

EGF receptor (human) Recombinant monoclonal antibody (528) (Mouse IgG2ak)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ErbB-1, HER1, Epidermal growth factor receptor, EGFR
Application	Flow Cytometry, IF, IHC (PS), IP, WB
Clone	528
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Partially purified EGFR from A431 cells.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P00533
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee