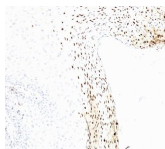


Human Papillomavirus 16 (HPV-16) Monoclonal Mouse Antibody (CAMVIR-1)



Product Description

Reacts with a protein of 57 kDa, identified as the L1 protein of human papilloma virus type 16 (HPV-16). It is the major capsid protein of HPV-16. Infection with specific types of HPV has been associated with an increased risk of developing cervical neoplasia. HPV types 6 and 11 have been associated with relatively benign diseases such as genital warts but types 16 and 18 are strongly associated with cervical, vaginal, and vulvar malignancies. The antibody reacts very strongly with formalin-fixed, paraffin-embedded tissues containing HPV-16 or -33; very weak reactions were occasionally observed with biopsy specimens or smears containing HPV-6 or HPV-11. It cross-reacts with HPV37.

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 0502, Anti-Human Papillomavirus-16 (CAMVIR-1)

Product attributes

Antibody number	#0502
Antibody reactivity (target)	Human Papillomavirus-16
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CAMVIR-1
Isotype	IgG2a, kappa
Molecular weight	57 kDa
Synonyms	HPV-16; HPV-16 capsid; HPV16 L1; HPV16 major capsid protein L1; Human papillomavirus type 16 L1; Human papillomavirus type 16 major capsid protein L1
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Human papilloma virus type 16, major capsid protein L1
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	HPV-16
Antibody application notes	Immunohistology formalin-paraffin 0.5-1.0 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Immunofluorescence 0.5-1 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	HPV-16 infected cells or cervical tissue
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	Infectious disease, Virology
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 20, 2025 at 07:59:55 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/monoclonal-anti-human-papillomavirus-16-hpv16-camvir-1/>