

VPS72 polyclonal antibody - Classic

Other name: YL1, TCFL1, CFL1

Cat. No. C15200224

Type: Monoclonal

Source: Mouse

Lot #: 001

Size: 50 µg /50 µl

Concentration: 1 µg/µl

Specificity: Human, mouse: positive / Other species: not tested

Purity: Affinity purified monoclonal antibody in PBS. Does not contain any preservative.

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 VPS72 (Vacuolar Protein Sorting 72 Homolog), using a recombinant protein.

Applications

Applications	Suggested dilution/amount	Results
Western blotting	1:1,000	Fig 1
Immunohistochemistry	1:400	Fig 2

Target description

VPS72 (UniProt/Swiss-Prot entry Q15906) is a subunit of two different multi-component protein complexes, the histone acetyltransferase complex TRRAP/TIP60 and the chromatin remodeling SRCAP-containing complex. The TRRAP/TIP60 complex acetylates nucleosomal histones playing an important role in transcriptional regulation, double strand DNA break repair and apoptosis. The SRCAP-containing complex catalyzes the exchange of histone H2A with the histone variant H2AFZ, which has a role in chromosome segregation. VPS72 may be responsible for binding H2AFZ.

Results

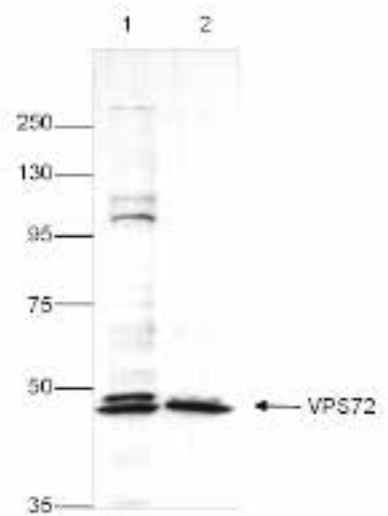


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against VPS72

Whole cell extracts from HeLa (lane 1) or MEF (lane 2) cells were analysed by Western blot using the Diagenode antibody against VPS72 (Cat. No. C15200224) diluted 1:1,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left

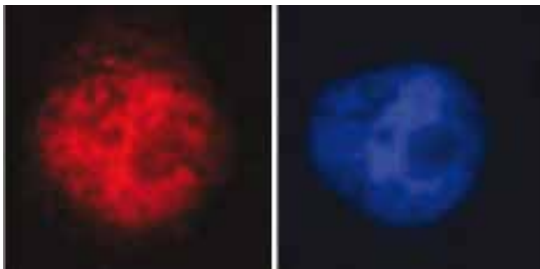


Figure 2. Immunofluorescence using the Diagenode monoclonal antibody directed against VPS72

HeLa cells were stained with the Diagenode antibody against VPS72 (Cat. No. C15200224) diluted 1:400 followed by an anti-mouse antibody conjugated to Alexa594 (left). The right figure shows staining of the nuclei with DAPI.

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: August 21, 2015