

# DNA-RNA immunoprecipitation (DRIP) protocol

Created July 7, 2019

## DRIP protocol

A step-by-step DRIP protocol, including R-loop preparation and associated reagents.

DNA-RNA immunoprecipitation (DRIP) uses the S9.6 anti-RNA-DNA hybrid antibody to capture RNA-DNA hybrids along chromosomes. DRIP is typically followed by mapping DNA fragments on a few loci or even across the whole genome with qPCR, microarray hybridization, or deep sequencing.

Thanks to [Professor Frédéric Chédin's lab at UC Davis](#) for providing us with this protocol.

## R-Loop preparation

### Reagents

25 mM rNTP stock (NEB N0466S) - dilute to 2.5 mM rNTP for experiment

T3 RNA Polymerase, 50 U/μL (NEB M0378S)

10 x RNAPol Reaction Buffer (NEB M0378S)

1M DTT (from frozen stock, made in house)

2.5% Tween-20 (diluted in water, made in house)

pCALM3\_2 plasmid (pCALM3\_2 carries an R-loop forming portion of the human *CALM3* gene)

RNase A, 10 mg/mL (DNase free) – dilute to 1.0 mg/mL for experiment

RNase H, 5 U/μL (NEB M0297S)

Proteinase K, 10 mg/mL (Sigma P2308)

DuRed (nucleic acid dye), 1000X (Brigen D009)

ApaLI, 2500 units (NEB R0507S)

## Protocol

1. Mix:

pCALM3\_2 2 µg  
10x buffer 10 µL  
1M DTT 2 µL  
2.5% Tween-20 2 µL  
2.5 mM rNTP 10 µL  
H<sub>2</sub>O to 99.4 µL total

2. Initiate reaction by adding 0.6 µL of T3 RNA Polymerase, mix gently, and split into two reactions (50 µL each). Incubate at 37°C for 10 minutes.
3. Inactivate enzyme by adding 1 µL of 10 mM EDTA.
4. To one sample, add 10 µL of 0.1 mg/mL RNase A, and the other (negative control) add 10 µL 0.1 mg/mL RNase A and 4 µL RNase H. Incubate for 30 minutes at 37°C.
5. Add 4 µL of Proteinase K and incubate for 30 minutes at 37°C.
6. Cleaned up on Axygen PCR purification kit and eluted in 50 µL ddH<sub>2</sub>O, separately.
7. Each sample (2 samples totally) split into two tubes, put one tube on ice, the other one was used for the following digestion.

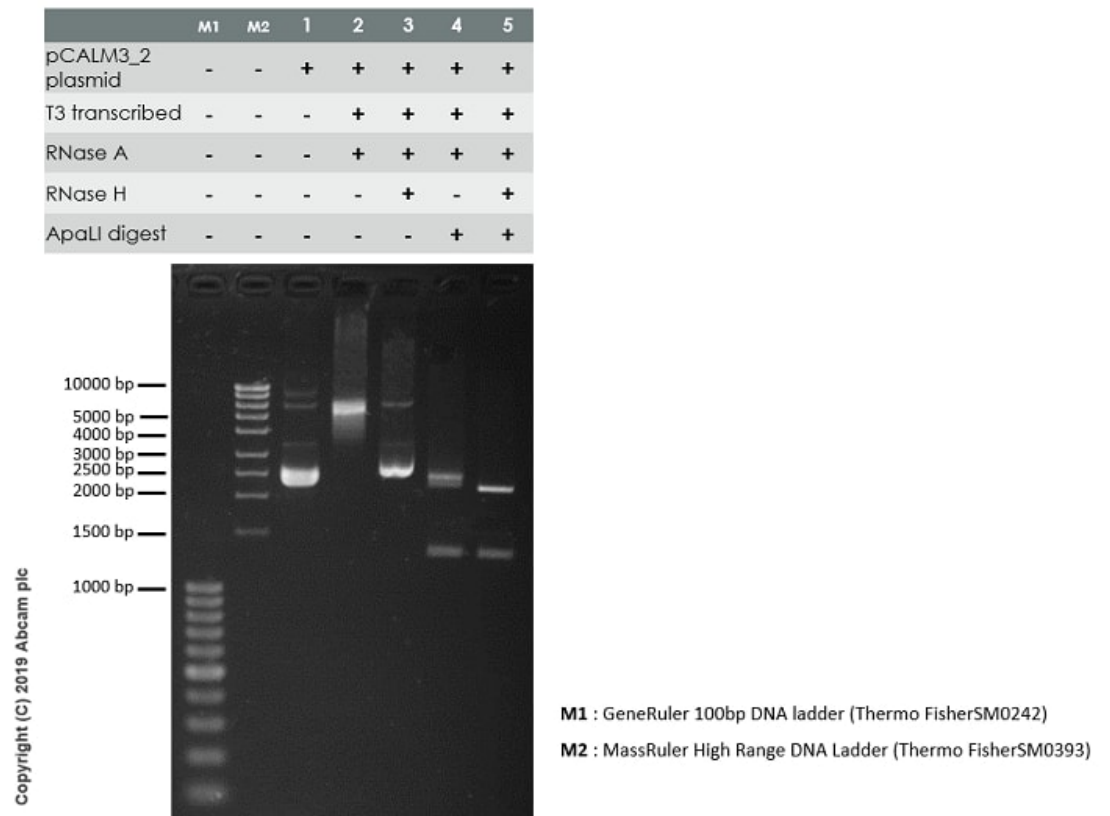
8. Digest:

25 µL DNA  
5 µL 10X Cutsmart buffer NEB  
1 µL LI ApaLI NEB  
19 µL ddH<sub>2</sub>O  
50 µL total

9. 37°C 2 hours.
10. Cleaned up on Axygen PCR purification kit and eluted in 25 µL ddH<sub>2</sub>O. Four samples (2–5) plus one pCALM3\_2 plasmid (1) were obtained for DRIP experiment.
11. In order to verify that R-loop formation did occur, load ~200 ng on a 0.9% 1x TBE gel **WITHOUT** nucleic acid dye and run at 90V for 60 min.

Use Glycerol at 10% final as a loading dye. Post-stain with DuRed (nucleic acid dye).

R-loop formation causes a characteristic shift in mobility compared to untranscribed or RNase H-treated samples. Results are shown below (Figure 1).



**Figure 1.** Transcription from pCALM3\_2 to generate R-loops. Each digestion reaction was run on an agarose gel. pCALM3\_2 carries a portion of the human *CALM3* gene that forms R-loops when transcribed with the T3 RNA polymerase. Treatment with RNase A (digests single-stranded RNA) does not affect the R-loops structure (lane 2) whereas treatment with RNase H (digests RNA in DNA-RNA hybrids) destroy R-loops structure (lane 3). pCALM3\_2 plasmid can be digested by ApaLI restriction enzyme without affecting the R-loop structures.

# DNA-RNA hybrid

## Immunoprecipitation by using antibodies pre-immobilized on beads

### Reagents

PBS (phosphate buffer)

Triton X100

ssDNA (Salmon sperm single strand DNA)

[Recombinant Anti-DNA:RNA hybrid antibody \[S9.6\] \(ab234957\)](#)

Isotype antibody (Mouse (G3A1) mAb IgG1 Isotype Control #5415, 2.5 mg/ml, Mouse IgG1, kappa)

EDTA (Ethylene Diamine Tetraacetic Acid)

2.5% Tween-20 (diluted in water, made in house)

ApaLI (NEB R0507S)

RNase A, 10 mg/mL (DNase free) – dilute to 1.0 mg/mL for experiment

RNase H, 5 U/μL (NEB M0297S)

Proteinase K, 10 mg/mL (Sigma P2308)

DuRed (nucleic acid dye), 1000X (Brigen D009)

Qiagen PCR purification kit

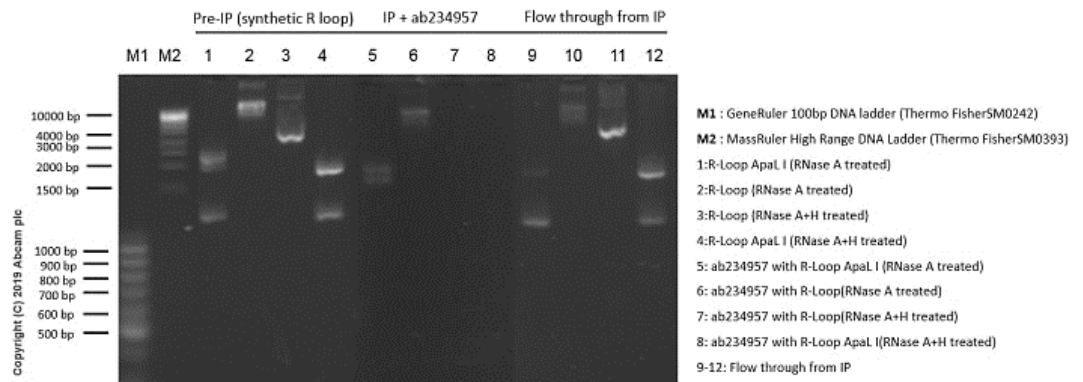
## Protocol

### Antibodies pre-immobilized on beads

1. Prepare eight tubes of Protein A beads.
2. 100  $\mu$ L protein A beads each tube washed twice in 1 mL of 1X PBS, 0.1% Triton X100, centrifuge 1 min at 3,600 rpm 4°C, carefully aspirate the supernatant.
3. Each tube resuspended with 1 mL 1X PBS, 0.1% Triton X100 and 7.5  $\mu$ g ssDNA (Salmon sperm single strand DNA/20  $\mu$ L beads), shake gently 10 mins at room temperature.  
  
Then centrifuge 1 min at 3,600 rpm at 4°C, aspirate the supernatant. Wash once in 1 mL of 1X PBS, 0.1% Triton X100, centrifuge 1 min at 3,600 rpm 4°C, carefully aspirate the supernatant.
4. 5  $\mu$ L S9.6 antibody (1mg/mL) (test antibody 5  $\mu$ g, add to protein A) added to four tubes as positive control and 5  $\mu$ L isotype antibody (5  $\mu$ g) added to rest of four tubes as a negative control, make up the samples to 1 mL with 1X PBS, 0.1% Triton X100.  
  
Shake gently 10 mins at room temperature.
5. Wash twice with 1 mL 1X PBS, 0.1% Triton X100, centrifuge 1min at 3,600 rpm 4°C, aspirate the supernatant.
6. Resuspend each tube in 100  $\mu$ L PBS, 0.1% Triton X100 and add 1  $\mu$ L 0.5M EDTA.

### Add DNA

7. Remove 5  $\mu$ L of each input R-loop (RNase A treated), R-loop (RNase A+H treated), ApaLI digested R-loop (RNase A treated), ApaLI digested R-loop (RNase A+H treated) for gel analysis as control.
8. Add 35  $\mu$ L of R-loop (RNase A treated), R-loop (RNase A+H treated), ApaLI digested R-loop (RNase A treated), ApaLI digested R-loop (RNase A+H treated) to two tubes (S9.6, isotype), respectively.
9. Rotate gently for 2 hours at 4°C.
10. Centrifuge 1 min at 3,600 rpm 4°C, remove all depleted supernatants and retain for electrophoresis.
11. Wash three times in 1X PBS, 0.1% Triton X100, centrifuge 1 min at 3,600 rpm 4°C and aspirate the supernatant.
12. Add 50  $\mu$ L elution buffer + 5  $\mu$ L proteinase K, then shake in 1,400 rpm 30 mins 50°C, centrifuge 1 min at 13,000 rpm at room temperature. Collect supernatant.
13. Clean up by Qiagen PCR purification kit and eluted in 30  $\mu$ L.



**Figure 2.** DNA-RNA hybrid Immunoprecipitation (DRIP) data. pCALM3\_2 was used to generate R-loops. S9.6 ([ab234957](#)) immunoprecipitates R-loops in the presence or absence of prior digestion by ApaLI, which does not affect R-loop structure. Prior treatment with RNase A (digests single-stranded RNA) does not affect the IP signal whereas prior treatment with RNase H (digests RNA in DNA-RNA hybrids) eliminates the signal.