-polyclonal-rabbit-antibody/>