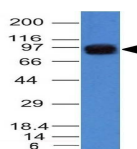


UACA / Nucling Monoclonal Mouse Antibody (AE-5)



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

Product Description

UACA (Uveal Autoantigen with Coiled-coil domains and Ankyrin repeats) is a 1,416 amino acid nuclear membrane protein. It was originally identified as an autoantigen in patients with panuveitis, a characteristic of Vogt-Koyanagi-Harada disease, and in patients with Graves' disease. UACA was also later identified as Nucling, an mRNA differentially expressed in F9 embryonal carcinoma cells during cardiac muscle differentiation. UACA appears to function as a pro-apoptotic protein that recruits the apaf-1-pro-caspase-9 complex for the induction of apoptosis to mediate the cell-death pathway.

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 0956, Anti-Nucling (AE-5)

Product attributes

Antibody number	#0956
Antibody reactivity (target)	Nucling, UACA
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	AE-5
Isotype	IgG1, kappa
Molecular weight of antigen	160 kDa
Synonyms	Nuclear membrane binding protein; Uveal autoantigen with coiled-coil domains and ankyrin repeats (UACA); NUCLING
Human gene symbol	UACA
Entrez gene ID	55075
SwissProt	Q9BZF9
Unigene	108049
Immunogen	Nuclei of myeloid leukemia biopsy cells
Verified antibody applications	WB (verified)
Antibody target cellular localization	Nuclear membrane
Species reactivity	Human, Mouse
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 0.5-1.0 ug/mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	HeLa or 293T cells. Highly expressed in skeletal muscle, heart, kidney and pancreas. Also expressed in epidermal melanocytes, eye muscles and thyroid follicular cells.
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	Apoptosis
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), 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 June 7, 2026 at 03:46:37 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-uaca-nucling-ae-5/>