

IL-6 monoclonal antibody (2C9)

IL6 is an immunoregulatory cytokine that activates a cell surface signaling assembly composed of IL6, IL6RA (IL6R; MIM 147880), and the shared signaling receptor gp130 (IL6ST; MIM 600694).

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

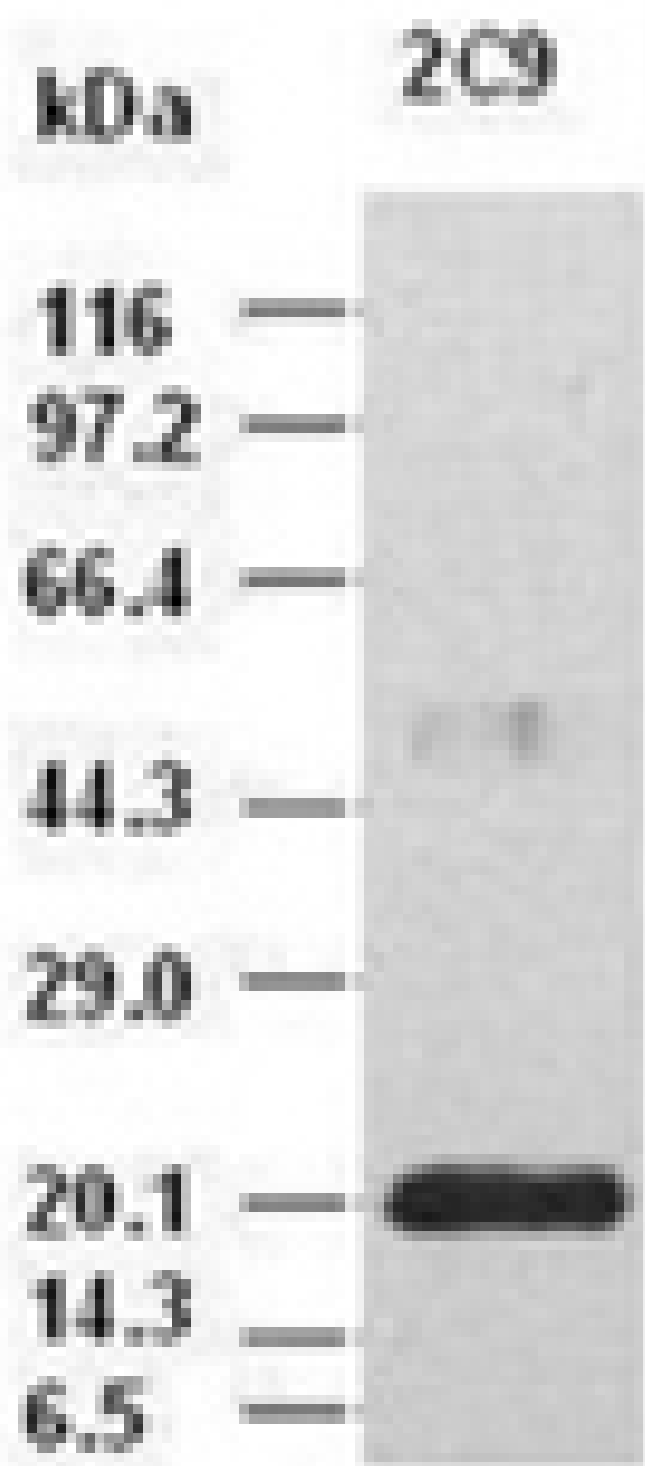
Ordering Information

[Order Online »](#)

ADI-905-884-100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Recombinant human IL-6, probed with IL-6 mAb (2C9) at 1:500000.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-6

Application WB

Application Notes Detects a band of ~24kDa (full length) by Western blot.

Clone 2C9

Formulation Liquid. In PBS, pH 7.3, containing 1% BSA, 50% glycerol and 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human IL-6 (aa 29-212).

Isotype IgG2b

Purity Detail Protein G affinity purified.

Recommendation
Dilutions/Conditions Western Blot (1:500,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P05231

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-
fr@enzolifesciences.com](mailto:info-fr@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)