

## Drosha antibody (pAb)

**Catalog Nos:** 39783, 39784

**RRID:** AB\_2793342

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 0.44  $\mu$ g/ $\mu$ l

**Molecular Weight:** 160 kDa

**Background:** Drosha is an RNase III enzyme that is involved in the process of generating microRNA (miRNA). miRNA interacts with the RNA-induced silencing complex (RISC) to cleave mRNA and thus downregulate gene expression. Drosha is a double stranded RNase and part of the Microprocessor complex. Other RNase III enzymes include Dicer and Argonaute. Drosha initiates processing of the miRNA precursor (pri-miRNA) in the nucleus, which is then exported to the cytoplasm and further cleaved by Dicer into a mature miRNA.

**Immunogen:** This Drosha antibody was raised against a peptide within the C-terminal region of human Drosha.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

IP: 10  $\mu$ l per IP

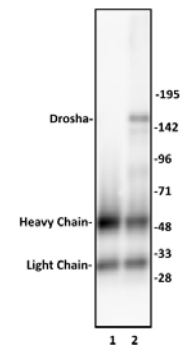
WB\*: 1:500 - 1:2,000 dilution

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.

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

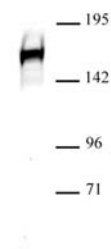
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not 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.



### Drosha antibody (pAb) tested by Immunoprecipitation.

10  $\mu$ l of Drosha antibody was used to immunoprecipitate Drosha from 250  $\mu$ g of HeLa nuclear cell extract (lane 2). 10  $\mu$ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the Drosha antibody at a dilution of 1:500.



### Drosha antibody (pAb) tested by Western blot.

HeLa cell nuclear extract (40  $\mu$ g) probed with Drosha antibody at a dilution of 1:500.