

Adiponectin polyclonal antibody

Adipose tissue is the source for a variety of bioactive molecules that regulate metabolism and immunity termed adipokines, most of which are elevated in the circulation of obese individuals. Adiponectin functions as an insulin sensitizing factor, resulting in reduced hepatic glucose production, increased glucose uptake in the muscle, and increased fatty acid oxidation in the liver and muscle. In contrast to most adipokines, adiponectin is detected at high levels in normal human plasma, but is often reduced in obese subjects and in those with increased insulin resistance and type 2 diabetes. Due to its suggested protective role in insulin resistance, adiponectin may be a useful pharmacological target for various metabolic diseases.

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

Citations: 1

[View Online »](#)

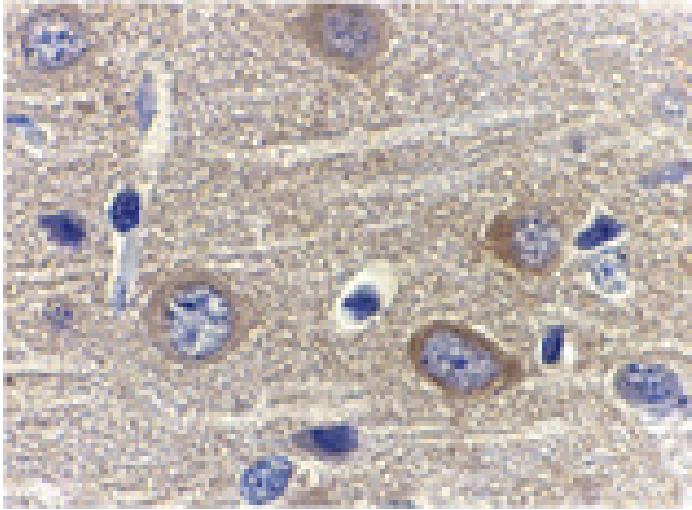
Ordering Information

[Order Online »](#)

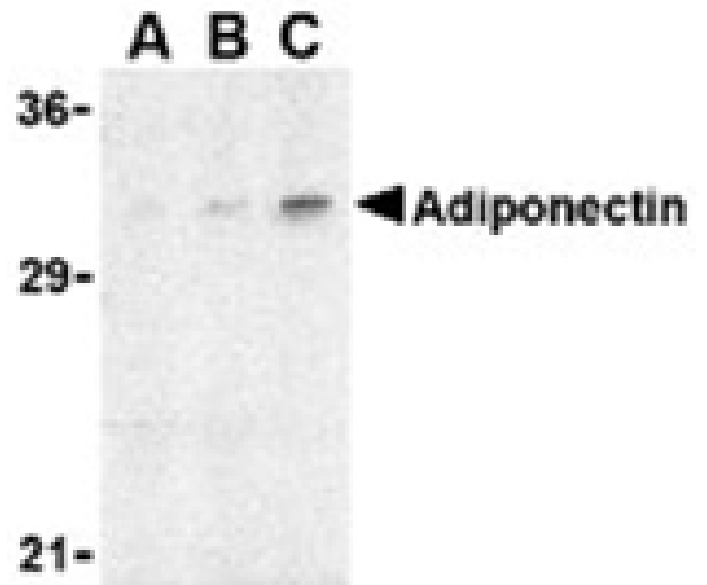
ADI-905-714-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of adiponectin in rat brain tissue with adiponectin antibody at 1 μ g/ml.



Western blot analysis of adiponectin in rat brain cell lysate with adiponectin antibody at (A) 1, (B) 2, and (C) 4 μ g/ml.

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ADIPOQ, ACDC, ACRP30, APM1, GBP28

Application IHC, WB

Application Notes Detects a band of ~32kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of human Adiponectin.

Purity Detail Peptide affinity purified.

Recommendation Immunohistochemistry (1µg/ml)Western Blot (1µg/ml)Suggested dilutions/conditions
Dilutions/Conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID Q15848

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)