

DPPA2 antibody (pAb)

Catalog Nos: 65619, 65919, 65620

RRID: AB_3216317

Application(s): WB

Reactivity: Human

Quantities: 100 µg, 50 µg, 10 µg

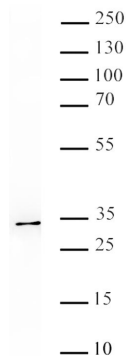
Purification: Protein A Chromatography

Host: Rabbit

Isotype: IgG

Molecular Weight: 32 kDa

Background: DPPA2 and DPPA4 (developmental pluripotency associated 2 and 4) are small heterodimerizing nuclear proteins linked in the biological processes of transcription and transcription regulation. The encoded proteins have a scaffold-attachment factor A/B, acinus and PIAS (SAP) domain that binds DNA and is thought to modify chromatin. The nuclear factor encoded by the gene is involved in the maintenance of pluripotency in stem cells and is found to be essential for embryogenesis. **DPPA2 and DPPA4** are required to maintain H3K4me3 and H3K27me3 at important bivalent promoters depicted in pluripotent cells by low H3K4me3 levels and initiating but not elongating RNA polymerase II. This gene is thought to play an important role in tumor progression because when mutated it is highly expressed in embryonal carcinomas, pluripotent germ cell tumors, and other cancers.



DPPA2 antibody (pAb) tested by Western blot.

Nterra nuclear extract (20 µg per lane) was probed with DPPA2 antibody (0.5 µg/mL).

Immunogen: This antibody was raised against a peptide corresponding to DPPA2.

Buffer: PBS containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB: 0.5µg/mL-2µg/mL

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.