

PRODUCT NAME		
<b>hRbAp46/48 monoclonal antibody</b>		
Other names: RBBP7/4, RBBP-7/4, RBAP46/48, NURF 55		
<b>Cat. No.</b> C15200033 (MAb-033-050)	<b>Type:</b> Monoclonal <b>Isotype:</b> IgG2b	<b>Size:</b> 50 µg/ 25 µl
<b>Lot #:</b> 001	<b>Source:</b> Mouse	<b>Concentration:</b> 2.0 µg/µl

**Description:** Monoclonal antibody raised in mouse against the full length human RbAp48 protein (Retinoblastoma-binding protein p48) fused to GST. The antibody recognizes both RbAp48 and RbAp46.

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

Applications	Suggested dilution	References
Western blotting	1:1,000	Fig 1

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

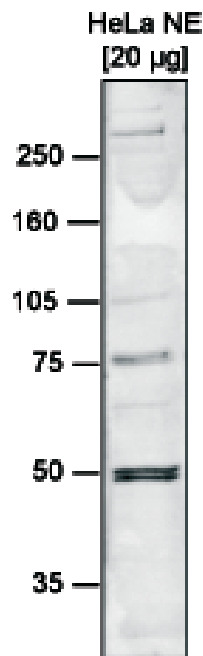
**Storage:** For long storage, store at -20°C/ -80°C. Avoid multiple freeze-thaw cycles.

**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

RbAp48 (UniProtKB/Swiss-Prot entry Q09028) is a component of several complexes which regulate chromatin metabolism. These include (1) the chromatin assembly factor 1 (CAF-1) complex, which is required for chromatin assembly following DNA replication and DNA repair; (2) the core histone deacetylase (HDAC) complex, which promotes histone deacetylation and consequent transcriptional repression; (3) the nucleosome remodelling and histone deacetylase complex (the NuRD complex), which promotes transcriptional repression by histone deacetylation and nucleosome remodeling; (4) the PRC2/EED-EZH2 complex, which promotes repression of homeotic genes during development; and (5) the NURF (nucleosome remodeling factor) complex. Therefore it may target chromatin assembly factors as well as chromatin remodeling factors and histone deacetylases to their histone substrates in a nucleosomal DNA dependent manner.



**Figure 1**

**Western blot analysis using the Diagenode monoclonal antibody directed against hRbAp46/48**

Western blot was performed using nuclear extracts from HeLa cells (HeLa NE, 20 µg) and the Diagenode monoclonal antibody against hRbAp46/48 (cat# MAb-033-050) diluted 1:1,000 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.