

TWIST antibody (mAb)

Catalog Nos: 61097, 61098

RRID: AB_2793505

Isotype: IgG1

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

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 21 kDa

Background: TWIST is a regulator of transcription that serves as a master regulator of **embryonic morphogenesis**. Through its bHLH motif it interacts with many different transcription factors to affect multiple pathways of gene expression. TWIST plays a pivotal role in **mesodermal**, myoblast, and osteoblast differentiation. TWIST is involved in the **epithelial–mesenchymal transition (EMT)**, a period during embryonic development characterized by loss of cell adhesion and increased cell mobility, essential events for mesoderm and neural tube formation. The EMT shares many features of metastatic cancer, and elevated levels of TWIST are observed in many types of cancer. Mutations in the gene encoding TWIST have been associated with Saethre-Chotzen syndrome, an inherited developmental disorder characterized by skull deformations, limb abnormalities, and facial dysmorphism.

Immunogen: This TWIST antibody was raised against a recombinant protein corresponding to full-length mouse TWIST.

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.