

Insulin receptor β monoclonal antibody (C18C4)

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](#).

Citations: 2

[View Online »](#)

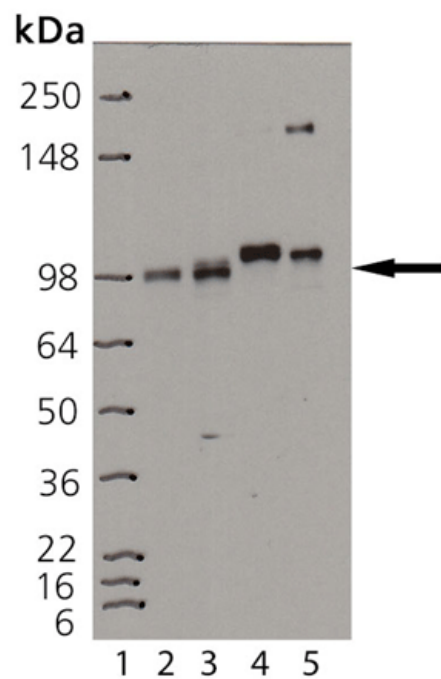
Ordering Information

[Order Online »](#)

ADI-905-683-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of ADI-905-683: Lane 1: MW Marker, Lane 2: HeLa lysate, Lane 3: MCF-7 lysate, Lane 4: PC-12 lysate, Lane 5: 3T3 lysate.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	INSR
Application	WB
Application Notes	Detects a band of ~98kDa by Western blot.
Clone	C18C4
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Recombinant cytoplasmic domain of human Insulin Receptor β .
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, colorimetric, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Bovine, Hamster, Human, Mouse, Porcine, Rabbit, Rat, Sheep
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)