

SNAI1 / Snail antibody (mAb)

Catalog Nos: 61367, 61368

RRID: AB_2793608

Clone: 9H2

Isotype: IgG2a

Application(s): IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 32 kDa

Background: **SNAI1 / Snail** (snail homolog 1 (Drosophila)) is involved in the epithelial to mesenchymal transition (EMT) and formation and maintenance of embryonic mesoderm (By similarity). Binds to 3 E-boxes of the E-cadherin gene promoter and represses its transcription.

Immunogen: This antibody was raised against a full-length recombinant protein corresponding to human SNAI1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91255.

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.