

| PRODUCT NAME                            |                                   |                          |
|-----------------------------------------|-----------------------------------|--------------------------|
| hSirT1 monoclonal antibody              |                                   |                          |
| Other names: hSIR2, SIR2alpha, SIR2L1   |                                   |                          |
| Cat. No. <b>C15200063</b> (MAb-063-050) | Type: Monoclonal<br>Isotype: IgG1 | Size: 50 µg/ 72 µl       |
| Lot #: 001                              | Source: Mouse                     | Concentration: 0.7 µg/µl |

**Description:** Monoclonal antibody raised in mouse against human SirT1 (NAD-dependent deacetylase sirtuin-1) using a synthetic peptide.

**Specificity:** Human: positive  
Other species: not tested

| Applications     | Suggested dilution | References |
|------------------|--------------------|------------|
| Western blotting | 1:500              | Fig 1      |

**Purity:** Protein G purified monoclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

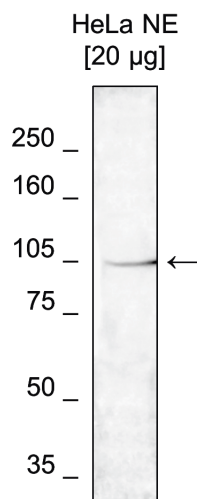
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** February 18, 2010

**Target description**

SirT1 (UniProtKB/Swiss-Prot entry Q96EB6) is a NAD-dependent deacetylase, which regulates processes such as apoptosis and muscle differentiation by the deacetylation of key regulatory proteins. Deacetylation of lysine 382 of p53/TP53 by Sirt1 impairs its ability to induce the proapoptotic program and modulate cell senescence, whereas deacetylation of TAF1B represses rDNA transcription by the RNA polymerase I. SirT1 is also involved in HES1- and HEY2-mediated transcriptional repression and inhibits skeletal muscle differentiation by deacetylating PCAF and MYOD1. SirT1 is recruited to the nuclear bodies via its interaction with PML.



**Figure 1**

**Western blot analysis using the Diagenode monoclonal antibody directed against hSirT1**

Western blot was performed using nuclear extracts from HeLa cells (HeLa NE, 20 µg) and the Diagenode monoclonal antibody against hSirT1 (cat# MAb-063-050) diluted 1:500 in TBS-Tween containing 5% skimmed milk. The molecular weight marker (in kDa) is shown on the left, the position of the protein of interest is shown on the right.