

ORC2 monoclonal antibody

Other name: ORC2L

Cat. No. C15200198

Type: Monoclonal

Isotype: IgG1k

Source: Mouse

Lot #: 001

Size: 50 µg/ 84 µl

Concentration: 0.6 µg/µl

Specificity: Human: positive
Other species: not tested

Purity: Protein G 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 human ORC2 (origin recognition complex, subunit 2), using two recombinant proteins containing the N-terminal and C-terminal part of the protein, respectively.

Applications

	Suggested dilution	Results
Western blotting	1:500	Fig 1

Target description

ORC2 (UniProt/Swiss-Prot entry Q13416) is one of the six subunits of the ORC (origin recognition complex) that binds origins of replication. The binding of ORC to DNA is ATP-dependent, but so far, specific DNA sequences that define origins of replication have not been identified. The ORC is required to assemble the pre-replication complex necessary to initiate DNA replication. The ORC is regulated in a cell cycle-dependent manner. It is assembled at the end of the anaphase of mitosis and disassembled when cells enter the S phase.

Results

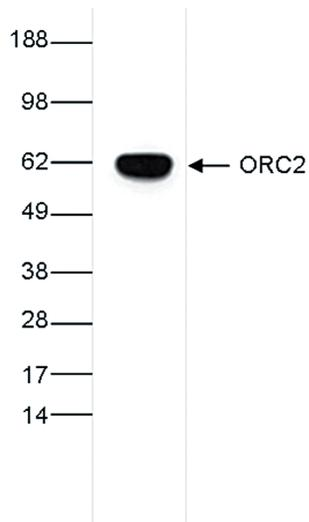


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against ORC2

RIPA cell lysates from HeLa cells (30 μ g) were analysed by Western blot using the Diagenode monoclonal antibody against ORC2 [Cat. No. C15200198] diluted 1:500 in PBS-Tween containing 10% milk. The position of the protein of interest is indicated on the right (expected size: 66 kDa); 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: October 26, 2012