

# Cullin 1 monoclonal antibody (AS97.1)

The human Cullin family of genes consists of Cul-1, Cul-2, Cul-3, Cul-4a, Cul-4b, and Cul-5. Among the 6 members, Cul-1 and Cul-2 are very well characterized. Cul-2 is homologous to the *Saccharomyces cerevisiae* cdc53 gene. Studies have shown that Cul-1 forms an E3 ubiquitin-ligase, called the SCF complex (Skp1 and F-box protein such as bTrCp or FWD1), which controls the proteolysis of G1 cell cycle regulatory proteins.

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

## Ordering Information

[Order Online »](#)

ADI-KAM-CC135-D	50µg
ADI-KAM-CC135-F	200µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CUL1

**Application** WB

**Application Notes** Detects a band of ~93kDa by Western blot.

**Clone** AS97.1

**Formulation** Liquid. In PBS, pH 7.2, containing 0.1mM PMSF and 50% glycerol.

**Host** Mouse

**Immunogen** Recombinant human cullin 1.

**Isotype** IgG2a

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1:250, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from mouse ascites.

**Species Reactivity** Bovine, Human, Monkey, Mouse, Rabbit, Rat

**UniProt ID** Q13616

**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)