

## HJURP monoclonal antibody

**Other names:** FAKTS, URLC9, HFLEG1, FLEG1

**Cat. No.** C15200228

**Type:** Monoclonal

**Isotype:** IgG1k

**Source:** Rabbit

**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 HJURP (Holliday Junction Recognition Protein), using a recombinant protein.

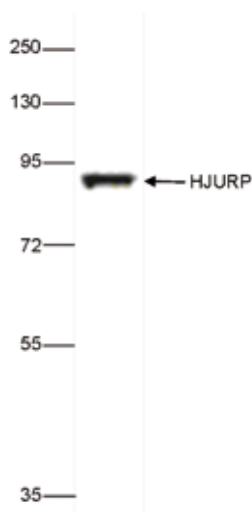
### Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1
IF	1:400	Fig 2

### Target description

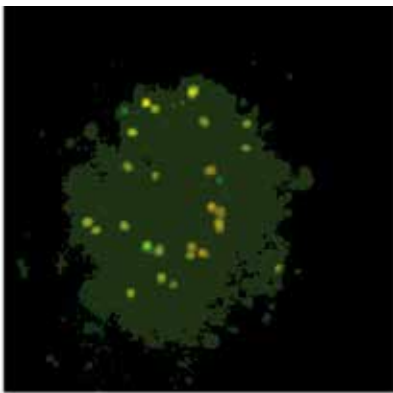
HJURP (UniProt/Swiss-Prot entry Q8NCD3) is a centromeric protein that plays a central role in the incorporation and maintenance of histone H3 variant CENPA at centromeres. It acts as a specific chaperone for CENPA and is required for the incorporation of newly synthesized CENPA molecules into the nucleosomes at replicated centromeres. HJURP prevents CENPA-H4 tetramerization and premature DNA binding by the CENPA-H4 tetramer.

## Results



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

HeLa whole cell lysates were analysed by Western blot using the Diagenode antibody against HJURP (cat. No. C15200228) 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 HJURP**

U2OS cells were stained with the Diagenode antibody against HJURP (cat. No. C15200228) diluted 1:400 and with a rabbit polyclonal CENPA antibody. Secondary antibodies were a goat anti-mouse antibody conjugated to Alexa 488 (green) and a goat anti-rabbit antibody conjugated to Alexa 647 (red).

### 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 17, 2015