

[pTyr¹³¹⁶]IGF-1 receptor monoclonal antibody (2B9)

Insulin-like growth factor-I (IGF-I) is a polypeptide hormone similar in structure to insulin. It is involved in regulation of neuronal growth and development in the central and peripheral nervous system. IGF-1 is a potent inhibitor of apoptosis in neurons, and activates the Akt signaling pathway via the IGF-1R and Insulin receptors. The IGF-1R acts as a receptor for IGF-1 and the related IGF-2, and is subject to autophosphorylation on multiple tyrosine residues that regulate receptor activation and assembly of signal transduction complexes.

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

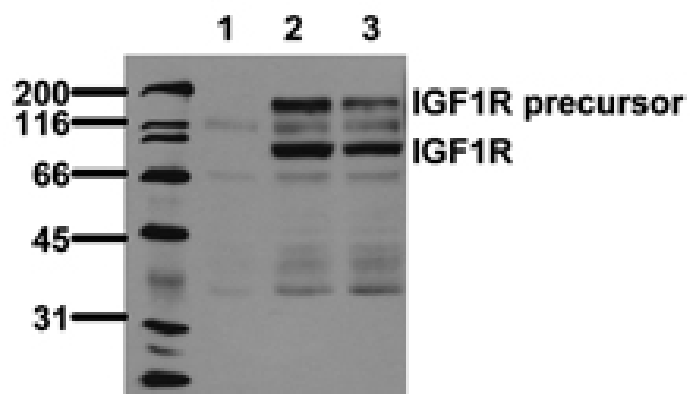
Ordering Information

[Order Online »](#)

ADI-905-643-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of untreated (1), Insulin-treated (2), or IGF1-treated (3) HeLa cell lysates, probed with IGF1 Receptor (pTyr1316) mAb (2B9).

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	Insulin-like growth factor-1 receptor, IGF-1R
Application	ELISA, WB
Application Notes	Detects a band of ~97kDa by Western blot.
Clone	2B9
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 human IGF-1 receptor phosphorylated at Tyr1316.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	ELISA (0.1µg/ml)Western Blot (1µg/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	Human
UniProt ID	P08069

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com