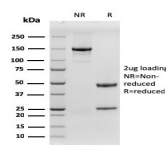


EGFR Recombinant Monoclonal Mouse Antibody (rGFR/1667)



Product Description

This MAb recognizes a protein of 170 kDa, identified as EGFR. EGFR is type I receptor tyrosine kinase with sequence homology to erbB-1, -2, -3 -4 or HER-1, -2, -3 -4. It binds to Epidermal Growth Factor (EGF), Transforming Growth Factor- α (TGF- α), Heparin-binding EGF (HB-EGF), amphiregulin, betacellulin and epiregulin. EGFR is overexpressed in tumors of breast, brain, bladder, lung, gastric, head neck, esophagus, cervix, vulva, ovary, and endometrium. It is predominantly present in squamous cell 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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 3605, Anti-EGFR (rGFR/1667)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #3605 |
| Antibody reactivity (target) | EGFR |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Recombinant Monoclonal |
| Clone | rGFR/1667 |
| Isotype | IgG1, kappa |
| Molecular weight | ~170 kDa (wild type) and ~145 kDa (vIII variant) |
| Synonyms | ErbB1; ERBB1; Errp; HER1; mENA; PIG61; Proto-oncogene c-ErbB-1; Receptor Tyrosine Protein Kinase; ErbB1; Urogastrone; wa2; Wa5 |
| Human gene symbol | EGFR |
| Entrez gene ID | 1956 |
| SwissProt | P00533 |
| Unigene | 488293 |
| Immunogen | Recombinant full-length human EGFR protein |
| Verified antibody applications | WB (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Positive control | A431 cells. Placenta, breast, colon or bladder cancer (IHC). |
| 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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |
| Tumor expression | Squamous cell carcinoma |

| 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 September 6, 2025 at 11:10:24 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/egfr-recombinant-monoclonal-mouse-antibody-rgr-1667/>