

## CHD1 antibody (mAb)

**Catalog Nos:** 39989, 39990

**RRID:** AB\_2793420

**Clone:** 2F11

**Isotype:** IgG2a

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

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Rat

**Concentration:** 1 µg/µl

**Molecular Weight:** 210 kDa

**Background:** CHD1 (Chromo-ATPase/Helicase-DNA-binding protein 1 ) is a member of the CHD family of ATP-dependent chromatin remodeling factors implicated in RNA polymerase II transcriptional regulation. They are characterized by the presence of a double chromodomains (Chromatin Organization Modifier) and SNF2-related helicase/ATPase domains. The CHD1/CHD2 sub-family of CHD proteins also contain a C-terminal DNA binding domain that recognizes AT-rich sequences. CHD proteins alter gene expression possibly by modification of chromatin structure thus altering access of the transcriptional apparatus to its chromosomal DNA template. CHD1 has been shown to interact with histone H3 trimethylated at lysine 4 (H3K4Me3) and also is required for the maintenance of stem cell pluripotency.

**Immunogen:** This CHD1 antibody was raised against a peptide corresponding to the C-terminus of mouse CHD1.

**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.