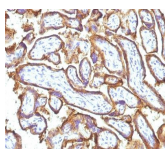


# EGFR Monoclonal Mouse Antibody (GFR1195 + 31G7)



## 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- $\alpha$  (TGF- $\alpha$ ), Heparin-binding EGF (HB-EGF), amphiregulin,  $\beta$ 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<sup>®</sup> dyes and other labels. CF<sup>®</sup> dyes offer exceptional brightness and photostability. See the [CF<sup>®</sup> Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF<sup>®</sup>405S and CF<sup>®</sup>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<sup>®</sup> 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 1200, Anti-EGFR (GFR1195 31G7)**

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

## Product attributes

<b>Antibody number</b>	#1200
<b>Antibody reactivity (target)</b>	EGFR
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	GFR1195 + 31G7
<b>Isotype</b>	IgG2a, kappa
<b>Molecular weight of antigen</b>	~170 kDa (wild type) and ~145 kDa (vIII variant)
<b>Synonyms</b>	ErbB1; ERBB1; Errp; HER1; mENA; PIG61; Proto-oncogene c-ErbB-1; Receptor Tyrosine Protein Kinase; ErbB1; Urogastrone; wa2; Wa5
<b>Human gene symbol</b>	EGFR
<b>Entrez gene ID</b>	1956
<b>SwissProt</b>	P00533
<b>Unigene</b>	488293
<b>Immunogen</b>	Human EGFR purified from A431 cells (31G7); Recombinant human EGFR protein (GFR1195)
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	Immunohistochemistry 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
<b>Positive control</b>	A431 cells. Placenta, Breast, Colon or Bladder cancer
<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, Signal transduction
<b>Product origin</b>	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.
<b>Tumor expression</b>	Squamous cell carcinoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF <sup>®</sup> 405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF<sup>®</sup>405S Features</a>
BNC88	CF <sup>®</sup> 488A	490/515	488	GFP, FITC	<a href="#">CF<sup>®</sup>488A Features</a>
BNC68	CF <sup>®</sup> 568	562/583	532, 561	RFP, TRITC	<a href="#">CF<sup>®</sup>568 Features</a>
BNC94	CF <sup>®</sup> 594	593/614	561	Texas Red <sup>®</sup>	<a href="#">CF<sup>®</sup>594 Features</a>
BNC40	CF <sup>®</sup> 640R	642/662	633-640	Cy <sup>®</sup> 5	<a href="#">CF<sup>®</sup>640R Features</a>
BNC47	CF <sup>®</sup> 647	650/665	633-640	Cy <sup>®</sup> 5	<a href="#">CF<sup>®</sup>647 Features</a>
BNC74	CF <sup>®</sup> 740	742/767	633-685	775/50	<a href="#">CF<sup>®</sup>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 9, 2026 at 12:50:04 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-egfr-gfr1195-31g7/>