

SMRT / NCoR2 antibody (pAb)

Catalog Nos: 39845, 39846

RRID: AB_2793365

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.89 μ g/ μ l

Molecular Weight: 460 kDa

Background: SMRT / NCoR2 (Silencing Mediator of Retinoic acid and Thyroid hormone receptor, Nuclear receptor CoRepressor 2, SMAP270, CTG26) is a co-repressor of transcription that associates with a subset of nuclear receptors including RAR, PR, GR and TR in the absence of ligand. It is part of a large co-repressor protein complex that includes SIN3A/B and histone deacetylases HDAC1 and HDAC2.

Immunogen: This SMRT / NCoR2 antibody was raised against a peptide within the N-terminal region of SMRT / NCoR2.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

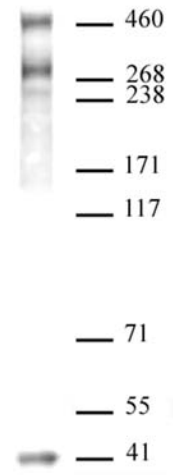
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



SMRT / NCoR2 antibody (pAb) tested by Western blot.

Whole cell extract of HeLa cells (40 μ g per lane) probed with SMRT / NCoR2 antibody (pAb) (1:500 dilution).