

## SMAD3 polyclonal antibody

**Other names:** MADH3, JV15-2, HMAD-3, HSMAD3, MAD3, LDS3, HSPC193, SMAD 3

**Cat. No. C15410328**

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 100 µg/ 86 µl

**Concentration:** 1.17 µg/µl

**Specificity:** Human: 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 SMAD3 (SMAD Family Member 3) using a KLH conjugated synthetic peptide from the C-terminal par of the protein.

### Applications

	Suggested dilution	Results
Western blotting	1:500	Fig 1
Immunohistochemistry	1:400	Fig 2

### Target description

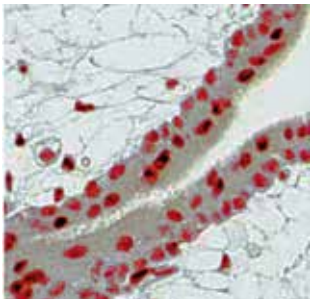
SMAD3 (UniProt/Swiss-Prot entry P84022) belongs to the SMAD ('mothers against decapentaplegic) family of proteins which are signal transducers and transcriptional modulators involved in multiple pathways. SMAD3 is activated by TGF-beta (transforming growth factor) and activin type 1 receptor kinases which phosphorylates the C-terminal serines 423 and 425. SMAD3 binds the TRE element in the promoter region of many genes that are regulated by TGF-beta and is involved in chondrogenesis, osteogenesis and wound healing. It is also thought to play a role in carcinogenesis.

## Results



**Figure 1. Western blot analysis using the Diagenode antibody directed against SMAD3**

Lysates from different cells treated with TGFb (+) or untreated control cells (-) were analysed by Western blot using the Diagenode antibody against SMAD3 (Cat. No. C15410328) diluted 1:500. The position of SMAD3 is indicated on the right. The antibody recognizes both the phosphorylated and the unphosphorylated SMAD3.



**Figure 2. Immunohistochemistry using the Diagenode antibody directed against SMAD3**

Formalin fixed paraffin embedded cancerous human liver tissue was stained with the Diagenode antibody against SMAD3 (Cat. No. C15410328) diluted 1:400, followed by a peroxidase labelled anti-rabbit antibody. The presence of the SMAD3 protein is shown by the grey color; red indicated the nuclei.

### 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