

# Insulin receptor $\beta$ 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](#).

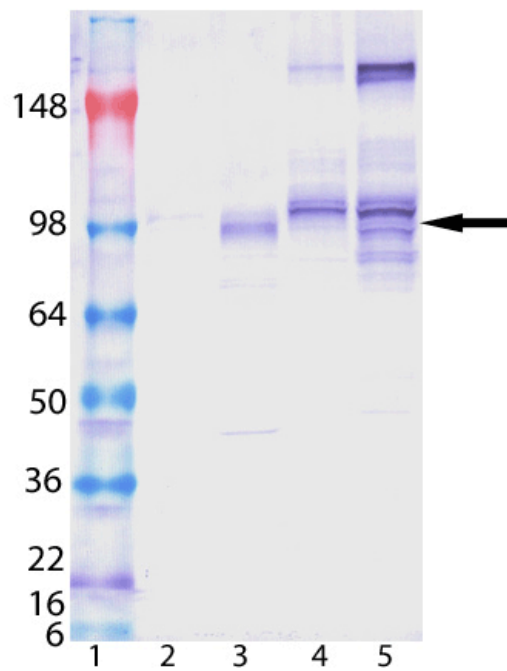
## Ordering Information

[Order Online »](#)

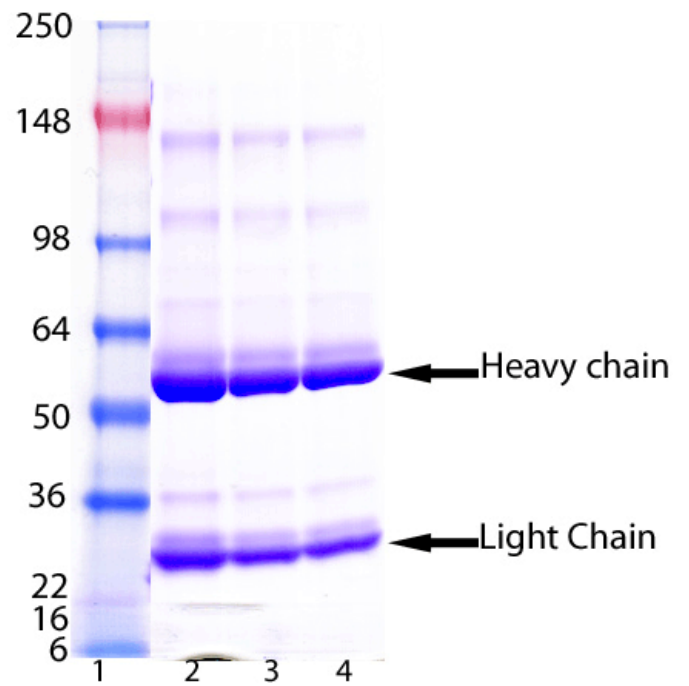
|                 |       |
|-----------------|-------|
| ENZ-ABS710-0100 | 100µg |
|-----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS710: Lane 1: MW Marker, Lane 2: HeLa lysate, Lane 3: MCF-7 lysate, Lane 4: PC-12 lysate, Lane 5: 3T3 lysate.



SDS-PAGE analysis of ENZ-ABS710: Lane 1: MW Marker, Lane 2: 10  $\mu$ g, Lane 3: 5 $\mu$ g, Lane 4: 2 $\mu$ g of Insulin receptor  $\beta$  monoclonal antibody (C18C4), resolved under reducing condition.

# 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 50% glycerol and 0.09% sodium azide.   |
| Host                                | Mouse   |
| Immunogen                           | Recombinant cytoplasmic domain of human Insulin Receptor $\beta$ .  |
| Isotype                             | IgG1  |
| Purity Detail                       | Protein G affinity purified.  |
| Recommendation Dilutions/Conditions | Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Source                              | Purified from hybridoma tissue culture supernatant.   |
| Species Reactivity                  | Bovine, Hamster, Human, Mouse, Porcine, Rabbit, Rat, Sheep  |
| UniProt ID                          | P06213  |
| Worry-free Guarantee                | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>  |



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)