

TFIIF polyclonal antibody - Classic

Cat. No. C15410256

Type: Polyclonal

Source: Rabbit

Lot #: 40723

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human,mouse.

Purity: Purified by peptide affinity chromatography.

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.

Applications

	Suggested dilution*	Results
ChIP assay	5 µg/ChIP	Figure 1
Immunoprecipitation	1:100-1:500	Figure 2
Western blot	1:500-1:3000	Figure 3,4

*Optimal dilutions/concentrations should be determined by the researcher.

Results

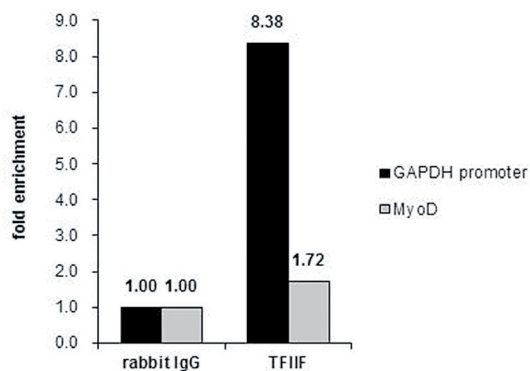


Figure 1. TFIIF ChIP results

ChIP was performed with HeLa chromatin extract and 5 µg of either control rabbit IgG or TFIIF antibody. The precipitated DNA was detected by PCR with primer set targeting to GAPDH promoter or MyoD.

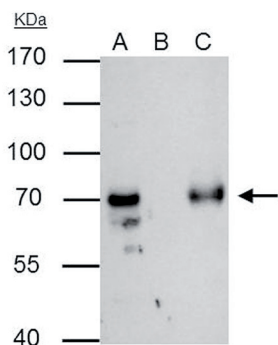


Figure 2. TFIIF IP results

TFIIF antibody immunoprecipitates TFIIF protein in IP experiments. IP Sample: 293T whole cell extract A: 30 µg whole cell extract of TFIIF protein expressing 293T cells B: Control with 2.5 µg of pre-immune rabbit IgG C: Immunoprecipitation of TFIIF with 2.5 µg of TFIIF antibody [C15410256] The immunoprecipitated TFIIF protein was detected by western blot with the TFIIF antibody diluted 1:1,000.

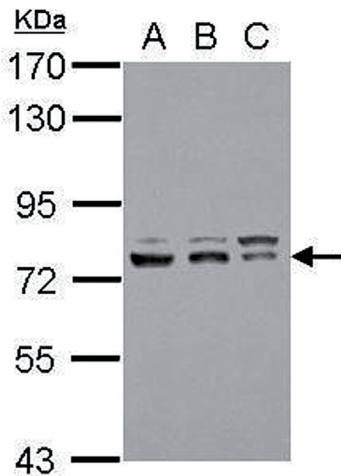


Figure 3. Western blot

Sample: 30 µg of whole cell lysate
A: 293T
B: A431
C: HeLa
7.5% SDS PAGE
TFIIIF antibody dilution 1:3,000

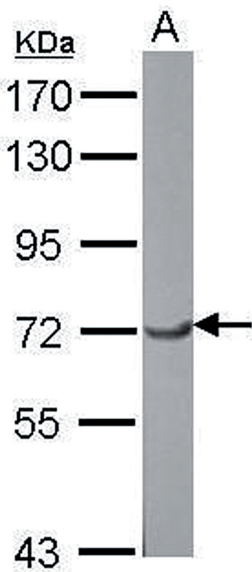


Figure 4. Western blot

Sample: 30 µg of NIH-3T3 whole cell lysate
7.5% SDS PAGE
TFIIIF antibody dilution 1:3,000

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: September 12, 2014