

EGF receptor monoclonal antibody (6F1)

The epidermal growth factor receptor (EGFR) is a 170-kDa membrane bound receptor consisting of an extracellular ligand-binding domain, a single hydrophobic transmembrane region, and a tyrosine kinase domain-containing cytoplasmic region. Ligand binding induces receptor homo- or heterodimerization, triggering autophosphorylation of cytoplasmic tyrosine residues that provide docking sites for Src homology 2 (SH2) domain-containing signaling molecules. Src-mediated phosphorylation of EGFR at Tyr845 may contribute to EGFR hyperactivity in cancer, while autophosphorylation of EGFR at Tyr1068 and/or Tyr1173 in response to ligand binding is known to recruit cofactors (e.g. SHP1, RasGAP, SOS1, Grb2) that activate pathways associated with cell growth, apoptosis, adhesion, receptor endocytosis, and protein degradation.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

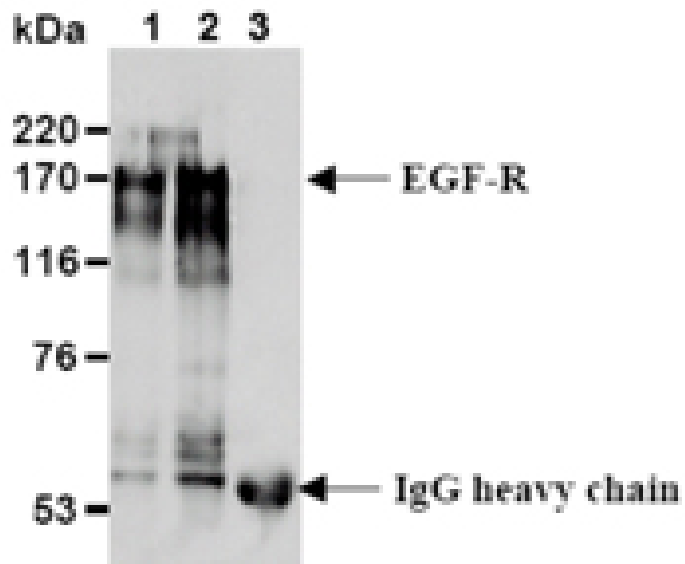
Ordering Information

[Order Online »](#)

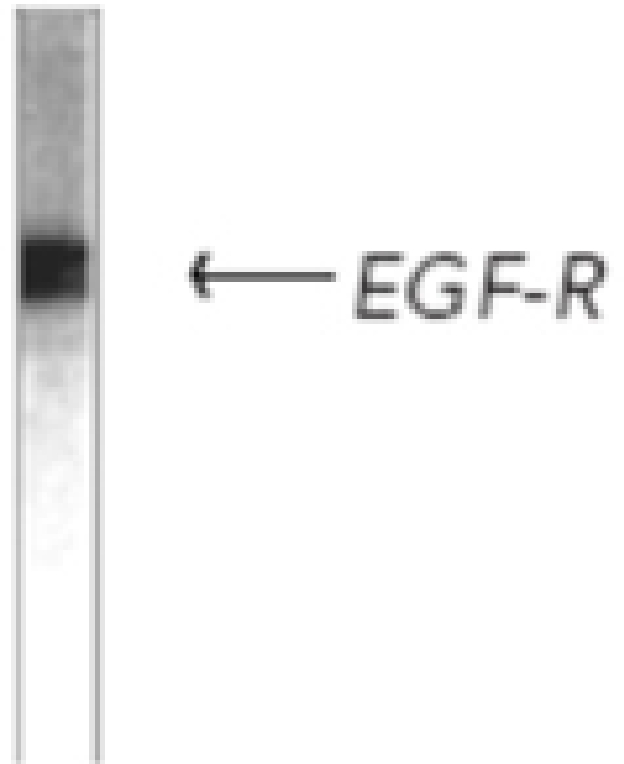
ADI-CSA-330-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunoprecipitation analysis of human EGF-R from A431 cells with 2µg (1) and 10µg (2) of EGFR mAb (6F1), and mouse IgG (3).



Western blot analysis of human A431 cell lysate probed with EGFR mAb (6F1).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	EGFR, epidermal growth factor receptor
Application	Flow Cytometry, ICC, IP, WB
Application Notes	Detects a band of ~170kDa by Western blot.
Clone	6F1
Formulation	Liquid. In PBS, pH 7.2 containing 50% glycerol.
Host	Mouse
Immunogen	Recombinant human EGF receptor (C-terminal fragment).
Isotype	IgG2b
Purity Detail	Ammonium sulfate precipitation purified.
Recommendation Dilutions/Conditions	Flow cytometry: 1:100 (final concentration)Immunocytochemistry: 1:100Immunoprecipitation: 2-5 µL/250 µL of cell extract from 2.5 x 10 ⁶ cellsWestern Blot: 1:1,000-1:2,000 for chemiluminescence detection systemSuggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P00533
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)