

NR0B1 antibody (pAb)

Catalog Nos: 61661, 61662

RRID: AB_2793724

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 60 kDa

Background: NR0B1 (Nuclear Receptor Subfamily 0, Group B, Member 1) is an orphan nuclear receptor and a component of a cascade required for the development of the hypothalamic-pituitary-adrenal-gonadal axis. NR0B1 inhibits the transcriptional activity of other nuclear receptors through heterodimeric interactions. It may also have a role in embryonic development and in the maintenance of embryonic stem cell pluripotency as it up-regulates Oct-4 expression ES cells. Defects in NR0B1 are a cause of dosage-sensitive sex reversal (DSSR) and X-linked adrenal hypoplasia congenital (XL-AHC). DSSR results from a duplication of the region of the X-chromosome containing NR0B1, resulting in males with two copies of NR0B1, which inhibits normal testis formation, and thus XY individuals develop as females. XL-AHC is a developmental disorder of the adrenal gland, characterized by a structural disorganization of the glands. Hypogonadotropic hypogonadism (HHG) is frequently associated with this disorder. HHG is a condition resulting from or characterized by abnormally decreased gonadal function, with retardation of growth and sexual development.

Immunogen: This antibody was raised against recombinant human NR0B1 protein.

Buffer: Purified IgG in PBS with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

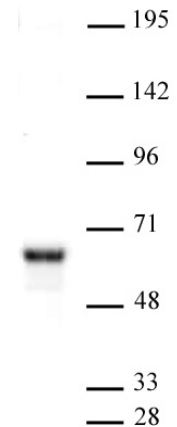
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 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.



NR0B1 antibody (pAb) tested by Western blot.

Detection of NR0B1 by Western blot. Jurkat whole cell extract (20 μ g) probed with NR0B1 antibody at a 1:500 dilution.