

# [pTyr<sup>1322</sup>]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](#).

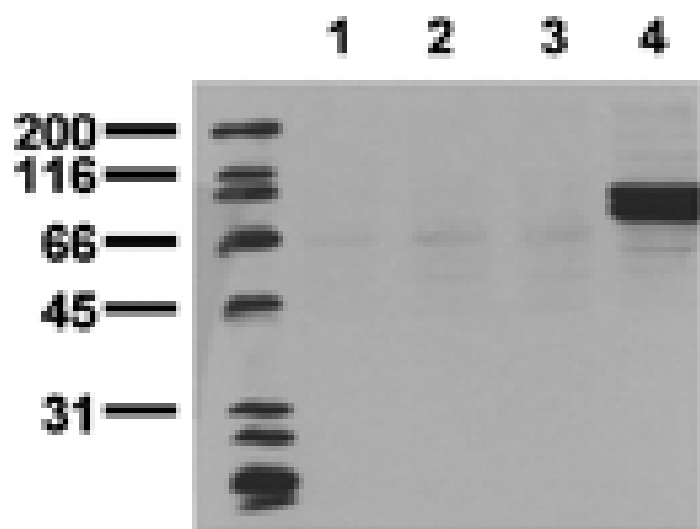
## Ordering Information

[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.

**Isotype** IgG1

**Purity Detail** Thiophilic adsorption and size exclusion chromatography purified.

**Recommendation** ELISA (0.1µg/ml)Western Blot (0.5µg/ml, ECL)Suggested dilutions/conditions may not  
**Dilutions/Conditions** 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



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

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +32 3 466 0420  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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

**Worry-free Guarantee**

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