## ORC1 Monoclonal Mouse Antibody (7F6/1)



## **Product Description**

The initiation of DNA replication is a multi-step process that depends on the formation of pre-replication complexes, which trigger initiation. Among the proteins required for establishing these complexes are the origin recognition complex (ORC) proteins. ORC proteins bind specifically to origins of replication where they serve as scaffold for the assembly of additional initiation factors. Human ORC subunits 1-6 are expressed in the nucleus of proliferating cells and tissues, such as the testis. ORC1 and ORC2 are both expressed at equivalent concentrations throughout the cell cycle; however, only ORC2 remains stably bound to chromatin. ORC4 and ORC6 are also expressed constantly throughout the cell cycle. ORC2, ORC3, ORC4 and ORC5 form a core complex upon which ORC6 and ORC1 assemble. The formation of this core complex suggests that ORC proteins play a crucial role in the G1-S transition in mammalian cells.

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3067, Anti-ORC1 (7F6/1)

## Product attributes

| Product attributes                    |   |  |
|---------------------------------------|---|--|
| Antibody number                       | #3067   |  |
| Antibody reactivity (target)          | ORC1  |  |
| Antibody type                         | Primary   |  |
| Host species                          | Mouse   |  |
| Clonality                             | Monoclonal  |  |
| Clone                                 | 7F6/1   |  |
| Isotype                               | IgG1, kappa   |  |
| Molecular weight                      | 120 kDa   |  |
| Synonyms                              | HSORC1; MmORC1; Origin Recognition Complex 1 (ORC1);<br>Origin recognition complex subunit 1 (yeast homolog) like<br>(ORC1L); Origin recognition complex subunit 1 like (S. cerevisia;<br>Origin recognition complex subunit 1 like; PARC1; Replication<br>control protein 1  |  |
| Human gene symbol                     | ORC1L   |  |
| Entrez gene ID                        | 4998  |  |
| SwissProt                             | Q13415  |  |
| Unigene                               | 17908   |  |
| Immunogen                             | Recombinant full-length human ORC1 protein  |  |
| Antibody target cellular localization | Nucleus   |  |
| Species reactivity                    | Human   |  |
| Antibody application notes            | For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |  |
| Positive control                      | HeLa 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   |  |
| 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.   |  |

Call us: 800-304-5357 Email: btinfo@biotium.com

| Antibody # prefix | Conjugation           | Ex/Em (nm) | Laser line | Detection channel           | Dye Features     |
|-------------------|-----------------------|------------|------------|-----------------------------|------------------|
| BNC04             | CF®405S               | 404/431    | 405        | DAPI (microscopy),<br>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,<br>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 November 4, 2025 at 11:00:47 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/orc1-monoclonal-mouse-antibody-7f6-1/">https://biotium.com/product/orc1-monoclonal-mouse-antibody-7f6-1/</a>