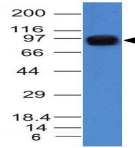


## UACA / Nucling Monoclonal Mouse Antibody (AE-5)

UACA (Uveal Autoantigen with Coiled-coil domains and Ankyrin repeats) is a 1,416 amino acid nuclear membrane protein.



### 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](mailto: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)**

| Antibody # | prefix                 | Conjugation | Ex/Em                | Concentration | Storage Buffer              |
|------------|------------------------|-------------|----------------------|---------------|-----------------------------|
| BNC04      | CF®405S                |             | 404/431 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC05      | CF®405M                |             | 408/452 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC06      | CF®405L                |             | 395/545 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC88      | CF®488A                |             | 490/515 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC14      | CF®514                 |             | 516/548 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC43      | CF®543                 |             | 541/560 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC55      | CF®555                 |             | 555/565 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC68      | CF®568                 |             | 562/583 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC94      | CF®594                 |             | 593/614 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC40      | CF®640R                |             | 642/662 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC47      | CF®647                 |             | 650/665 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC60      | CF®660C                |             | 667/685 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC61      | CF®660R                |             | 663/682 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC80      | CF®680                 |             | 681/698 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC81      | CF®680R                |             | 680/701 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC00      | CF®700                 |             | 695/720 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNC70      | CF®770                 |             | 770/797 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCR       | R-PE (PE)              |             | 496, 546, 565/578 nm | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCA       | APC                    |             | 650/660 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCP       | PerCP                  |             | 482/677 nm           | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCB       | Biotin                 | N/A         |                      | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCAP      | Alkaline Phosphatase   | N/A         |                      | 0.1 mg/mL     | PBS, 0.1% BSA, 0.05% azide  |
| BNCH       | Horseradish Peroxidase | N/A         |                      | 0.1 mg/mL     | PBS, 0.05% BSA, no azide    |
| BNUB       | Purified, with BSA     | N/A         |                      | 0.2 mg/mL     | PBS, 0.05% BSA, 0.05% azide |
| BNUM       | Purified, BSA-free     | N/A         |                      | 1 mg/mL       | PBS, no BSA, no azide       |

### References

Yamada, K., et al. 2001. Identification of a novel autoantigen UACA in patients with panuveitis. *Biochem. Biophys. Res. Commun.* 280: 1169-1176. | Ohkura, T., et al. 2004. Detection of the novel autoantibody (anti-UACA antibody) in patients with Graves disease. *Biochem. Biophys. Res. Commun.* 321: 432-440

### Product attributes

|                         |   |
|-------------------------|---|
| Antibody number         | 0956  |
| Reactivity (target)     | Nucling, UACA   |
| Antibody type           | Primary   |
| Host species            | Mouse   |
| Clonality               | Monoclonal  |
| Clone                   | AE-5  |
| Isotype                 | IgG1, kappa   |
| Molecular weight        | 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   |
| Cellular localization   | Nuclear membrane  |
| Species reactivity      | Human, Mouse  |
| Applications            | Immunofluorescence, Flow cytometry, Western   |
| Application notes       | 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)   |
| Supplied As             | 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, BSA-free: 1 mg/mL in PBS without azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide |
| Antibody research areas | Apoptosis   |