# [pTyr1322]Insulin receptor monoclonal antibody (21G12)

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.

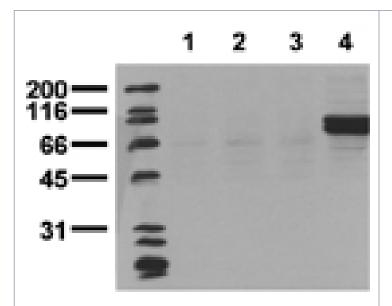
# **Ordering Information**

Order Online »

**ADI-905-646-100** 100μg

Manuals, SDS & CofA

**View Online** »



Western blot analysis of MDA-MB-231 cell lysates, untreated (1), Insulin treated (2), IGF-1 treated (3), or pervanadate treated (4), probed with Insulin Receptor (pTyr1322) mAb (21G12).

## **Handling & Storage**

**Use/Stability** Stable at -80°C up to 1 year, at 4°C up to 3 months.

Long Term Storage -20°C

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name INSR

**Application** ELISA, WB

**Application Notes**Detects a band of ~97kDa by Western blot.

**Clone** 21G12

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium

azide, PEG, and sucrose.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to a portion of insulin

receptor phosphorylated at Tyr1322.

lgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

Recommendation Dilutions/Conditions ELISA (0.1µg/ml)Western Blot (0.5µg/ml, ECL)Suggested

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

individually for each application.

**Reconstitution** Reconstitute with 1ml water (15 minutes at room

temperature).

Species Reactivity Dog, Human, Mouse

UniProt ID P06213

Last modified: May 29, 2024



info-

eu@enzolifesciences.com