

DAX-1 / NR0B1 antibody (mAb)

Catalog No: 39983

RRID: AB_2793417

Clone: 1DA-2F4

Isotype: IgG1, k

Application(s): ICC, IF, IHC, WB

Reactivity: Human, Mouse, Rat

Quantity: 100 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 52 kDa

Background: DAX-1 (DSS-AHC critical region on the X chromosome protein 1, 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. DAX-1 / 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 DAX-1 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 DAX-1, resulting in males with two copies of DAX-1, 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 DAX-1 / NR0B1 antibody was raised against a peptide corresponding to amino acids 135-166 of human DAX-1 / NR0B1.

Buffer: Purified IgG in 60 mM Sodium Citrate, 105 mM Tris-HCl, 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.