

[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](#).

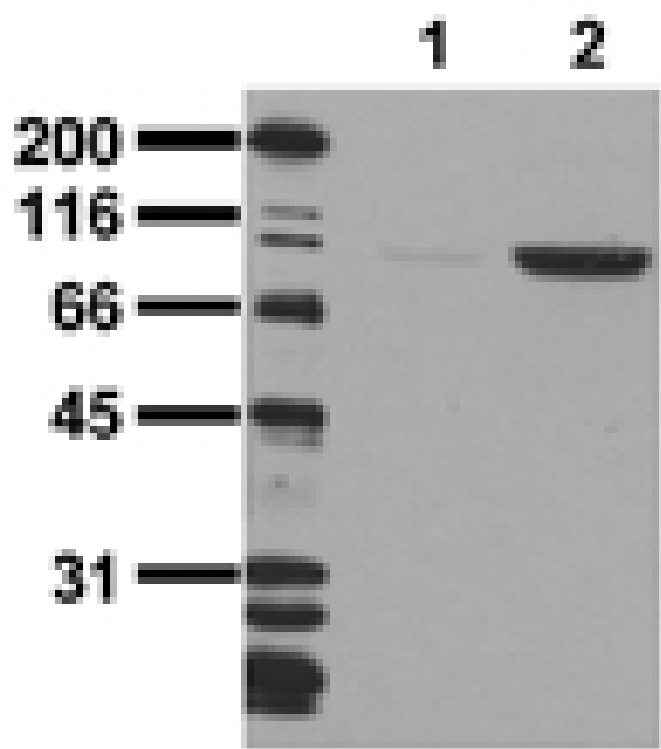
Ordering Information

[Order Online »](#)

ADI-905-645-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of untreated (1) or insulin-stimulated (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.

Isotype IgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography purified.

Recommendation 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

Worry-free Guarantee

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



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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