

5-fC polyclonal antibody

Full name: 5-formylcytosine polyclonal antibody

Cat. No. C15310200

Type: Polyclonal

Source: Rabbit

Lot #: A1941-001

Size: 100 µl

Concentration: not determined

Specificity: Human, mouse, other (wide range): positive

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 5-formylcytosine (5-fC) conjugated to KLH.

Applications

| | Suggested dilution | Results |
|-------|--------------------|---------|
| DIP | 3 µl per IP | Fig 1 |
| ELISA | 1:10,000 | Fig 2 |

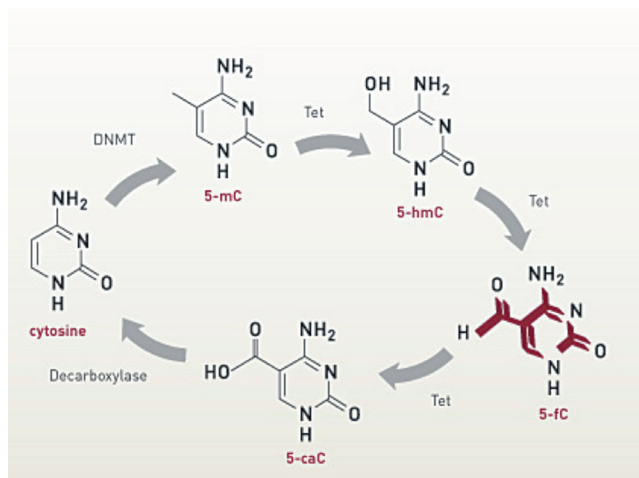
* Please note that the optimal antibody amount per ChIP should be determined by the end-user. We recommend testing 1-10 µl per IP.

Target description

Until a few years ago, 5-methylcytosine (5-mC) was the only known modification of DNA for epigenetic regulation. In 2009, however, a second methylated cytosine, 5-hydroxymethylcytosine (5-hmC) was discovered. This new modified base is generated by enzymatic conversion of 5-mC into 5-hmC by the TET family of oxygenases.

Recent results indicate that 5-hmC plays important roles distinct from 5-mC. Although its precise role has still to be shown, early evidence suggests that 5-hmC may well represent a new pathway to demethylate DNA involving a repair mechanism converting 5-hmC to cytosine. As such it may play a role in the regulation of gene activity. This pathway includes further oxidation of the hydroxymethyl group to a formyl or carboxyl group, both catalyzed by TET oxygenases. The formyl and carboxyl groups of 5-Formylcytosine (5-fC) and 5-Carboxylcytosine (5-caC) can be enzymatically removed without excision of the base.

Due to their structural similarity, the different modified cytosine analogues are difficult to discriminate. The development of highly specific affinity-based reagents, such as antibodies, appears to be the most powerful way to differentially and specifically enrich 5-mC and 5-hmC sequences. We previously released highly specific antibodies directed against 5-mC, 5-hmC and 5-caC. Now, we also present a unique rabbit polyclonal antibody against 5-fC.



Results

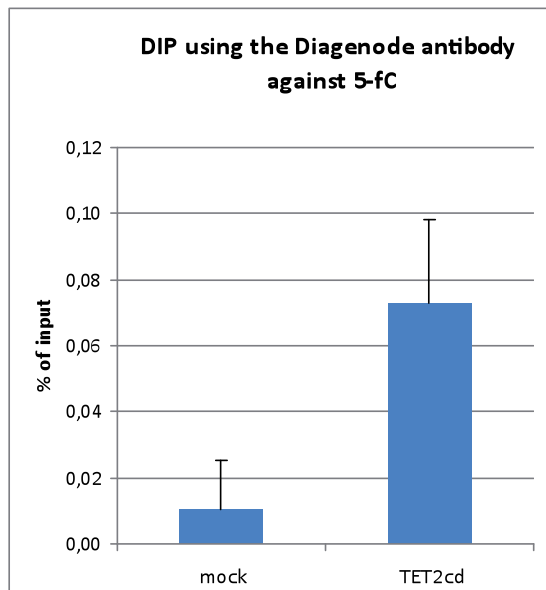


Figure 1. DIP results obtained with the Diagenode antibody directed against 5-fC

HEK293 cells were transfected with a reporter gene and hydroxymethylated in vitro with either a pCAG expression vector containing the TET2 catalytic domain (TET2cd) or a negative control pCAG vector. DIP assays were performed on 4 µg of sheared and denatured DNA using 3 µl of the Diagenode antibody against 5-fC (Cat. No. C15310200) in a total of 500 µl IP buffer. QPCR was performed with primers specific for the reporter gene. Figure 1 shows the recovery, expressed as a % of input (mean +standard deviation of 3 different experiments).

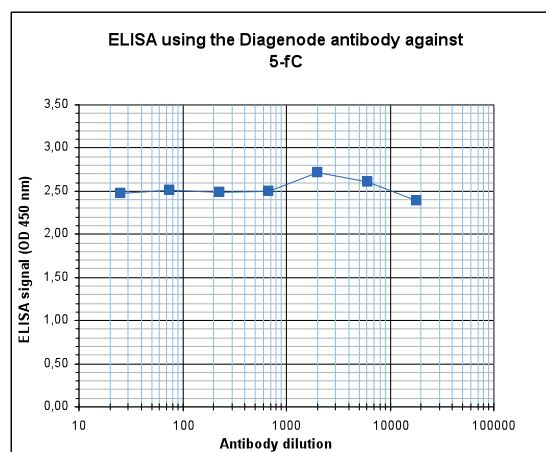


Figure 2. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against 5-fC (Cat. No. C15310200). The plates were coated with the immunogen. By plotting the absorbance against the antibody dilution (Figure 2), the titer of the antibody was estimated to be >1:100,000.

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: January 15, 2015