

## SMRT / NCoR2 antibody (mAb)

**Catalog Nos:** 61105, 61106

**RRID:** AB\_2615017

**Clone:** 5.10

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 300 kDa

**Background:** **SMRT / NCoR2** (Nuclear receptor corepressor 2, CTG repeat protein 26, SMAP270, Silencing mediator of retinoic acid and thyroid hormone receptor, Thyroid and retinoic-acid-receptor-associated corepressor, TRAC) is a transcriptional corepressor for nuclear receptors. **SMRT / NCoR2** has been shown to play a role in regulating immune response, mediating DNA repair, and in neuronal differentiation and cardiac development. **SMRT / NCoR2** is part of a large corepressor complex that contains SIN3A/B and histone deacetylases HDAC1 and HDAC2. It also includes TBL1, TBLR1, and GPS2 that help stabilize the overall structure of the complex. SMRT / NCoR2 mediates the transcriptional repression activity of nuclear receptors by promoting histone deacetylation and chromatin condensation, thus preventing access of the basal transcription machinery.

**Immunogen:** This SMRT / NCoR2 antibody was raised against a recombinant protein corresponding to amino acids 2004-2310 of human 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:

**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.