

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.

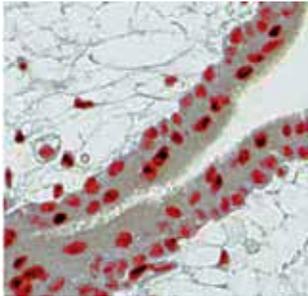


Figure 2. Immunohistochemistry using the Diagenode antibody directed against SMAD3

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