EGF receptor (human) Recombinant monoclonal antibody (C225) (Rabbit IgGκ)

The epidermal growth factor receptor is the cell-surface receptor for members of the epidermal growth factor family (EGF-family) of extracellular protein ligands.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS485-0200

200µg

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, FUNC

Clone C225 (Cetuximab)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Human EGFR.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original human IgG1 format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P00533

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com