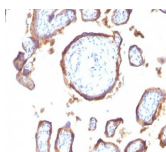


EGFR Monoclonal Mouse Antibody (GFR1195)



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, β cellulin 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 1195, Anti-EGFR (GFR1195)

Product attributes

| | |
|---------------------------------------|---|
| Antibody number | #1195 |
| Antibody reactivity (target) | EGFR |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | GFR1195 |
| Isotype | IgG2a, 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 human EGFR protein |
| Verified antibody applications | IHC (FPPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Antibody application notes | Immunohistology formalin-fixed 2-4 ug/mL, Digest formalin-fixed tissues with Protease Sigma Cat. #P-6911 at 1 mg/mL PBS, pH 7.4 for 10 min at 37C, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Optimal dilution for a specific application should be determined by user |
| Positive control | A431 cells, Placenta, Breast, Colon or Bladder cancer |
| 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 | Cancer, Signal transduction |
| 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. |
| 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 December 23, 2024 at 07:26:23 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-egfr-gfr1195/>