

On-Column DNase Digestion Kit

(Cat# K2066-50, 50 preparations, Store at -20 °C)

I. Introduction:

BioVision's On-Column DNase Digestion Kit provides the user a quick and efficient way to remove contaminating genomic DNA from RNA preparations. The kit is designed to be used in conjunction with BioVision's Mammalian RNA Isolation Kit (Cat# K2065) or One Prep DNA/RNA Purification Kit (Cat# K2058) to clean-up or eliminate DNA from purified RNA or from enzymatic reactions. The purified RNA is suitable for numerous downstream applications including RT-PCR. Additionally, the digestion protocol can be used with other commercially available RNA extraction kits.

II. Application:

- Removal of trace amounts of DNA from RNA being purified using spin-column technology.

III. Key Features:

- On-column digestion of DNA
- DNA digestion prior to RNA clean-up

IV. Sample Types:

- Up to 1×10^7 (cultured cells) or 20 mg (tissue)
- Enzymatic RNA reaction

V. Kit Contents:

Components	50 preparations	Cap Code	Part Number
RNase-free DNase	100 μ l	Red	K2066-50-1
10X RNase-free DNase Buffer	500 μ l	Yellow	K2066-50-2
RNase-free Water	5 ml	NM	K2066-50-3

VI. User Supplied Reagents and Equipment:

- Ethanol (96-100%)
- Barrier pipette tips
- 1.5 ml RNase free microcentrifuge tubes
- Benchtop centrifuge
- One Prep DNA/RNA Purification Kit (**BioVision Cat# K2058**)
- Mammalian RNA Isolation Kit (**BioVision Cat# K2065**)

VII. Shipment and Storage:

The kit is shipped at 4 °C but should be stored at -20 °C. Aliquots of RNase-free DNase should be stored at -20 °C but can be stored at 4 °C for up to 1 month. Avoid repeated freeze-thaw cycles.

VIII. DNA Digestion Protocols:

On-Column DNA Digestion (after step 4 of Total RNA Purification Protocol under section VIII of BV Cat# K2058 or # K2065))

- Prepare the **DNA digestion master mix** containing 5 μ l of 10X RNase-free DNase Buffer, 43 μ l of RNase-free water and 2 μ l of RNase-free DNase. Mix well. **Note:** When performing multiple digestions, mix enough reagents for the number of digestions.
- Add 350 μ l of RW1 Buffer to the RNA Column and centrifuge at 10,000 x g and RT for 30 sec. Discard the flow through.
- Load 50 μ l of DNA digestion master mix in the center of the RNA Column. Incubate at room temperature (RT) for 15 min to remove any trace amounts of genomic DNA.
- Add 350 μ l RW1 Buffer to the RNA Column and centrifuge at 10,000 x g and RT for 30 sec. Discard the flow-through
- Return back to steps 6-10 of the **Total RNA Purification Protocol** (BV # K2058 or BV # K2065).

RNA Clean-Up Protocol (use either with BV # K2058 or BV # K2065)

- Add RNase-free water to the enzymatic reaction to adjust the final volume to 100 μ l.
- Add 350 μ l of RNA-Lysis buffer to the diluted RNA and mix well.
- Add 250 μ l of Ethanol (96-100%) and mix well.
- Load the 700 μ l of the solution onto an RNA column and centrifuge at 10,000 x g and RT for 30 sec. Discard the flow-through.
Perform Optional On-Column DNA digestion, if required.
- Add 500 μ l RW2 (or W2) Buffer (*add Ethanol to RW2 or W2 Buffer before use) to the column and centrifuge at 10,000 x g and RT for 30 sec. Discard the flow-through.
- Add another 500 μ l RW2 (or W2) Buffer to the column and centrifuge at 10,000 x g at RT for 1 min.
- Put the column in a new collection tube. Centrifuge the column at max speed for 1 min.
- Transfer the column to an RNase-free 1.5 ml tube and add 50-100 μ l RNase free water to the center of the column. Centrifuge at 10,000 x g and RT for 1 min to elute the RNA.
- Store the RNA solution at -80 °C or -20 °C for short term storage. To quantify using spectroscopy, RNA should be diluted in 10 mM Tris-HCl, pH 7.5.

Alternate RNA Clean-Up and DNA Digestion Protocol

1. Add 10 μ l 10X RNase-free DNase Buffer and 2 μ l RNase-free DNase to the RNA enzymatic reaction. Adjust the volume to 100 μ l with RNase-free water and mix well.
2. Incubate at 37 $^{\circ}$ C for 15-30 min.
3. Add 350 μ l of RNA-Lysis Buffer and mix well.
4. Add 250 μ l of Ethanol (96-100%) and mix well.
5. Load 700 μ l of the solution onto an RNA Column and centrifuge at 10,000 x g and RT for 30 sec. Discard the flow-through.
6. Return to steps 6-10 the **Total RNA Purification Protocol** (BV # K2065 or BV # K2058) or steps 5-9 of the **RNA Clean-Up Protocol** above.

I. General Troubleshooting Guide:

Low A_{260}/A_{280} ratio	• Protein contamination	• Perform all the wash steps, particularly with RW1 buffer.
	• Low yield	• Ensure that the sample is diluted in 10 mM Tris, pH 7.5, and the dilution solution is used as blank.
Low Yield	• Degraded sample	• Purification should be performed using fresh samples or well-preserved samples.
	• Poor column binding	• Make sure Ethanol is added and the sample is mixed well before binding to the column.
	• Exceeding the column capacity	• It is best to split the samples when the RNA yield is > 50 μ g.
	• RNase contamination	• Be careful in preventing RNase contamination.

IX. Functional Test Data:

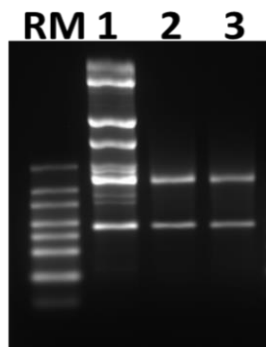


Figure 1: 5 μ g Jurkat cell RNA mixed with Hind III-digested λ DNA was purified. 1. RNA without DNase treatment. 2. DNase treatment prior to the RNA Clean-Up Protocol. 3. RNA Clean-Up Protocol with optional On-Column DNA digestion step. RM: Riboruler marker. All samples were incubated at RT for 30 min for DNase treatment. 10 μ l was loaded per lane.

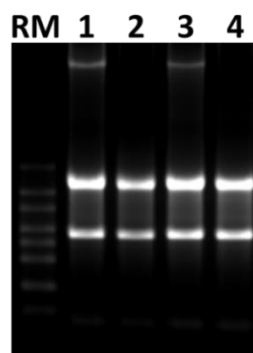


Figure 2: RNA was purified from Jurkat cells using BV Cat# K2065. Lanes 1 & 3. RNA without DNase treatment. Lanes 2 & 4. RNA with DNase treatment. RM: Riboruler marker. All samples were incubated at RT for 15 min for DNase treatment. 10 μ l of elution was loaded per lane.

X. Related Products:

- One Prep DNA/RNA Purification Kit (Cat# K2058)
- Mammalian RNA Isolation Kit (Cat# K2065)
- RNAkeepTM Solution (Cat# M1241)
- EasyRNATM Bacterial RNA Kit (Cat# K1351)
- EasyRNATM Blood RNA Mini Kit (Cat# K1373)
- EasyRNATM Plant RNA Mini Kit (Cat# K1374)
- Yeast RNA Mini Kit (Cat# K1418)
- EasyRNATM Fungal RNA Kit (Cat# K1419)

FOR RESEARCH USE ONLY! Not to be used on humans.