

TRIM33 recombinant protein

Other names: TIF1G, TIF1gamma, TF1G, RFG7, ECTO, PTC7

Cat. No. C23010110

Source: SF9 cells

Lot #: 001

Size: 10 µg/ 20 µl

Concentration: 0.5 µg/µl

Specificity: Human

Purity: Purified using FPLC, >95% purity as determined by SDS-PAGE.

Storage buffer: 20 mM Tris-Cl pH 7.6, 1 mM EDTA, 0.15 M NaCl, 10% glycerol, 0.5 mM PMSF and 1 mM DTT.

Storage: Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Full length recombinant human N-terminal Flag tagged TRIM33 (Tripartite Motif Containing 33) protein, produced in SF9 insect cells.

Protein description

TRIM33 (UniProt/Swiss-Prot entry Q9UPN9) is a member of the tripartite motif (TRIM) family. The TRIM motif includes three zinc-binding domains - a RING, a B-box type 1 and a B-box type 2 - and a coiled-coil region. TRIM33 is an E3 ubiquitin-protein ligase, which promotes SMAD4 ubiquitination, nuclear exclusion and degradation via the ubiquitin proteasome pathway. It may act as a transcriptional corepressor and inhibits the transcriptional response to the TGF-beta/BMP signaling cascade. However, molecules that interact with this protein have not yet been identified..TRIM33 also plays a role in the control of cell proliferation. Its association with SMAD2 and SMAD3 stimulates erythroid differentiation of hematopoietic stem/progenitor. Defects in TRIM24 cause thyroid papillary carcinoma.

Quality control

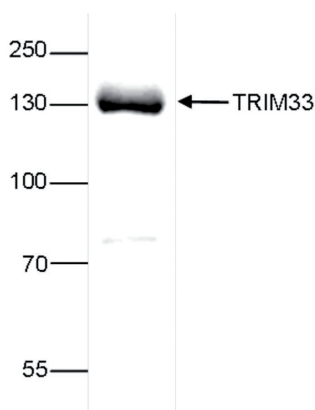


Figure 1.

SDS page of the TRIM33 recombinant protein. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

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: August 28, 2015