# 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.

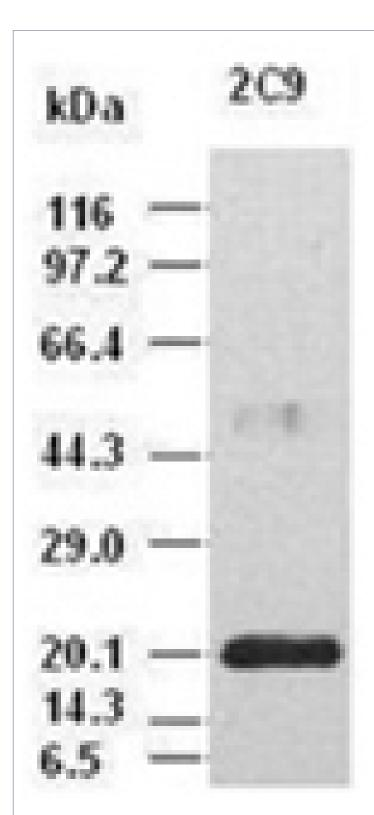
# **Ordering Information**

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).

lsotype lgG2b

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

