

Geminin antibody (pAb)

Catalog Nos: 61135, 61136

RRID: AB_2793521

Application(s): IF, IP, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Isotype: Serum

Molecular Weight: 33 kDa

Background: Geminin is a crucial cell cycle regulatory protein that controls DNA replication and centrosome function. Geminin inhibits DNA replication by preventing the binding of the MCM complex into the pre-replication complex. Geminin accumulates during S, G2, and M phases, and disappears at the time of the metaphase-anaphase transition. Geminin functions as a homotetramer and interacts with DNA replication licensing factor CDT1. The interaction between Geminin and CDT1 is crucial to restrict DNA replication to once per cell cycle. Geminin is required for the formation of pluripotent cells in the developing embryo and essential for pluripotency in the embryonic stem cells. In ESCs, Geminin remains during G1 and sustains core pluripotency factors (Oct4, Sox2, and Nanog) by antagonizing the action of the chromatin remodeling protein BRG1.

Immunogen: This Geminin antibody was raised against full-length recombinant human Geminin protein.

Buffer: Rabbit serum containing 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.