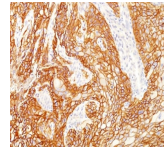


EGFRvIII Recombinant Monoclonal Rabbit Antibody (GFR/2600R)



Product Description

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 2600, Anti-EGFRvIII (GFR/2600R)**

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #2600 |
| Antibody reactivity (target) | EGFRvIII |
| Antibody type | Primary |
| Host species | Rabbit |
| Clonality | Recombinant Monoclonal |
| Clone | GFR/2600R |
| Isotype | IgG |
| Molecular weight | ~145 kDa (vIII variant) |
| Synonyms | Erb1 variant III; ERBB1 variant III; HER1 variant III; Proto-oncogene c-ErbB-1 variant III; Receptor Tyrosine Protein Kinase variant III; ErbB1 variant III |
| Human gene symbol | EGFR |
| Entrez gene ID | 1956 |
| SwissProt | P00533 |
| Unigene | 488293 |
| Immunogen | KLH conjugate of EGFRvIII-specific, 14-amino acid peptide from the predicted amino acid sequence at the fusion junction (LEEKKNYVVDHDC) |
| Antibody target cellular localization | Plasma membrane |
| Verified antibody applications | IHC (FFPE) (verified) |
| Species reactivity | Human |
| Antibody application notes | 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. Optimal dilution for a specific application should be determined by user |
| Positive control | A431 cells, Placenta, Breast, Colon or Bladder Carcinoma |
| 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 |
| 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 |
| 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, LI-COR Bioscience.