

# EGFR Monoclonal Mouse Antibody (B1D8)

## Product Description

This antibody reacts with the extracellular domain of EGFR and blocks the EGF/TGF induced activation. It also blocks tumor growth in vivo. It is excellent for purification of 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. Primary antibodies are available purified, or with a selection of fluorescent CF $\oplus$  dyes and other labels. CF $\oplus$  dyes offer exceptional brightness and photostability. See the [CF \$\oplus\$  Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF $\oplus$ 405S and CF $\oplus$ 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 $\oplus$  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 0643, Anti-EGFR (B1D8)**

## References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. J Biol Chem (1984) 259(19): 11895-11900. (IP, functional studies)
2. J Immunol Meth (2004) 287: 147-158. (Flow; epitope mapping)

### Product attributes

Antibody number	#0643
Antibody reactivity (target)	EGFR
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	B1D8
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	Microsomes from A431 cells.
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IP (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Does not react with mouse or rat, others not known, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Affinity Purification, for coupling, order Ab without BSA, Immunoprecipitation 1-2 ug/500 ug protein lysate. Optimal dilution for a specific application should be determined by user
Positive control	A431 cells, Breast 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
Tumor expression	Squamous cell carcinoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF $\oplus$ 405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF<math>\oplus</math>405S Features</a>
BNC88	CF $\oplus$ 488A	490/515	488	GFP, FITC	<a href="#">CF<math>\oplus</math>488A Features</a>
BNC68	CF $\oplus$ 568	562/583	532, 561	RFP, TRITC	<a href="#">CF<math>\oplus</math>568 Features</a>
BNC94	CF $\oplus$ 594	593/614	561	Texas Red $\oplus$	<a href="#">CF<math>\oplus</math>594 Features</a>
BNC40	CF $\oplus$ 640R	642/662	633-640	Cy $\oplus$ 5	<a href="#">CF<math>\oplus</math>640R Features</a>
BNC47	CF $\oplus$ 647	650/665	633-640	Cy $\oplus$ 5	<a href="#">CF<math>\oplus</math>647 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, LI-COR Bioscience.