

LRWD1 polyclonal antibody

Other names: ORCA, CENP-33

Cat. No. C15310232 (CS-PA012-100)

Type: Polyclonal

Source: Rabbit

Lot #: A1016-001

Size: 100 µl

Concentration: not determined

Specificity: Human: positive / Other species: not tested

Purity: Whole antiserum from rabbit 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: Polyclonal antibody raised in rabbit against human LRWD1 (leucine-rich repeats and WD repeat domain containing 1) using two KLH-conjugated synthetic peptides containing a sequence from the N-terminal and the central region of the protein, respectively.

Applications

	Suggested dilution	Results
ELISA	1:1,000	Fig 1

*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:1,000 dilution

Target description

LRWD1 (UniProtKB/Swiss-Prot entry Q9UFC0) is a component of the origin recognition complex (ORC) required to recruit and stabilize the ORC complex to chromatin. LRWD1 is probably required for the association of ORC on chromatin during G1 to establish pre-replication complex (preRC) and to heterochromatic sites in post-replicated cells. It binds histone repressive methylation repressive marks on heterochromatin H3K9me3, H3K27me3 and H4K20me3 in a cooperative manner with DNA methylation. It is unclear whether it recognizes and binds these repressive marks by itself or needs additional factors.

Results

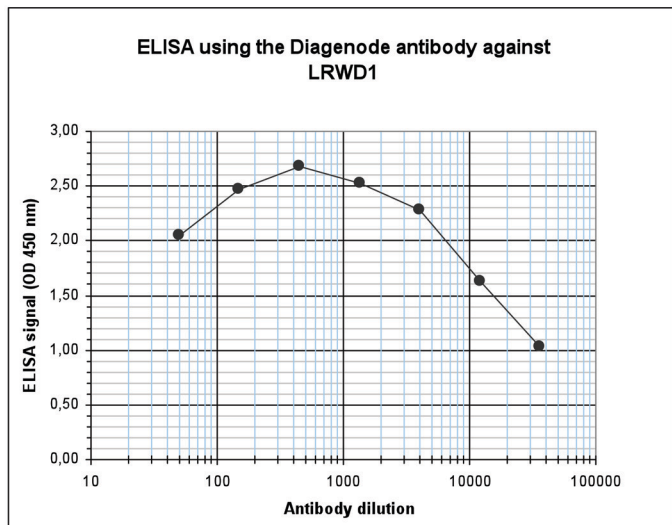


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human LRWD1 (cat. No. CS-PA012-100). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:24,900.

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: September 4, 2012