

TRIM24 recombinant protein

Other names: TIF1A, TIF1, TF1A, HTIF1, RNF82, PTC6

Cat. No. C23010109

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 TRIM24 (Tripartite Motif Containing 24) protein, produced in SF9 insect cells.

Protein description

TRIM24 (UniProt/Swiss-Prot entry O15164) 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. It acts as a transcriptional coactivator through its interaction with several nuclear receptors, including the estrogen, retinoic acid, and vitamin D3 receptors, and coactivators thereby modulating the transcription of target genes. TRIM24 is thought to associate with chromatin. This association depends on histone H3 modifications with the highest affinity for histone H3 that is both unmodified at 'Lys-4' and acetylated at 'Lys-23'. Defects in TRIM24 cause thyroid papillary carcinoma.

Quality control

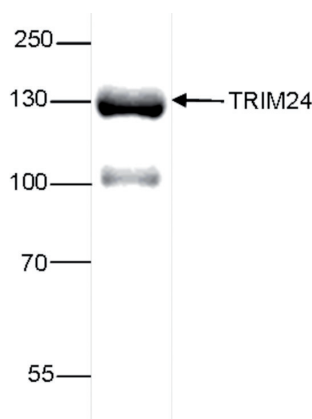


Figure 1.

SDS page of the TRIM24 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 25, 2015