

## Ty1 Antibody – ChIP Grade

**Cat. No. C15200054**

Type: Monoclonal	Specificity:
Size: 50 µg	Isotype: IgG1
Concentration: 2.2 µg/µl	Host: Mouse
Lot No.: 007	Purity: Protein A purified monoclonal antibody.
Storage buffer: PBS containing 0.05% azide	Storage conditions: 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: July 30, 2021

## Description

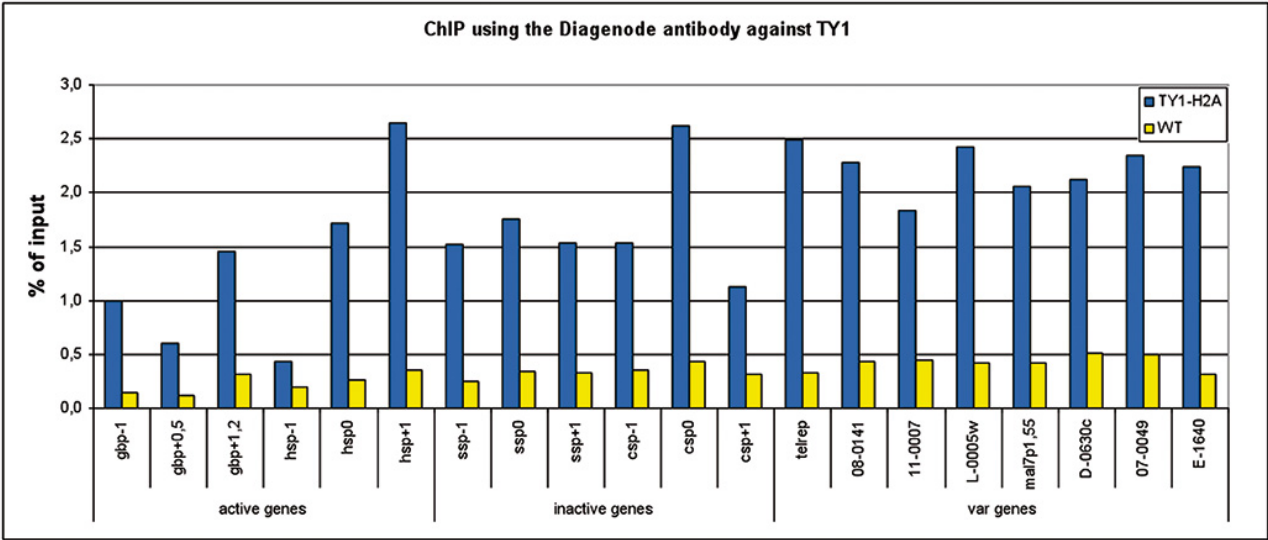
Monoclonal antibody raised in mouse against the **Ty1 tag** (amino acid sequence EVHTNQDPLD).

## Applications

Applications	Suggested dilution	References
ChIP *	2 µg/ChIP	Fig 1
Western Blotting	1:1,000 - 1:2,000	Fig 2, 3

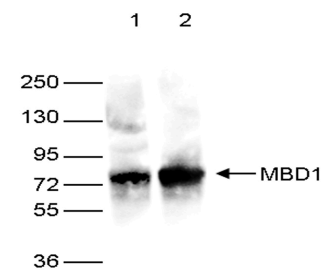
\* Please note that of the optimal antibody amount per IP should be determined by the end-user. We recommend testing 1-5 µg per IP.

Validation Data



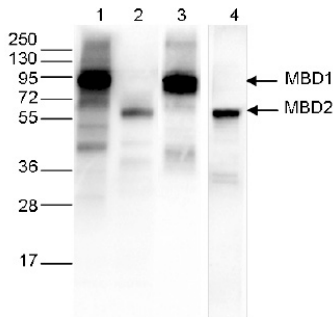
**Figure 1. ChIP results obtained with the Diagenode monoclonal antibody against the Ty1 tag**

ChIP assays were performed using *Plasmodium falciparum* cells stably expressing Ty1-H2A (blue bars) and wild type *Plasmodium falciparum* cells (brown 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 Ty1-tagged MBD1. The WB was performed with the Diagenode antibody against the Ty1 tag (cat. No. C15200054) diluted 1:2,000 in TBS-Tween containing 5% skimmed milk (lane 2) or with an MBD1 antibody (C15410078, lane 1). 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.



**Figure 3. 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.