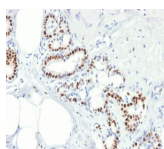


# Estrogen Receptor, Alpha Recombinant Monoclonal Mouse Antibody (rESR1/1935)



## Product Description

This MAb is specific to ER alpha and shows minimal cross-reaction with other members of the family. ER is an important regulator of growth and differentiation in the mammary gland. Presence of ER in breast tumors indicates an increased likelihood of response to anti-estrogen (e.g. tamoxifen) therapy. This MAb is excellent for staining of formalin-fixed, paraffin-embedded breast carcinomas.

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 2300, Anti-Estrogen Receptor, alpha (rESR1/1935)**

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

## Product attributes

|  |  |
|--|--|
| <b>Antibody number</b>                       | #2300  |
| <b>Antibody reactivity (target)</b>          | Estrogen Receptor, alpha   |
| <b>Antibody type</b>                         | Primary  |
| <b>Host species</b>                          | Mouse  |
| <b>Clonality</b>                             | Monoclonal, Recombinant Monoclonal   |
| <b>Clone</b>                                 | rESR1/1935   |
| <b>Isotype</b>                               | IgG1   |
| <b>Molecular weight of antigen</b>           | 67 kDa   |
| <b>Synonyms</b>                              | Estrogen Receptor alpha delta 4*5,6,7/654 isoform; Estrogen Receptor alpha delta 4 +49 isoform; Nuclear receptor subfamily 3 group A member 1  |
| <b>Human gene symbol</b>                     | ESR1   |
| <b>Entrez gene ID</b>                        | 2099   |
| <b>SwissProt</b>                             | P03372   |
| <b>Unigene</b>                               | 208124   |
| <b>Immunogen</b>                             | Recombinant full-length human Estrogen Receptor alpha protein  |
| <b>Antibody target cellular localization</b> | Nucleus  |
| <b>Verified antibody applications</b>        | IHC (FFPE) (verified), WB (verified)   |
| <b>Species reactivity</b>                    | Human  |
| <b>Antibody application notes</b>            | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Western blotting 1-2 ug/mL. Optimal dilution for a specific application should be determined by user |
| <b>Positive control</b>                      | MCF-7 Cells or Breast Cancers.   |
| <b>Shipping condition</b>                    | Room temperature   |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended  |
| <b>Regulatory status</b>                     | For research use only (RUO)  |
| <b>Antibody/conjugate formulation</b>        | 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  |
| <b>Antibody research areas</b>               | Cancer, Endocrinology, Nuclear receptors   |
| <b>Product origin</b>                        | 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.  |
| <b>Tumor expression</b>                      | Breast cancer  |

| Antibody # prefix | Conjugation        | Ex/Em (nm) | Laser line | Detection channel        | Dye Features                     |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04             | CF®405S            | 404/431    | 405        | DAPI (microscopy), AF405 | <a href="#">CF®405S Features</a> |
| BNC88             | CF®488A            | 490/515    | 488        | GFP, FITC                | <a href="#">CF®488A Features</a> |
| BNC68             | CF®568             | 562/583    | 532, 561   | RFP, TRITC               | <a href="#">CF®568 Features</a>  |
| BNC94             | CF®594             | 593/614    | 561        | Texas Red®               | <a href="#">CF®594 Features</a>  |
| BNC40             | CF®640R            | 642/662    | 633-640    | Cy®5                     | <a href="#">CF®640R Features</a> |
| BNC47             | CF®647             | 650/665    | 633-640    | Cy®5                     | <a href="#">CF®647 Features</a>  |
| BNC74             | CF®740             | 742/767    | 633-685    | 775/50                   | <a href="#">CF®740 Features</a>  |
| 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 4, 2026 at 01:23:16 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/estrogen-receptor-alpha-recombinant-monoclonal-mouse-antibody-resr1-1935/>