

# 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
-----------------	------

## 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 <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)