

DNA-RNA Hybrid antibody (mAb)

Catalog Nos: 65683, 65983, 65684**Clone:** S9.6**Application(s):** ChIP, ChIP-Seq, DB, EMSA, ICC, IF, IP**Reactivity:** Wide Range Predicted**Quantities:** 100 µg, 50 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Isotype:** IgG2a, k**Concentration:** 1 µg/µl**Molecular Weight:** N/A

Background: DNA-RNA hybrids are a natural occurrence within eukaryotic cells, with levels of these hybrids increasing at sites with high transcriptional activity, such as during transcription initiation, repression, and elongation. Because RNA-DNA hybrids influence genomic instability, the S9.6 antibody is a useful reagent to help study the consequences of R-loops and lesions formed by these hybrids during DNA replication or other cellular processes. In addition, the S9.6 antibody is effective in recognizing RNA-DNA hybridization for microarray studies.

Immunogen:

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

Application Notes:

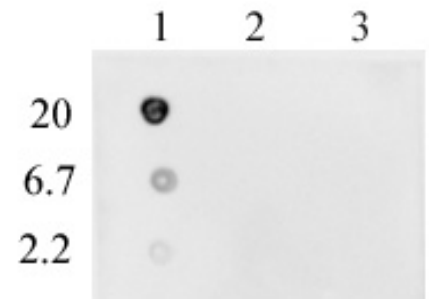
Applications Validated by Active Motif:

DB: 1 µg/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.

**DNA-RNA Hybrid antibody tested by dot blot.**

23-mer DNA, RNA or DNA-RNA Hybrid samples were spotted (indicated in ng on the left) on to a positively charged nylon membrane and blotted with DNA-RNA Hybrid antibody at 1 µg/ml dilution. Lane 1: Double-stranded DNA-RNA hybrid. Lane 2: Single-stranded DNA oligo. Lane 3: Single-stranded RNA oligo.