

OCT4 antibody (other name: POU5F1, OTF4, OCT3, OTF3)

Cat. No. C15410305

Type: Polyclonal

Source: Rabbit

Lot: A2089P

Size: 50 µg

Concentration: 1.5 µg/µl

Specificity: Human, mouse: positive.

Other species: not tested.

Purity: Affinity purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

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 OCT4 (Octamer-Binding Protein 4), using 4 KLH-conjugated synthetic peptides containing sequences from different parts of the protein.

Applications

Applications	Suggested dilution	References
ChIP*	5 µg/ChIP	Fig 1, 2
ELISA	1:1,000-10,000	Fig 3
Western blotting	1:1,000	Fig 4

Target description

OCT4 (UniProtKB/Swiss-Prot entry Q01860) is a transcription factor that plays a key role in embryonic development and stem cell pluripotency. During embryonic development OCT4 forms a trimeric complex with SOX2 and controls the expression of a number of genes important for embryogenesis such as YES1, FGF4, UTF1 and ZFP206. Aberrant expression of OCT4 in adult tissues is associated with tumorigenesis. A translocation with the Ewing's sarcoma gene on chromosome 21 also leads to tumor formation.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. 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, 2015

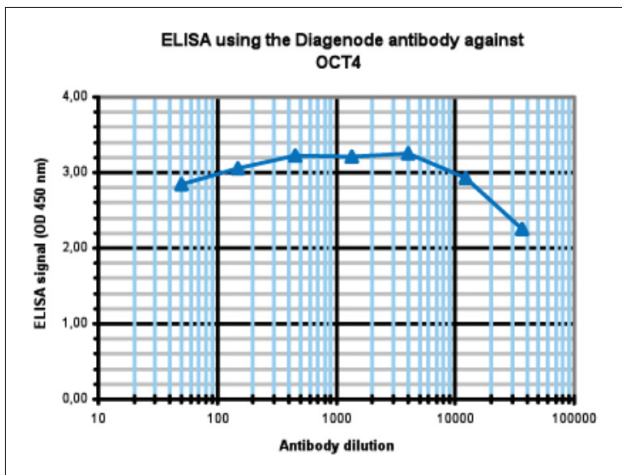


Figure 3. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against OCT4 (Cat. No. C15410305). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 2), the titer of the antibody was estimated to be 1:100,000.

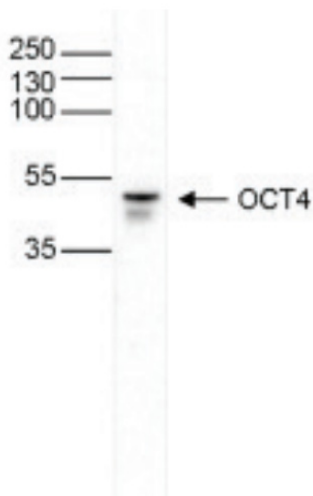


Figure 4. Western blot analysis using the Diagenode antibody directed against OCT4

Nuclear extracts (15 µg) from human embryonic stem cells were analysed by Western blot using the Diagenode antibody against OCT4 (Cat. No. C15410305) diluted 1:1,000 in TBSTween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.