

STAT2Y690p polyclonal antibody

Other names: STAT113, P113, ISGF-3

Cat. No. C15410330

Type: Polyclonal

Source: Rabbit

Lot #: 001

Size: 100 µg/ 87 µl

Concentration: 1.15 µg/µl

Specificity: Human, mouse, rat., dog, pig, horse, bovine, monkey: 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 STAT2 (Signal Transducer And Activator Of Transcription 2) phosphorylated at Tyr690 using a KLH conjugated synthetic peptide.

Applications

	Suggested dilution	Results
Western blotting	1:1,000	Fig 1

Target description

STAT2 (UniProt/Swiss-Prot entry P52630) 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. STAT2 binds to the IFN stimulated response element (ISRE) to activate the transcription of interferon stimulated genes and mediates antiproliferation and antiviral responses of the cell.

Results

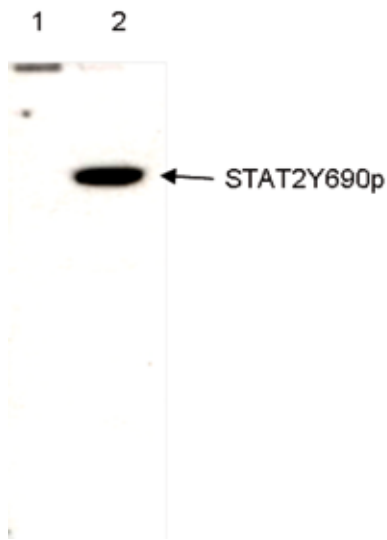


Figure 1. Western blot analysis using the Diagenode antibody directed against STAT2Y690p

Lysates from Jurkat cells treated with IFN- α (lane 2) or untreated control cells (lane 1) were analysed by Western blot using the Diagenode antibody against STAT2Y690p (cat. No. C15410330) diluted 1:1,000. The position of the phosphorylated STAT2 is indicated on the right. The antibody recognizes the phosphorylated STAT2 but does not react with the unphosphorylated protein.

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