

SMAD3S423/425p polyclonal antibody

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

Cat. No. C15410327

Type: Polyclonal

Source: Rabbit

Lot #: 001

Size: 100 µg/ 100 µl

Concentration: 1 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) phosphorylated at Ser423 and Ser425, using a KLH conjugated synthetic peptide.

Applications

| | Suggested dilution | Results |
|----------------------|--------------------|---------|
| Western blotting | 1:2,000 | 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

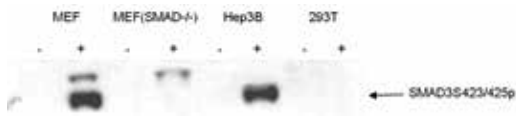


Figure 1. Western blot analysis using the Diagenode antibody directed against SMAD3S423/425p

Lysates from different cells treated with TGFb (+) or untreated control cells (-) were analysed by Western blot using the Diagenode antibody against SMAD3S423/425p (cat. No. C15410327) diluted 1:2,000. The position of the phosphorylated SMAD3 is indicated on the right. Figure 1 clearly shows the antibody recognizes the phosphorylated SMAD3 whereas it has no reactivity towards the unphosphorylated SMAD3.

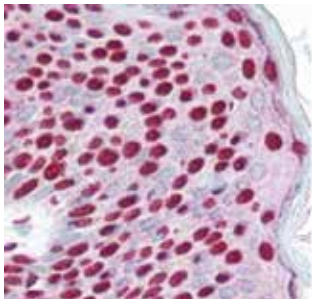


Figure 2. Immunohistochemistry using the Diagenode antibody directed against SMAD3S423/425p

Formalin fixed paraffin embedded epidermal keratinocytes were stained with the Diagenode antibody against SMAD3S423/425p (cat. No. C15410327) diluted 1:400, followed by a peroxidase labelled anti-rabbit antibody. Figure 2 shows nuclear staining of SMAD3 (red).

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