

Goosecoid antibody (pAb)

Catalog Nos: 61121, 61122

RRID: AB_2793515

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 40 kDa

Background: Goosecoid (Gsc) is a homeodomain transcription factor important for vertebrate development and organogenesis. First identified in *Xenopus*, it is expressed in the area of the Spemann organizer, a region within the embryo that control development and differentiation. Goosecoid regulates cell migration from the Spemann organizer and is overexpressed in a majority of human breast tumors. It is also involved in craniofacial and ribcage development.

Immunogen: This Goosecoid antibody was raised against a peptide in the N-terminal region of human homeobox protein Goosecoid.

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

Application Notes:

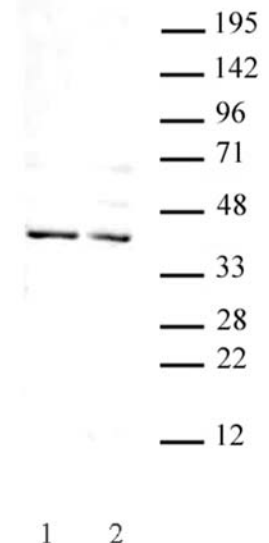
Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

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.



Goosecoid antibody (pAb) tested by Western blot.

Detection of Goosecoid by Western blot analysis. Lane 1: Nuclear extract of HT-29 cells (30 µg). Lane 2: Nuclear extract of SW48 cells (30 µg). Both were probed with Goosecoid antibody (pAb) at a 1:500 dilution.