

PureLink[™] Pro 96 Viral RNA/DNA Kit

For rapid, efficient purification of viral nucleic acids from cell-free samples

Catalog no. 12280-096A

Rev. Date: 7 May 2009 Part no. 100007167

MAN0001672

User Manual

Contents

Contents	iii
Experienced Users Procedure	v
Kit Contents and Storage	vii
Accessory Products	viii
Introduction	1
About the Kit	1
Methods	4
Before Starting	4
Experimental Overview	
General Guidelines	٩
Preparing Lysates	
	10
Preparing Lysates	10 11
Preparing Lysates Purification Using Centrifugation	10 11 12
Preparing Lysates Purification Using Centrifugation Purification Using a Vacuum Manifold	10 11 12 14
Preparing Lysates Purification Using Centrifugation Purification Using a Vacuum Manifold Example of Expected Results	10 11 12 14 15
Preparing Lysates Purification Using Centrifugation Purification Using a Vacuum Manifold Example of Expected Results Troubleshooting	
Preparing Lysates Purification Using Centrifugation Purification Using a Vacuum Manifold Example of Expected Results Troubleshooting Appendix	

Experienced Users Procedure

Introduction

This quick reference sheet is included for experienced users of the PureLink[™] Pro 96 Viral RNA/DNA Kit. If you are a first time user, follow the detailed protocol in this manual.

Step	Action
Preparing Lysates	The lysate preparation protocol is described below for $200 \ \mu l$ of starting material.
	1. Add 25 μl Proteinase K to each well.
	2. Add 200 μ l of cell-free sample to the well.
	Note: If processing <200 µl sample, adjust the final volume of the sample to 200 µl using PBS (phosphate buffered saline) or 0.9% NaCl.
	 Add 200 μl Lysis Buffer (containing 5.6 μg Carrier RNA). Mix by pipetting up and down 3–5 times.
	4. Incubate at 56℃ for 15 minutes. Briefly centrifuge.
	5. Add 250 µl 96–100% ethanol to each well and mix by pipetting up and down 3–5 times.
	6. Incubate the lysate for 5 minutes at room temperature.
Purification	Perform all centrifuge steps at 25°C.
Procedure with	 Transfer each lysate to the PureLink[™] 96 Well Viral Filter Plate. Cover unused wells with foil tape.
Centrifu- gation	 Place the Viral Filter Plate onto a 96-well, Deep-Well Block and centrifuge at ≥2,100 × g for 1–2 minutes.
	3. Remove the flow-through and place the Viral Filter Plate back on the Deep-Well Block.
	4. Add 500 μl Wash Buffer II with ethanol.
	5. Centrifuge 5 minutes at ≥2,100 × g until liquid passes through the Viral Filter Plate.
	6. Remove the flow-through and place the Viral Filter Plate back on the Deep-Well Block.
	7. Repeat Steps 4–6 above.
	8. Centrifuge the Viral Filter Plate at $\geq 2,100 \times g$ for 10 minutes.
	 Add 100 µl RNase-free water and incubate the plate for 1 minute at room temperature.
	 Place the Viral Filter Plate onto a clean Deep-Well Block. Centrifuge at ≥2,100 × g for 1–2 minutes.
	11. Use the RNA/DNA for the desired downstream application. To store the purified RNA/DNA, cover the wells with foil tape, and store at -80°C.

Experienced Users Procedure, Continued

Step	Action
Purification Procedure	Prepare lysate as described on previous page. Perform all steps at 25°C.
with Vacuum	 Place a PureLink[™] 96 Well Viral RNA/DNA Filter Plate on top of the manifold.
	2. Transfer lysates from Preparing Lysates, Step 6 to the Viral Filter Plate. Cover unused wells with foil tape.
	3. Apply vacuum for 2 minutes. Release the vacuum.
	4. Add 500 μl Wash Buffer II with ethanol.
	5. Apply vacuum 2 minutes. Release the vacuum.
	6. Repeat Steps 4 and 5 above.
	7. Tap the Viral Filter Plate on a stack of paper towels and pat dry.
	8. Return the Viral Filter Plate to the vacuum manifold and apply vacuum for 10 minutes. Release the vacuum.
	9. Replace the waste collection tray with a clean Deep-Well Block and place the Viral Filter Plate on top of the Deep-Well Block.
	 Add 150 μl RNase-free water and incubate the plate for 1 minute at room temperature.
	11. Apply vacuum for 2 minutes. Release vacuum.
	12. The RNA/DNA is eluted into the Deep-Well Block in a volume of 110–130 μl. Use the RNA/DNA for the desired downstream application. To store the purified RNA/DNA, cover the wells with foil tape, and store at –80°C.

Kit Contents and Storage

Shipping and Storage		All components of the PureLink [™] Pro 96 Viral RNA/DNA Kit is shipped at room temperature. Upon receipt, store all kit components except Carrier RNA at room temperature.		
		Store Carrier RNA at -20°C.		
	Note: The Proteinase K solution is stable for 1 ye room temperature. For long-term storage (>1 ye temperature is >25°C, store the Proteinase K solu		if room	
		The components and amounts included in the I Pro 96 Viral RNA/DNA Kit is listed below.	PureLink™	
		Sufficient reagents are in the kit to perform 384 reactions.	(4 × 96)	
	Component		Amount	
	PureLink [™] P	ro 96 Viral Lysis Buffer (L22)	100 ml	
	PureLink [™] P	ro 96 Wash Buffer II (5X)	2 × 87.5 ml	
	Proteinase K	(20 mg/ml) in storage buffer (proprietary)	10 ml	
	Carrier RNA	(lyophilized)	2.2 mg	
	PureLink [™] P	ro 96 RNase-free Water	75 ml	
	PureLink [™] 96	5 Well Viral RNA/DNA Filter Plate	4	
	96 Deep-Well Block		2×6/bag	
	Foil Tape		$3 \times 4/bag$	
			Ū.	

Accessory Products

Additional Products

The following products are also available from Invitrogen. For details on these products, visit www.invitrogen.com or contact Technical Support (page 17).

Product	Quantity	Catalog no.
RNase AWAY®	250 ml	10328-011
RNaseZAP [®]	250 ml	AM9780
PureLink™ Foil Tape	50 pieces	12261-012
Viral Lysis Buffer (L22)	500 ml	12282-500
PureLink [™] 96 RNA Wash Buffer II (5X)	2 x 1 L	12173-033
EveryPrep [™] Universal Vacuum Manifold	1 manifold	K211101
UltraPure [™] DEPC-treated Water	1 L	750023
UltraPure [™] DNase/RNase-Free Distilled Water	500 ml	10977-015
Yeast tRNA	25 mg	15401-011
Phosphate Buffered Saline (PBS), 1X	500 ml	10010-023
SuperScript [™] III One-Step RT-PCR System with Platinum [®] <i>Taq</i> DNA Polymerase	100 reactions	12574-026
RNA UltraSense™ One-Step Quantitative RT-PCR System	100 reactions	11732-927
Platinum [®] Quantitative PCR SuperMix- UDG w/ROX	100 reactions	11743-100

Introduction

About the Kit

Introduction	The PureLink [™] Pro 96 Viral RNA/DNA Kit provides a rapid and efficient method to simultaneously purify viral RNA/DNA from fresh or frozen cell-free biological fluids (plasma, serum, cerebrospinal fluid) and cell culture supernatants.
	The PureLink [™] Pro 96 Viral Kit is specifically designed for obtaining the highest yield with samples of low titer. The kit allows high throughput isolation of viral nucleic acid using Tecan or Beckman instruments.
	The purified viral RNA/DNA is devoid of proteins and nucleases, and is suitable for use in downstream applications that allow viral detection and genotyping.
System Overview	The kit allows efficient lysis of viral particles at elevated temperatures using Proteinase K and selective binding of viral nucleic acids to the silica matrix under highly denaturing conditions.
	The viral particles in the cell-free samples are lysed using Proteinase K and Lysis Buffer (L22) containing Carrier RNA at 56°C. The Lysis Buffer (L22) is specifically formulated to allow lysis of different types of viral particles.
	Ethanol is added to the lysate to a final concentration of 37% and the sample is loaded onto a PureLink [™] 96 Well Viral RNA/DNA Filter Plate. The viral RNA/DNA molecules bind to the silica-based media and impurities such as proteins and nucleases are removed by thorough washing with Wash Buffer. The RNA/DNA is then eluted in sterile, RNase free water.

About the Kit, Continued

Advantages	 The PureLink[™] Pro 96 Viral RNA/DNA Kit provides the following advantages: Highest sensitivity for samples with very low viral titer High-throughput processing of samples Automation using Tecan or Beckman robotic systems Purified nucleic acid free of contaminants such as proteins and nucleases
	Reliable performance of the purified viral nucleic acids in downstream applications
Carrier RNA	The Carrier RNA included with the PureLink [™] Pro 96 Viral RNA/DNA Kit is yeast tRNA (page viii). The presence of an excess amount of Carrier RNA as compared to viral nucleic acids during lysate preparation and purification:
	 Increases the binding of viral nucleic acids to the silica matrix
	 Reduces any viral nucleic acid degradation from nucleases present in the sample
	The purification protocol recommends using 5.6 µg Carrier RNA for 200 µl of sample. Most of the Carrier RNA is removed during the purification process because it is < 200 bp and any remaining Carrier RNA does not interfere with downstream applications. However,

depending on your application, you may validate the assay using less Carrier RNA.

About the Kit, Continued

Proteinase K	The Proteinase K is used for efficient lysis of viral particles. Proteinase K is active in the highly denaturing conditions of the lysis step.			
Downstream Applications	 The purified viral RNA and DNA is suitable for use in RT-PCR, qRT-PCR, and qPCR, and can be used for: Viral load monitoring Viral detection Viral genotyping 			
Kit Specifications	Dimensions:	Standard SBS (Society for Biomolecular Screening) footprint		
	Starting Material:	Up to 200 µl cell-free sample		
	Binding Capacity:	At least 40 μg nucleic acid		
	Filter Reservoir Capacity:	1.0 ml		
	Deep-Well Block Capacity:	1.09 ml		
	Centrifuge Compatibility:	Capable of centrifuging at ≥2,100 × g		
		Bucket depth 7 cm		
	Elution Volume:	100–150 μl		

Methods

Before Starting

Review the information in this section before starting. Guidelines are included for the recommended amount of starting material for use and to obtain high-quality RNA.

CAUTION

The PureLink[™] Viral RNA/DNA Kit buffers contain guanidine isothiocyanate. Always wear a laboratory coat, disposable gloves, and eye protection when handling buffers.

Do not add bleach or acidic solutions directly to solutions containing guanidine isothiocyanate or sample preparation waste as it forms reactive compounds and toxic gases when mixed with bleach or acids.



If there is any precipitate present in the buffers, warm the buffer up to 25–37°C to dissolve the precipitate before use.

General Handling of RNA Observe the following guidelines to prevent RNase contamination:

- Use disposable, individually wrapped, sterile plastic ware
- Use only sterile, new pipette tips (aerosol-barrier pipet tips recommended) and microcentrifuge tubes
- Wear latex gloves while handling reagents and RNA samples to prevent RNase contamination
- Always use proper microbiological aseptic techniques when working with RNA
- Use RNase AWAY[®] Reagent (page viii) to remove RNase contamination from surfaces (including from non-disposable equipment like homogenizers)

Before Starting, Continued

Starting Material	 The PureLink[™] Pro 96 Viral RNA/DNA Kit is designed to isolate viral nucleic acid from cell-free biological fluids such as plasma, serum, and CSF (cerebrospinal fluid) as well as cell culture supernatant using fresh or frozen samples. To obtain high yield of viral nucleic acids and minimize any degradation, follow these guidelines: Collect the sample (e.g. plasma or serum) and proceed immediately to purification. If desired, store the sample at 4°C for short-term storage (up to 4 hours) or freeze the sample at -20°C or -80°C for long-term storage. Do not freeze-thaw plasma or serum samples more than once. Remove any visible cryoprecipitates from samples by centrifugation at ~7,000 × g for 2–3 minutes. Use the clear supernatant immediately for purification. If you need to concentrate the cell culture supernatant use appropriate centrifugal concentrators.
Sample Volume	The PureLink [™] Pro 96 Viral RNA/DNA Kit can process sample volumes of up to 200 µl without preparation of any additional buffers or carry over of any impurities.
Preparing Wash Buffer	Add 350 ml of 96–100% ethanol to 87.5 ml PureLink [™] Pro 96 Wash Buffer II included in the kit. Place a check in the box on the Wash Buffer II label to indicate the ethanol is added. Store the Wash Buffer II with ethanol at room temperature.

Before Starting, Continued

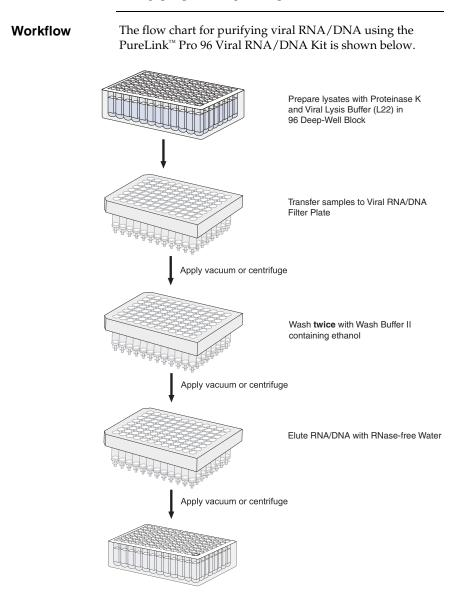
Preparing Carrier RNA	RN lesa am	The recommended purification protocol uses 5.6 µg Carrier RNA per sample (for up to 200 µl sample). If you wish to use less Carrier RNA per sample, you need to validate the amount of Carrier RNA needed for each sample type and downstream application.		
	То	prepare Carrier RNA (5.6 µg/sample):		
	1.	Add 310 μ l RNase-free Water (included with the kit) to 310 μ g lyophilized Carrier RNA (supplied with the kit) to obtain a 1 μ g/ μ l Carrier RNA stock solution.		
	2.	Mix thoroughly and divide the solution into smaller aliquots. Store the aliquots at –20°C. Avoid repeated freezing and thawing.		
	3.	Calculate the volume of Lysis Buffer/Carrier RNA mix required to process the desired number of samples simultaneously using the formula below:		
		$N \times 0.21$ ml (volume of Lysis Buffer/reaction) = A ml		
		A ml × 28 μ l/ml = B μ l		
		where		
		N = number of samples		
		A = calculated volume of Lysis Buffer (L22)		
		B = calculated volume of $1 \mu g/\mu l$ Carrier RNA stock solution to add to Lysis Buffer (L22)		
	4.	Thaw the required amount of $1 \mu g/\mu l$ Carrier RNA stock solution.		
	5.	In a sterile tube, add the volume of Carrier RNA stock solution (B, calculated as above) to the volume of Lysis Buffer (A, calculated as above). Mix gently by pipetting up and down. Avoid vortexing as it generates foam.		
	6.	Store at 4°C until use. Use the buffer within 1 hour.		
	Exa	ample:		
	and	e example below shows the amount of Lysis Buffer (L22) d Carrier RNA stock solution required to process samples using the above formula:		
	10	\times 0.21 ml (volume of Lysis Buffer/reaction) = 2.1 ml		
	2.1	$ml \times 28 \ \mu l/ml = 58.8 \ \mu l$		
	pro	prepare Lysis Buffer containing Carrier RNA for ocessing 10 samples, mix 58.8 µl Carrier RNA stock ution with 2.1 ml Lysis Buffer (L22).		

Before Starting, Continued

Instrument Compatibility	The PureLink [™] 96 Well Viral RNA/DNA Filter Plates are compatible with the following instruments:
for 96-Well Plates	• Vacuum Manifold: The manifold must accommodate the PureLink [™] 96 Well Plate (half-skirted filter plate) and be capable of collecting the filtrate (e.g. EveryPrep [™] Universal Vacuum Manifold from Invitrogen).
	• Centrifuge: The centrifuge must be capable of centrifuging 96-well plates at ≥2,100 × g, and accommodate a 7.0 cm microtiter plate stack.
	• Automated Liquid Handling Workstation: The workstation must be equipped with a vacuum manifold and a vacuum source, and accommodate the PureLink [™] 96 Well Plate (half-skirted filter plate).
Calibrating Vacuum for	Use a vacuum pressure of -12 to -15 in. Hg to obtain the best results.
Use with 96- Well Plates	Using higher vacuum pressure than the recommended pressure may cause sample splattering or inefficient DNA binding, while using lower vacuum pressure will affect the elution resulting in lower recovery.
	To check the vacuum pressure:
	1. Place an unused PureLink [™] 96 Well Viral RNA/DNA Filter Plate on top of the vacuum manifold.
	2. Apply vacuum and check the vacuum pressure on the vacuum regulator (usually attached to the manifold or a vacuum pump).
	3. Adjust the vacuum pressure on the regulator to obtain the recommended pressure of -12 to -15 in. Hg.
	Note: During purification the vacuum pressure may exceed the recommended value.

Experimental Overview

The viral nucleic acid purification procedure described below uses plate-based centrifugation or vacuum for highthroughput processing of samples.



General Guidelines



Follow the recommendations below to obtain the best results:

- Perform all steps at room temperature
- Be sure to add ethanol to Wash Buffer II before use (page 5)
- Perform the recommended wash steps to obtain highquality RNA
- Pipet water in the center of the silica matrix and perform a 1 minute incubation before elution

Handle all viruses in compliance with established institutional guidelines. Since safety requirements for mportant use and handling of viruses may vary at individual institutions, we recommend consulting the health and safety guidelines and/or officers at your institution. Be sure to take the appropriate precautions (wear a ٠ laboratory coat, disposable gloves, and eye protection) when handling viral samples. The eluates collected during wash steps contain biohazardous waste. Dispose the eluate and collection tubes appropriately as biohazardous waste. Elution The PureLink[™] Pro 96 Viral RNA/DNA Kit elutes viral Volume nucleic acid in $100-150 \mu$ l for downstream applications. Note: Using larger elution volume decreases the viral nucleic acid concentration in the eluate. Automation For automation scripts, contact Technical Support page 17. Scripts *Continued on next page*

Preparing Lysates

Materials	• Cell-free samples (such as plasma or serum samples))
Needed	 Appropriate amount of Lysis Buffer (L22) containing Carrier RNA (see page 6) 	5
	 Heat block for 96-well plate, water bath, or incubator to 56°C 	r set
	• 96–100% ethanol	
	• 96-well Deep-Well Blocks (supplied with kit)	
	• Sterile, RNase-free PBS or 0.9% NaCl	
	• Centrifuge capable of ≥2,100 × g with buckets that ca accommodate 96-well plates that have a plate height clearance of 7.0 cm	
Preparing Lysate	The lysate preparation protocol is described below for 20 of starting material.	
	 Add 25 µl Proteinase K into each well of a Deep-Wel Block that contain samples. 	1
	2. Add 200 µl of cell-free sample (equilibrated to room	
	temperature) to each well. Note: If sample is less than 200 μl, adjust the final volume of sample to 200 μl using sterile PBS or 0.9% NaCl.	of the
	 Add 200 µl Lysis Buffer (containing 5.6 µg Carrier Rl Mix by pipetting up and down 3–5 times. 	NA).
	 Seal the plate with the foil tape and incubate at 56°C 15 minutes. 	for
	5. Centrifuge the plate briefly to collect the liquid in the bottom of the well.	9
	 Add 250 μl 96–100% ethanol to each well to obtain a ethanol concentration of 37% and mix by pipetting u and down 3–5 times. 	
	Note: If you are processing multiple samples, you may add ethanol to all wells and then mix.	l
	7. Incubate the lysate with ethanol for 5 minutes at root temperature.	m
	 Proceed to Purification Using Centrifugation (page Purification Using a Vacuum Manifold (page 12), o Purification Using the EveryPrep[™] Universal Vacuu Manifold (page 17). 	r

Purification Using Centrifugation

Materials Needed	• • • •	Viral lysate from cell-free samples (see previous page) Centrifuge capable of ≥2,100 × g with buckets that can accommodate 96-well plates that have a plate height clearance of 7.0 cm PureLink [™] Pro 96 Wash Buffer II with ethanol (page 5) PureLink [™] Pro 96 RNase-free Water (supplied with kit) PureLink [™] 96 Well Viral RNA/DNA Filter Plate (supplied with kit) 96-well Deep-Well Blocks (supplied with kit) Foil tape (supplied with kit)
Purification	Perform all steps at room temperature:	
Procedure Using Centrifugation	1.	Transfer each lysate sample to a well of the PureLink [™] 96 Well Viral RNA/DNA Filter Plate. Cover any unused wells with foil tape.
	2.	Place the Viral Filter Plate onto a new or used 96-well Deep-Well Block and centrifuge at ≥2,100 × g for 1–2 minutes.
	3.	Discard the flow-through and place the Viral Filter Plate back on the Deep-Well Block.
	4.	Add 500 μl Wash Buffer II (with ethanol) into each well of the Viral Filter Plate.
	5.	Centrifuge at ≥2,100 × g for 1–2 minutes, until all liquid passes through the Viral Filter Plate.
	6.	Discard the flow-through and place the Viral Filter Plate back on the Deep-Well Block.
	7.	Repeat Steps 4–6 above.
	8.	Centrifuge the Viral Filter Plate at $\ge 2,100 \times g$ for 10 minutes to dry the plate completely.
	9.	Add 100 μ l RNase-free water to the center of each well and incubate the plate for 1 minute at room temperature.
	10.	Place the Viral Filter Plate onto a new Deep-Well Block (supplied in the kit). Centrifuge at $\geq 2,100 \times g$ for 1–2 minutes.
	11.	Use the RNA/DNA for the desired downstream application. To store the purified RNA/DNA, cover the wells with foil tape, and store at -80°C.

Purification Using a Vacuum Manifold

Materials Needed	• • • •	Viral lysate from cell-free samples (see previous page) Vacuum manifold and a vacuum pump (producing pressure of -12 to -15 in. Hg) or automated liquid handling workstation PureLink [™] Pro 96 Wash Buffer II with ethanol (page 5) PureLink [™] Pro 96 RNase-free Water (supplied with kit) PureLink [™] 96 Well Viral RNA/DNA Filter Plate (supplied with kit) 96-well Deep-Well Blocks (supplied with kit) Foil tape (supplied with kit)
Binding and Washing RNA/DNA	ma	semble the vacuum manifold according to the nufacturer's instructions. For a protocol using the eryPrep [™] Universal Vacuum Manifold, see page 17. Place a PureLink [™] 96 Well Viral RNA/DNA Filter Plate on top of the vacuum manifold. Transfer the lysates from the Deep-Well Block to the Viral Filter Plate. Cover unused wells with foil tape. Apply vacuum for 2 minutes. Release vacuum. Add 500 µl Wash Buffer II (with ethanol) into each well of the Viral Filter Plate. Apply vacuum for 2 minutes. Release vacuum. Repeat Steps 4 and 5 above. Place the Viral Filter Plate on the vacuum manifold and apply vacuum for 10 minutes. Release vacuum. Alternatively, centrifuge the Viral Filter Plate at ≥2,100 × g for 10 minutes to dry the plate completely. Place the Viral Filter Plate with the filter side down on a stack of paper towels, and pat firmly to blot any residual liquid. Proceed to elution using vacuum manifold (Steps 10–12, next page), or using centrifugation (Steps 9–10, previous page).

Purification Using a Vacuum Manifold,

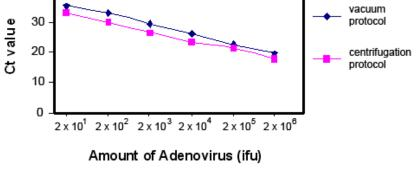
Continued

Eluting RNA/DNA	10.	Place a clean Deep-Well Block (included in the kit) in the vacuum manifold in place of the waste collection tray and place the Viral Filter Plate on top of the Deep- Well Block.
	11.	Add 150 μ l of RNase-free water (included in the kit) to the center of the membrane in each well of the Viral Filter Plate and incubate for 1 minute at room temperature.
	12.	Apply vacuum for 2 minutes. Release vacuum. The RNA/DNA is eluted into the Deep-Well Block in a volume of $110-130 \mu$ l.
	13.	Use the RNA/DNA for the desired downstream

volume of 110–130 μl.
13. Use the RNA/DNA for the desired downstream application. To store the purified RNA/DNA, cover the wells with foil tape, and store at -80°C.

Example of Expected Results

Analyzing Viral RNA/DNA	Since the amount of viral RNA/DNA present in cell-free body fluids is low we recommend that you do not use UV absorbance to determine yield. Use qRT-PCR or RT-PCR for RNA virus, and qPCR and PCR for DNA virus using appropriate viral-specific probes to determine yield or presence of viral nucleic acid.		
	To analyze viral nucleic acid size, use agarose gel electrophoresis followed by hybridization using viral specific labeled probes and autoradiography.		
Results	Examples of results obtained after purification of viral RNA/DNA using the PureLink [™] Pro 96 Viral RNA/DNA Kit are shown below.		
	Plasma samples (200 µl) were spiked with Adenovirus at the indicated ifu (infectious unit). Viral DNA were purified using the PureLink [™] 96 Viral RNA/DNA Kit as described in this manual.		
	qPCR was performed on the resulting samples using the Platinum [®] quantitative PCR Supermix-UDG Kit with TaqMan [®] primers in an ABI 7900HT instrument.		
	Consistent recovery of viral DNA is reflected by linear performance of the PureLink [™] Pro 96 Viral RNA/DNA Kit when analyzed by qPCR.		
40 -			
ou 30 -			



Troubleshooting

Problem	Cause	Solution
Low nucleic acid yield	Incomplete lysis or the column or	Be sure to use the appropriate volumes of reagents during lysate preparation.
	matrix is clogged	Do not use more than 200 μ l sample.
		If cryoprecipitates are visible in frozen viral samples, remove the cryoprecipitates by centrifugation (page 5) to avoid clogging of the column or matrix.
		Precipitates in buffers may affect lysis. Dissolve precipitates by warming the buffers at 25–37°C.
	Poor quality of sample material	Avoid repeated freezing and thawing of samples. Use fresh samples and process immediately after collection or use samples thawed only once for best results.
		Check the quality of the RNA in the original samples using qRT-PCR or RT-PCR.
	Lysis Buffer without Carrier RNA used or	To prepare lysates from cell-free samples, use Lysis Buffer (L22) with Carrier RNA (page 6).
	Carrier RNA inactivated	Once the Carrier RNA is reconstituted in water, aliquot the Carrier RNA and store at -20° C. Do not perform multiple freeze-thaw cycles.
	Incorrect binding conditions	For efficient binding of viral nucleic acids, always add ethanol to the lysate to a final concentration of 37% prior to loading the lysate onto the column or matrix.
	Ethanol not added to Wash Buffer II	Be sure to add 96–100% ethanol to Wash Buffer II as described on page 5. Do not use denatured 95% ethanol.
	Incorrect elution conditions	Add water to the center of the silica matrix and perform incubation for 1 minute with water before eluting.

Troubleshooting, Continued

Problem	Cause	Solution
Low nucleic acid yield, continued	Quantitation performed using UV absorbance	Since viral nucleic acids are present in low amounts in cell-free samples, do not use UV absorbance for quantitation. Analyze viral nucleic acids using qRT- PCR, RT-PCR, qPCR, or PCR.
RNA degraded	RNA contaminated with RNase	Follow the guidelines on page 4 to prevent RNase contamination.
	Poor quality of samples	Always use fresh samples or samples frozen at –80°C. For lysis, process the sample quickly to avoid degradation.
Poor performance of nucleic acids in downstream enzymatic reactions	Presence of ethanol or use of denatured 95% ethanol in	Traces of ethanol from the Wash Buffer II can inhibit downstream enzymatic reactions.
	purified nucleic acids	To remove residual buffer, discard Wash Buffer II flow through. Completely dry the Viral Filter Plate, and always use a new Deep-Well Block for elution.
		Use only 96–100% ethanol. Do not use denatured 95% ethanol.
	Assay may be sensitive to Carrier RNA concentration	You may need to optimize the amount of Carrier RNA that is required for optimal purification and is suitable for your downstream applications.
	Reagents for enzymatic reactions inactive	Ensure that the enzymes and reagents used for performing downstream applications have not expired or inactivated. Repeat the reaction with fresh enzyme and reagents.
	Viral nucleic acid eluate too dilute	Optimize the amount of viral nucleic acid eluate required for your specific application and perform elution using the desired kit and elution volume (10–150 µl).
Carrier RNA not enough to process samples	Incorrect Carrier RNA amount used per sample	We recommend using a maximum of 5.6 µg Carrier RNA per sample when processing up to 200 µl sample volume.

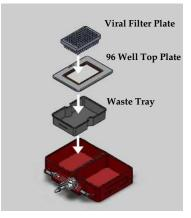
Appendix

Purification Using the EveryPrep[™] Universal Vacuum Manifold

Introduction	Instructions are provided below to purify DNA using the EveryPrep [™] Universal Vacuum Manifold (see page viii). Refer to the manual for the EveryPrep [™] Universal Vacuum Manifold for detailed instructions on operation with the 96 Well Top Plate. All steps are performed at room temperature.	
Materials Needed	 Viral lysate from cell-free samples (see previous page) Vacuum manifold and vacuum pump (producing pressure of 12–15 in. Hg) PureLink[™] Pro 96 Wash Buffer II with ethanol (page 5) PureLink[™] Pro 96 RNase-free Water (supplied with kit) PureLink[™] 96 Well Viral RNA/DNA Filter Plate (supplied with kit) 96-well Deep-Well Blocks (supplied with kit) Foil tape (supplied with kit) 	

Purification Using the EveryPrep[™] Universal Vacuum Manifold, Continued

- EveryPrep[™] Universal Vacuum Manifold Assembly
- Assemble the EveryPrep[™] Universal Vacuum Manifold: Place the Waste Tray in the Binding Chamber, cover the top with the 96 Well Top Plate, and place the PureLink[™] Viral RNA/DNA Filter Plate over the Top Plate.



2. Proceed to Binding RNA/DNA, below.

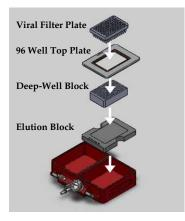
Binding and Washing RNA/DNA

- 1. Transfer the lysates from the Deep-Well Block to the Viral Filter Plate. Cover unused wells with foil tape.
- 2. Apply vacuum for 2 minutes. Release vacuum.
- 3. Add 500 μl Wash Buffer II (with ethanol) into each well of the Viral Filter Plate.
- 4. Apply vacuum for 2 minutes. Release vacuum.
- 5. Repeat Steps 4 and 5 above.
- 6. Apply vacuum for an additional 10 minutes. Release vacuum.
- Place the Viral Filter Plate with the filter side down on a stack of paper towels, and pat firmly to blot any residual liquid.
- 8. Proceed to Eluting RNA/DNA, next page.

Purification Using the EveryPrep[™] Universal Vacuum Manifold, Continued

Eluting RNA/DNA

 Prepare the EveryPrep[™] Universal Vacuum Manifold for elution: Place the Elution Block and a clean Deep-Well Block in the Elution Chamber, cover the top with the 96 Well Top Plate, and place the PureLink[™] Viral RNA/DNA Filter Plate over the Top Plate.



- 2. Add 150 µl of RNase-free water (included in the kit) to the center of the membrane in each well of the Viral Filter Plate and incubate for 1 minute at room temperature.
- 3. Apply vacuum for 2 minutes. Release vacuum. The RNA/DNA is eluted into the Deep-Well Block in a volume of 110–130 μ l.
- 4. Use the RNA/DNA for the desired downstream application. To store the purified RNA/DNA, cover the wells with foil tape, and store at -80°C.

Technical Support



Visit the Invitrogen website at <u>www.invitrogen.com</u> for:

- Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.
- Complete technical support contact information
- Access to the Invitrogen Online Catalog

Additional product information and special offers

Contact Us

For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our website (<u>www.invitrogen.com</u>).

Corporate Headquarters:

5791 Van Allen Way Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 602 6500 E-mail: <u>tech_support@invitrogen.com</u>

Japanese Headquarters: LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519 E-mail: jpinfo@invitrogen.com

European Headquarters:

Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: +44 (0) 141 814 6100 Tech Fax: +44 (0) 141 814 6117 E-mail: <u>eurotech@invitrogen.com</u>

Certificate of Analysis The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to <u>www.invitrogen.com/support</u> and search for the Certificate of Analysis by product lot number, which is printed on the box.

Purchaser Notification

Limited Use Label License No. 5: Invitrogen Technology The purchase of this product conveys to the buyer the nontransferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008 or outlicensing@lifetech.com

Purchaser Notification, Continued

Disclaimer	Any scripts or methods for robotic systems have been developed with minimal involvement from the manufacturer and are provided only as a convenience to customers to use at their own risk. Invitrogen assumes no responsibility or liability for any special, incidental, direct, or consequential loss or damages whatsoever. Invitrogen provides no warranties of any kinds, express or implied, including without limitation implied warranties or merchantability or fitness for a particular purpose.
Limited Warranty	Invitrogen (a part of Life Technologies Corporation) is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Support Representatives. All Invitrogen products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. This warranty limits the Company's liability to only the price of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. The Company reserves the right to select the method(s) used to analyze a product unless the Company gargees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Support Representatives. Life Technologies Corporation shall have no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.
	RNase $AWAY^{\otimes}$ is a registered trademark of Molecular Bio-Products, Inc. TaqMan [®] is a registered trademark of Roche Molecular Systems, Inc. BioRobot [®] is a registered trademark of QIAGEN Corp. Freedom EVO [™] is a trademark of Tecan. Biomek [®] is a registered trademark of Beckman Coulter, Inc.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

invitrogen

Corporate Headquarters Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech_support@invitrogen.com

For country-specific contact information, visit our web site at www.invitrogen.com