

WDR4 antibody (pAb)

Catalog Nos: 65113, 69113, 65114**RRID:** AB_3216374**Application(s):** WB**Quantities:** 100 µg, 50 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 45 kDa

Background: **WDR4 (WD Repeat-Containing Protein 4)**, also known as **TRM82**, is a member of the WD repeat protein family. WD repeats are minimally conserved regions of approximately 40 amino acids typically bracketed by gly-his and trp-aspartate (GH-WD), which may facilitate formation of heterotrimeric or multiprotein complexes. **WDR4** is the non-catalytic subunit of N7-methylguanosine (m7G). **WDR4** and METTL1 form a complex required for tRNA (guanine(46)-N(7))-methyltransferase activity. This complex catalyzes tRNA m7G methylation to enhance translation efficiency.

Immunogen: This antibody was raised against a human peptide containing WD repeat domain 4.

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

Application Notes:

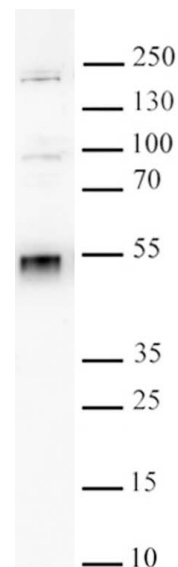
Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml

*Note: For optimal results, we recommend the addition of 0.05% Tween-20 to all blocking solutions to reduce background in Western blot.

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.

**WDR4 antibody (pAb) tested by Western blot**

20 µg of HeLa whole cell extract was run on SDS-PAGE and probed with antibody at a dilution of 2 µg/mL.