

## Ty1 monoclonal antibody

**Cat. No.** C15200054  
**Type:** Monoclonal ChIP-grade  
**Isotype:** IgG1  
**Source:** Mouse  
**Lot #:** 006  
**Size:** 50 µg/ 28 µl  
**Concentration:** 1.8 µg/µl

**Purity:** Protein A purified monoclonal antibody in PBS containing 0.05% azide.  
**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.

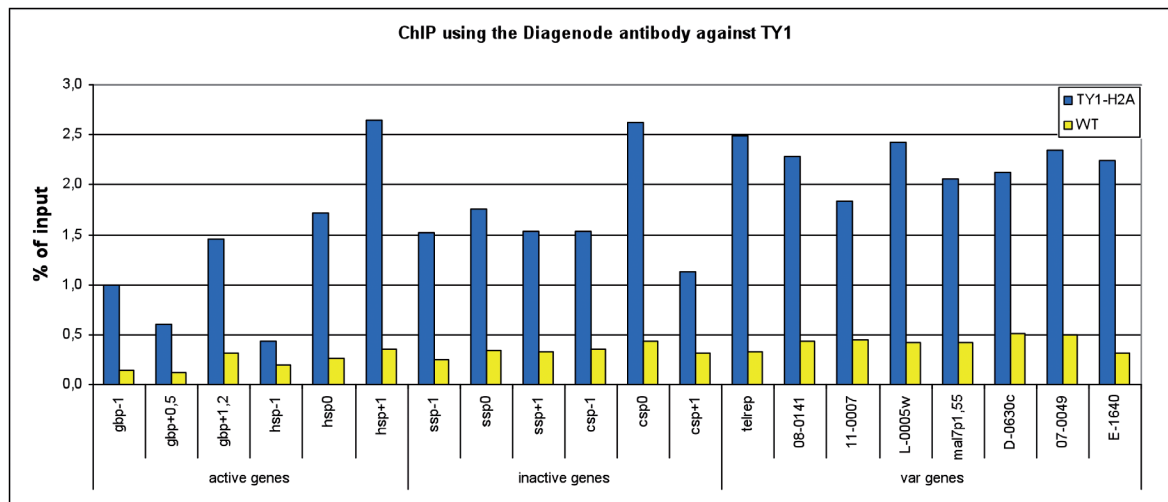
**Description:** Monoclonal antibody raised in mouse against the Ty1 tag (amino acid sequence EVHTNQDPLD).

### Applications

	Suggested dilution	Results
ChIP *	2 µg/ChIP	Fig 1
Western blotting	1:1,000 - 1:2,000	Fig 2

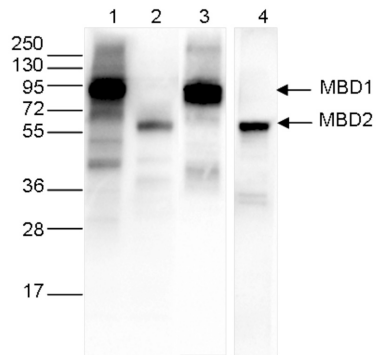
\* Please note that of the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 µg per IP.

### Results



**Figure 1. ChIP results obtained with the Diagenode monoclonal antibody against the Ty1 tag**

ChIP assays were performed using Plasmodium falciparum cells stably expressing Ty1-H2A (blue bars) and wild type Plasmodium falciparum cells (yellow bars). Primer pairs were used for active genes, inactive genes, and variable genes. Sheared chromatin from 1x10<sup>6</sup> cells and 2 µg of antibody against Ty1 (Cat. No. C15200054) were used per ChIP experiment.



**Figure 2. Western blot analysis using the Diagenode monoclonal antibody against the Ty1 tag**

Western blot was performed on 40 µg whole cell extracts from HeLa cells transfected with either Ty1-tagged MBD1 (lanes 1 and 3) or Ty1-tagged MBD2 (lane 2 and 4). The WB was performed with the Diagenode antibody against the Ty1 tag (Cat. No. C15200054) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk (lane 1 and 2) or with an MBD1 antibody (C15410078, lane 3) or an MBD2 antibody (C15410312, lane 4). The position of the Ty1-tagged MBD1 and Ty1-tagged MBD2 are shown on the right, the marker (in kDa) is shown on the left.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
 Rue Bois Saint-Jean, 3  
 4102 Seraing (Ougrée) - Belgium  
 Tel: +32 4 364 20 50  
 Fax: +32 4 364 20 51  
 orders@diagenode.com  
 info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
 Denville, NJ 07834 - USA  
 Tel: +1 862 209-4680  
 Fax: +1 862 209-4681  
 orders.na@diagenode.com  
 info.na@diagenode.com

**Last update:** November 24, 2017