

## STAT5aY694p polyclonal antibody

**Other names:** STAT5, MGF

**Cat. No.** C15410329

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 100 µg/ 91 µl

**Concentration:** 1.1 µg/µl

**Specificity:** Human, mouse, rat: positive.

Other species: not tested

**Purity:** Affinity purified antibody in PBS containing 0.01% sodium 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 mouse STAT5a (Signal Transducer And Activator Of Transcription 5A) phosphorylated at Tyr694 using a KLH conjugated synthetic peptide.

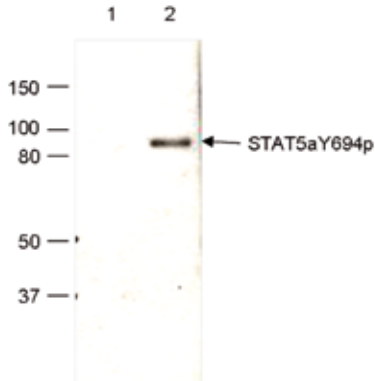
### Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1
Immunohistochemistry	1:50	Fig 2

### Target description

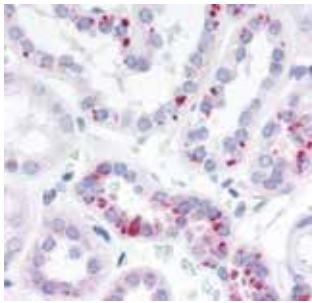
STAT5a (UniProt/Swiss-Prot entry P42229) is a member of the STAT family of transcription factors. These transcription factors are activated by phosphorylation in response to cytokines and other growth factors. Once phosphorylated the proteins are translocated from the cytoplasm to the nucleus where they activate their target genes. As such STAT proteins act as both signal transducers and transcriptional activators. STAT5 regulates the expression of milk proteins during lactation. Further, STAT5 is associated with different tumors and activation of STAT5 in myeloma and lymphoma has been shown to be essential for tumorigenesis.

## Results



**Figure 1. Western blot analysis using the Diagenode antibody directed against STAT5aY694p**

Lysates from NK92 cells treated with IL-2 (lane 2) or untreated control cells (lane 1) were analysed by Western blot using the Diagenode antibody against STAT5aY694p [Cat. No. C15410329] diluted 1:1,000. The position of the phosphorylated STAT5a is indicated on the right, the marker (in kDa) is shown on the left. The antibody recognizes the phosphorylated STAT5a but does not react with the unphosphorylated protein.



**Figure 2. Immunohistochemistry using the Diagenode antibody directed against STAT5aY694p**

Formalin fixed paraffin embedded human kidney tissue was stained with the Diagenode antibody against STAT5aY694p [Cat. No. C15410329] diluted 1:50, followed by a peroxidase labelled anti-rabbit antibody.

### 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: May 27, 2016