

Ty1 monoclonal antibody

Cat. No. C15200054

Type: Monoclonal ChIP-grade

Isotype: IgG1

Source: Mouse

Lot #: 005

Size: 50 µg/ 29 µl

Concentration: 1.73 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

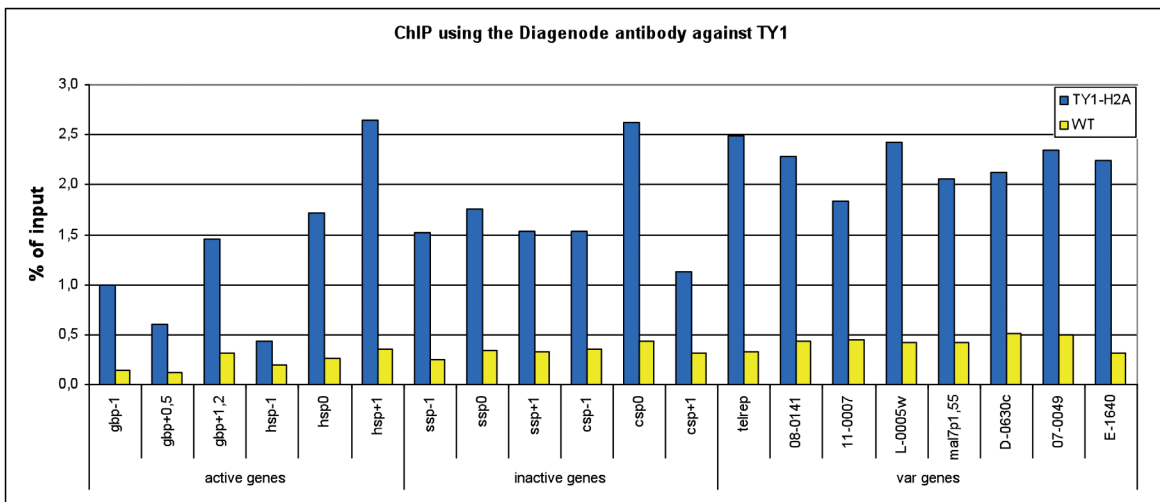


Figure 1. ChIP results obtained with the Diagenode monoclonal antibody against the Ty1 tag

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 1x10⁶ cells and 2 µg of antibody against Ty1 (Cat. No. C15200054) were used per ChIP experiment.

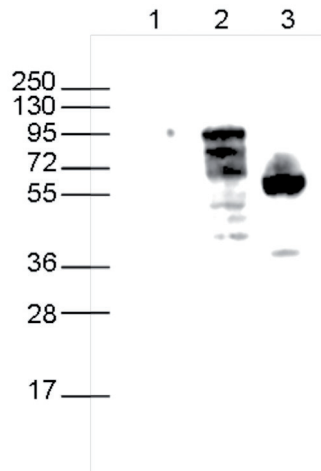


Figure 2. Western blot analysis using the Diagenode monoclonal antibody against the Ty1 tag

Western blot was performed on 25 µg whole cell extracts from HeLa cells using the antibody against the Ty1 tag (cat. No. C15200054) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. Figure 2 shows the results for untransfected cells, used as a negative control (lane 1) and for cells transfected with a Ty1-tagged MBD1 (lane 2) and a Ty1-tagged MBD2 (lane 3). 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: January 5, 2015