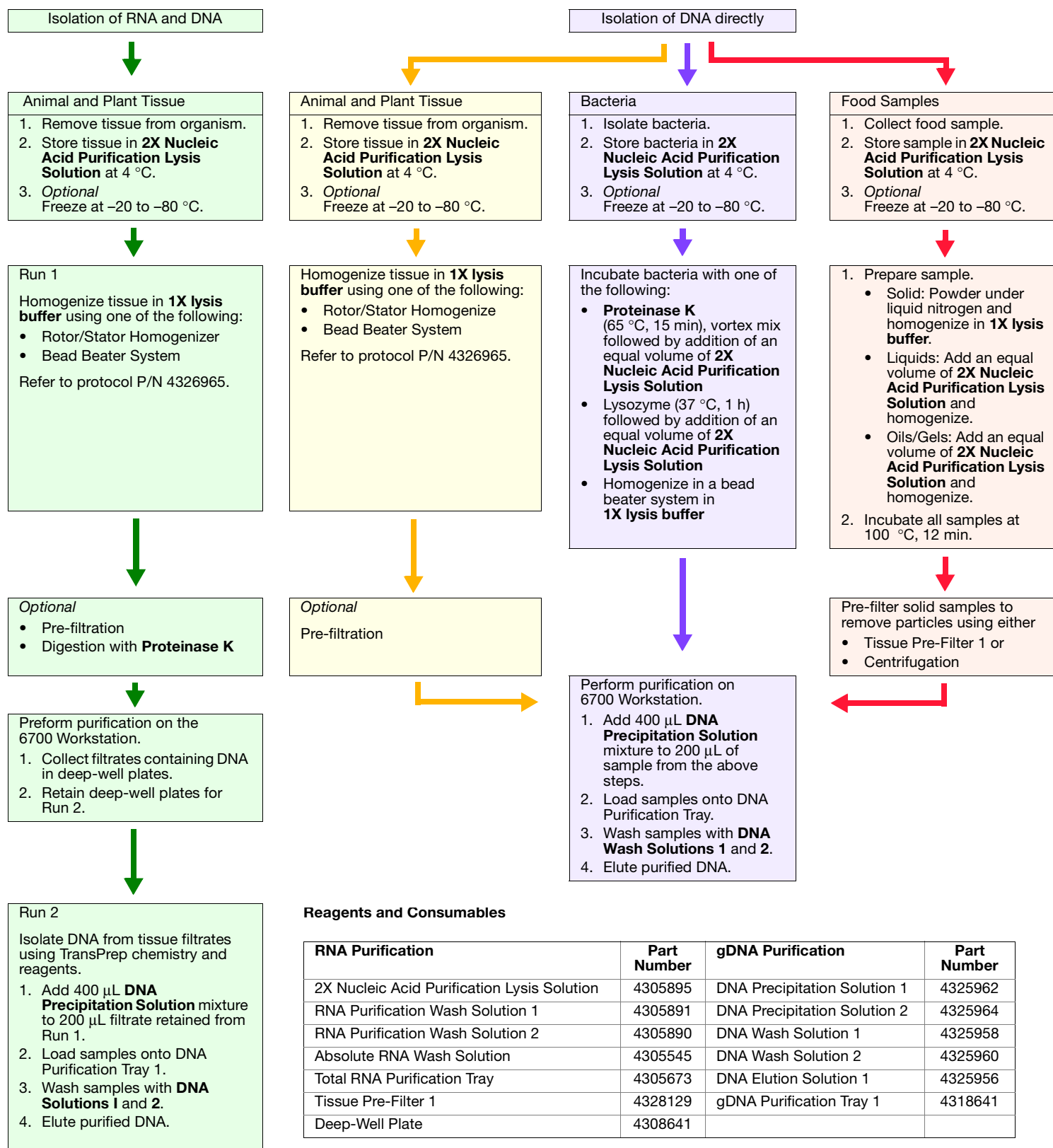


⚠ WARNING Biological samples have the potential to transmit infectious disease. For safety and biohazard guidelines, please refer to the “Safety” section in the *TransPrep Chemistry Purification of gDNA from Filtrates Obtained After the Isolation of RNA from Homogenized Animal or Plant Tissue Samples Protocol*, P/N 4326965. Follow specific safety practices when using this instrument. For all chemicals in **bold** type below, please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.



⚠ WARNING Biological samples have the potential to transmit infectious disease. For safety and biohazard guidelines, please refer to the "Safety" section in the *TransPrep Chemistry Purification of gDNA from Filtrates Obtained After the Isolation of RNA from Homogenized Animal or Plant Tissue Samples Protocol*, P/N 4326965. Follow specific safety practices when using this instrument. For all chemicals in **bold** type below, please read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

STEP	ACTION																																																																																									
1	Remove and store tissue	<div>a. Remove tissue from organism.</div> <div>b. Store tissue in 2X Nucleic Acid Purification Lysis Solution or freeze at −20 to −80 °C.</div> <div>For other types of samples see the <i>TransPrep Chemistry Purification of gDNA from Filtrates Obtained After the Isolation of RNA from Homogenized Animal or Plant Tissue Samples Protocol</i>.</div>																																																																																								
2	Prepare to perform Run 1	<div>a. Create the following new protocols:<div><div>• RNA Archive protocol, if performing Run 1</div><div>• DNA Precipitation protocol</div><div>• DNA Archive protocol</div></div></div> <div>b. Lyse and homogenize the tissue.</div> <div>c. Perform pre-filtration, if the tissue sample needs it.</div> <div>d. Pipette 250 μL of each lysed and homogenized tissue sample into wells of a 300-μL flat-bottom cell-culture plate.</div>																																																																																								
		<div><div><div><div>New RNA/DNA Archive Protocol</div><div><div>Protocol Name: <div>DNA Archive Protocol</div><div><input checked="" type="checkbox"/> In Use</div></div><div>Conditions for Transferring Samples to the Purification Tray</div><div><div>Lysis/DNA Precipitation Input</div><div><div>Deep-Well Plate</div><div>First Transfer: <div>None</div><div>550</div><div>3</div><div>600</div></div><div>Second Transfer: <div>0</div><div>0</div><div>0</div></div><div><input type="checkbox"/> High Viscosity Sample</div></div><div>Filtration Conditions</div><div><div><input type="checkbox"/> Create Deep-Well Filtrate Plate</div><div>Incubation Time: <div>0</div> (min.)</div><div>Vacuum Time: <div>120</div> (sec.)</div><div>Vacuum Pressure: <div>20</div> %</div></div><div>Wash Conditions</div><table><thead><tr><th>Step</th><th>Add</th><th>Volume (μL)</th><th>Temp. (°C)</th><th>Incubation (min)</th><th>Vacuum (sec)</th><th>Repeat (count)</th><th>Vacuum (%)</th></tr></thead><tbody><tr><td>1. <input checked="" type="checkbox"/></td><td>Wash Solution 1</td><td>650</td><td></td><td>0</td><td>90</td><td>1</td><td>20</td></tr><tr><td>2. <input checked="" type="checkbox"/></td><td>Wash Solution 2</td><td>650</td><td></td><td>0</td><td>90</td><td>1</td><td>20</td></tr><tr><td>3. <input type="checkbox"/></td><td></td><td>300</td><td></td><td>0</td><td>120</td><td>2</td><td>20</td></tr><tr><td>4. <input type="checkbox"/></td><td></td><td>300</td><td></td><td>0</td><td>120</td><td>1</td><td>20</td></tr><tr><td>5. <input type="checkbox"/></td><td></td><td>300</td><td></td><td>0</td><td>120</td><td>1</td><td>20</td></tr><tr><td>6. <input type="checkbox"/></td><td></td><td>300</td><td></td><td>0</td><td>120</td><td>1</td><td>20</td></tr><tr><td>7. <input type="checkbox"/></td><td></td><td>300</td><td></td><td>0</td><td>120</td><td>1</td><td>20</td></tr><tr><td colspan="4">Pre-Elution Vacuum</td><td></td><td>30</td><td></td><td>30</td></tr><tr><td colspan="2">Elution Solution</td><td>150</td><td>-</td><td>2</td><td>120</td><td>1</td><td>20</td></tr><tr><td colspan="2"><input type="checkbox"/> Final Addition Fluid</td><td></td><td></td><td></td><td></td><td></td><td></td></tr></tbody></table><div><div>Cancel</div><div>OK</div></div></div></div></div><div>New RNA/DNA Archive Protocol</div></div></div>	Step	Add	Volume (μL)	Temp. (°C)	Incubation (min)	Vacuum (sec)	Repeat (count)	Vacuum (%)	1. <input checked="" type="checkbox"/>	Wash Solution 1	650		0	90	1	20	2. <input checked="" type="checkbox"/>	Wash Solution 2	650		0	90	1	20	3. <input type="checkbox"/>		300		0	120	2	20	4. <input type="checkbox"/>		300		0	120	1	20	5. <input type="checkbox"/>		300		0	120	1	20	6. <input type="checkbox"/>		300		0	120	1	20	7. <input type="checkbox"/>		300		0	120	1	20	Pre-Elution Vacuum					30		30	Elution Solution		150	-	2	120	1	20	<input type="checkbox"/> Final Addition Fluid							
Step	Add	Volume (μL)	Temp. (°C)	Incubation (min)	Vacuum (sec)	Repeat (count)	Vacuum (%)																																																																																			
1. <input checked="" type="checkbox"/>	Wash Solution 1	650		0	90	1	20																																																																																			
2. <input checked="" type="checkbox"/>	Wash Solution 2	650		0	90	1	20																																																																																			
3. <input type="checkbox"/>		300		0	120	2	20																																																																																			
4. <input type="checkbox"/>		300		0	120	1	20																																																																																			
5. <input type="checkbox"/>		300		0	120	1	20																																																																																			
6. <input type="checkbox"/>		300		0	120	1	20																																																																																			
7. <input type="checkbox"/>		300		0	120	1	20																																																																																			
Pre-Elution Vacuum					30		30																																																																																			
Elution Solution		150	-	2	120	1	20																																																																																			
<input type="checkbox"/> Final Addition Fluid																																																																																										
3	Perform Run 1	<div>Perform Run 1 if RNA is required. Go to step 4 if only DNA is required.</div> <div>a. Make selections on the protocol tab.</div> <div>b. Set up the deckspace.</div> <div>c. Start the run.</div> <div>d. Finish the run.</div>																																																																																								
4	Prepare to perform Run 2	<div>a. Prepare DNA Precipitation solution mixture. Pipette the components listed below into a 120-mL reagent reservoir.</div> <div>IMPORTANT! These reagents should be mixed immediately before use and should not be stored as a mixture.</div> <table><thead><tr><th>Component</th><th>Volume for One Well (μL)*</th><th>Volume for 96 Wells (mL)*</th></tr></thead><tbody><tr><td>DNA Precipitation Solution 1</td><td>100</td><td>9.6</td></tr><tr><td>DNA Precipitation Solution 2</td><td>300</td><td>28.8</td></tr><tr><td>Total Volume</td><td>400</td><td>38.4</td></tr></tbody></table> <div>* Volumes given are for 200 μL of filtrate per well collected in a deep-well plate during the first run on the 6700 Workstation.</div> <div>b. Load the reagent reservoir on the 6700 Workstation.</div> <div>c. Remove the deep-well plate from Run 1 and place the plate in the secondary input station.</div>	Component	Volume for One Well (μL)*	Volume for 96 Wells (mL)*	DNA Precipitation Solution 1	100	9.6	DNA Precipitation Solution 2	300	28.8	Total Volume	400	38.4																																																																												
Component	Volume for One Well (μL)*	Volume for 96 Wells (mL)*																																																																																								
DNA Precipitation Solution 1	100	9.6																																																																																								
DNA Precipitation Solution 2	300	28.8																																																																																								
Total Volume	400	38.4																																																																																								
5	Perform Run 2	<div>a. Run both the new DNA Precipitation protocol and the new DNA Archive protocol.</div> <div>b. Make selections on the protocol tab.</div> <div>c. Set up the deckspace.</div> <div>d. Start the run.</div> <div>e. Finish the run.</div>																																																																																								