[pTyr¹¹⁵⁰/Tyr¹¹⁵¹]Insulin receptor monoclonal antibody (10C3)

The Insulin Receptor (InsR) is a heterodimeric receptor tyrosine kinase with an extracellular alpha-chain, a transmembrane domain and an intracellular beta-chain. InsR is activated upon binding of the peptide hormone insulin, leading to autophosphorylation of tyrosine residues 1146, 1150, and 1151 in the activation loop of the beta-chain. Additional autophosphorylation sites such as tyrosine residues 960, 972, 1316, and 1322 regulate the assembly of signal transduction complexes.

This antibody is covered by our Worry-Free Guarantee.

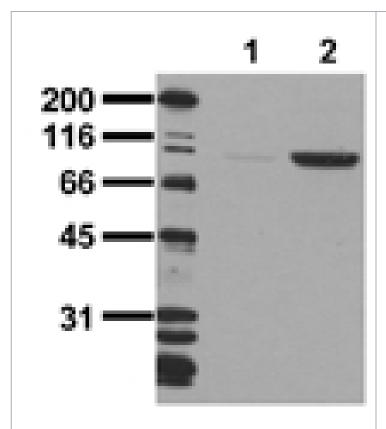
Ordering Information

Order Online »

ADI-905-645-100 100μg

Manuals, SDS & CofA

View Online »



Western blot analysis of untreated (1) or insulinstimulated (2) MCA-MB-213 cancer cells probed with Insulin Receptor (pTyr1150/1151) mAb (10C3).

Handling & Storage

Use/Stability Stable at -20°C up to 1 year.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name INSR

Application ELISA, WB

Clone 10C3

Crossreactivity Cross-reacts with IGF1 receptor.

Formulation Liquid. In PBS containing 50% glycerol, PEG, sucrose,

and 0.09% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of human

Insulin Receptor phosphorylated at Tyr1150/1151,

conjugated to KLH.

lsotype lgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

Recommendation Dilutions/Conditions ELISA (0.1µg/ml)Western Blot (0.2-0.5µg/ml,

ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P06213

Last modified: May 29, 2024



info-

eu@enzolifesciences.com