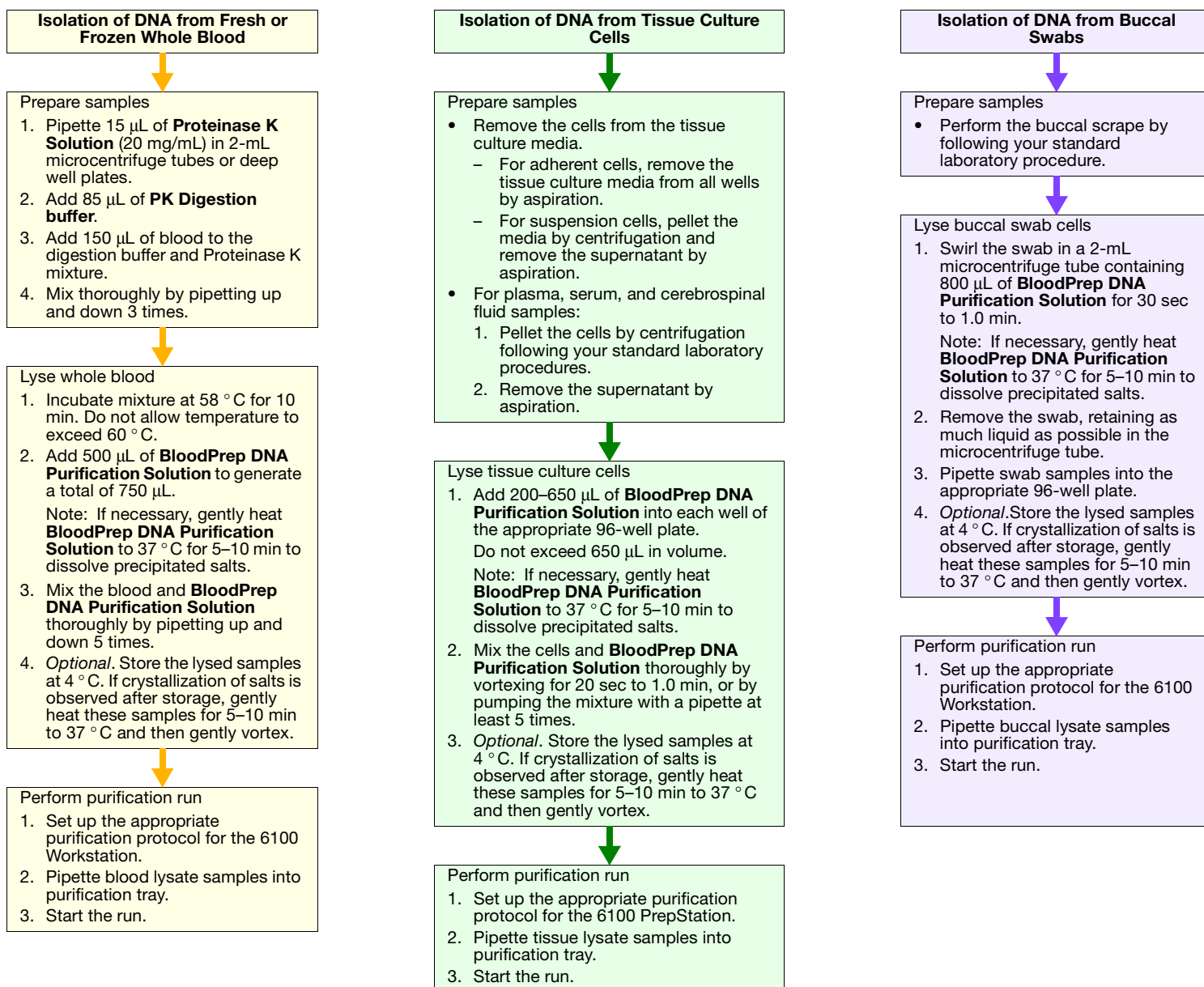


DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs on the ABI PRISM™ 6100 Nucleic Acid PrepStation

⚠ WARNING Biological samples have the potential to transmit infectious disease. For safety and biohazard guidelines, please refer to the “Safety” section in the *DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs Protocol*, P/N 4343586. 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.



Reagents and Consumables

Item	Part Number	Item	Part Number	Item	Part Number
96-Well Deep-well plate	4308641	BloodPrep™ DNA Purification Solution	4342775	Proteinase K Solution (20 mg/mL)	4333793
96-Well Optical Reaction Plate with Barcode	4306737	BloodPrep™ DNA Wash Solution	4342949	Splash Guard	4311758
Archive Tray Covers	4306286	BloodPrep™ PK Digestion Buffer	4342777	300-µL, flat-bottom, 96-well cell culture plate	MLS ^a
BloodPrep™ DNA Elution Solution 1	4342951	Genomic DNA Purification Tray II	4330172	Ethanol, 70%	MLS
BloodPrep™ DNA Elution Solution 2	4342950	Microcentrifuge tubes, 2 mL	4305936	Water, molecular biology grade	MLS

a. Major Laboratory Supplier

DNA Isolation from Fresh and Frozen Blood, Tissue Culture

Cells, and Buccal Swabs on the ABI PRISM™ 6100 Nucleic Acid PrepStation (continued)

⚠ WARNING Biological samples have the potential to transmit infectious disease. For safety and biohazard guidelines, please refer to the “Safety” section in the *DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs Protocol*, P/N 4343586. 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	Lysate Preparation from Fresh or Frozen Whole Blood
1	Prepare samples	<p>a. Pipette 15 µL of Proteinase K Solution (20 mg/mL) into 2-mL microcentrifuge tubes or deep well plates. Note: If blood is <24 hrs old, add the samples directly to tubes with Proteinase K Solution and proceed to step 2b.</p> <p>b. Add 85 µL of PK Digestion buffer.</p> <p>c. Add 150 µL of blood to the digestion buffer and Proteinase K mixture. Note: If using animal blood, use suggested volumes in Protocol, P/N 4343586.</p> <p>d. Mix thoroughly by pipetting up and down 3 times.</p>
2	Lyse whole blood	<p>a. Incubate mixture at 58 °C for 10 min. IMPORTANT! Temperatures in excess of 60 °C may cause degradation of isolated DNA and reduce activity of the Proteinase K.</p> <p>b. Add 500 µL of BloodPrep DNA Purification Solution to generate a total of 750 µL. IMPORTANT! The BloodPrep DNA Purification Solution may require gentle heating to 37 °C for 5–10 min to dissolve precipitated salts.</p> <p>c. Mix the blood and BloodPrep DNA Purification Solution thoroughly by pipetting up and down 5 times. IMPORTANT! The incubated blood sample mixture and purification solution must be thoroughly mixed with the pipette before proceeding, because vortexing does not effectively mix the viscous samples and the Proteinase K solution, no matter how long you vortex.</p> <p>d. <i>Optional</i>. If you cannot proceed with purification immediately, store the lysed samples at 4 °C. If crystallization of salts is observed after storage, gently heat these samples for 5–10 min to 37 °C and then gently vortex.</p>
STEP	ACTION	Lysate Preparation from Tissue Culture Cells
1	Prepare samples	<p>Remove the cells from the tissue culture media.</p> <ul style="list-style-type: none"> For adherent cells, remove the tissue culture media from all wells by aspiration. For suspension cells, pellet the media by centrifugation and remove the supernatant by aspiration. <p>Note: For plasma, serum, and cerebrospinal fluid samples, pellet the cells by centrifugation following your standard laboratory procedures, and remove the supernatant by aspiration.</p>
2	Lyse cells	<p>a. Add 200–650 µL of BloodPrep DNA Purification Solution into each well of the appropriate 96-well plate. Note: Do not exceed 650 µL in volume unless the sample is extremely viscous. If viscous, reduce the sample input volume or increase the volume of BloodPrep DNA Purification Solution appropriately. IMPORTANT! The BloodPrep DNA Purification Solution may require gentle heating to 37 °C for 5–10 min to dissolve precipitated salts.</p> <p>b. Mix the cells and BloodPrep DNA Purification Solution thoroughly by vortexing for 20 sec to 1.0 min, or by pumping the mixture with a pipette at least 5 times.</p> <p>c. <i>Optional</i>. If you cannot proceed with purification immediately, store the lysed samples at 4 °C. If crystallization of salts is observed after storage, gently heat these samples for 5–10 min to 37 °C and then gently vortex.</p>
STEP	ACTION	Lysate Preparation from Buccal Swab Cells
1	Prepare samples	Perform the buccal scrape by following your standard laboratory procedure.
2	Lyse cells	<p>a. Swirl the swab in a 2-mL microcentrifuge tube containing 800 µL of BloodPrep DNA Purification Solution for 30 sec to 1.0 min. IMPORTANT! The BloodPrep DNA Purification Solution may require gentle heating to 37 °C for 5–10 min to dissolve precipitated salts.</p> <p>b. Remove the swab, retaining as much liquid as possible in the microcentrifuge tube.</p> <p>c. Pipette swab samples into the purification tray.</p> <p>d. <i>Optional</i>. If you cannot proceed with purification immediately, store the lysed samples at 4 °C. If crystallization of salts is observed after storage, gently heat these samples for 5–10 min to 37 °C and then gently vortex.</p>

STEP	ACTION	Total DNA Purification of Blood, Tissue Culture Cells, or Buccal Cells Using the 6100 PrepStation																																																																																																																																																																																																																										
1	Set up the 6100 PrepStation for blood samples	Set up the DNA purification parameters for the blood sample on the protocol tab of the 6100 PrepStation software:																																																																																																																																																																																																																										
		<table><thead><tr><th>Step</th><th>Description</th><th>Volume (μL)</th><th>Position</th><th>Incubation (sec)</th><th>Vacuum (%)</th><th>Time (sec)</th></tr></thead><tbody><tr><td colspan="7">Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the 6100 waste station.</td></tr><tr><td>1</td><td>Load Samples^a</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>2</td><td>Add BloodPrep DNA Purification Solution^a</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>400</td></tr><tr><td>3</td><td>Add BloodPrep DNA Wash Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>4</td><td>Add BloodPrep DNA Wash Solution</td><td>600</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>5</td><td>Add BloodPrep DNA Wash Solution</td><td>300</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>6</td><td>Pre-Elution Vacuum</td><td>–</td><td>Waste</td><td>0</td><td>100</td><td>120</td></tr><tr><td>7</td><td>Touch Off</td><td>–</td><td>Waste</td><td>–</td><td>–</td><td>–</td></tr><tr><td>8</td><td>BloodPrep DNA Elution Solution 1^b</td><td>100</td><td>Collection</td><td>180</td><td>60</td><td>120</td></tr><tr><td>9</td><td>BloodPrep DNA Elution Solution 2^c</td><td>100</td><td>Collection</td><td>0</td><td>60</td><td>120</td></tr><tr><td>10</td><td>Touch Off</td><td>–</td><td>Collection</td><td>–</td><td>–</td><td>–</td></tr><tr><td colspan="7">a.If the well evacuates slowly, increase the vacuum setting to 100% and repeat the step.</td></tr><tr><td colspan="7">b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. You may lower the total elution volume to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.</td></tr><tr><td colspan="7">c.It is very important that you use the elution solutions in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.</td></tr><tr><td rowspan="10"></td><td rowspan="10">Set up the 6100 PrepStation for tissue culture or buccal swab samples</td><td>Set up the DNA purification parameters for the blood sample on the protocol tab of the 6100 PrepStation software:</td></tr><tr><td><table><thead><tr><th>Step</th><th>Description</th><th>Volume (μL)</th><th>Position</th><th>Incubation (sec)</th><th>Vacuum (%)</th><th>Time (sec)</th></tr></thead><tbody><tr><td colspan="7">Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..</td></tr><tr><td>1</td><td>Load Samples^a</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>2</td><td>Add BloodPrep DNA Purification Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>3</td><td>Add BloodPrep DNA Wash Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>4</td><td>Add BloodPrep DNA Wash Solution</td><td>600</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>5</td><td>Add BloodPrep DNA Wash Solution</td><td>300</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>9</td><td>Pre-Elution Vacuum</td><td>–</td><td>Waste</td><td>0</td><td>100</td><td>120</td></tr><tr><td>10</td><td>Touch Off</td><td>–</td><td>Waste</td><td>–</td><td>–</td><td>–</td></tr><tr><td>11</td><td>BloodPrep DNA Elution Solution 1^b</td><td>100</td><td>Collection</td><td>180</td><td>60</td><td>120</td></tr><tr><td>12</td><td>BloodPrep DNA Elution Solution 2^c</td><td>100</td><td>Collection</td><td>0</td><td>60</td><td>120</td></tr><tr><td>13</td><td>Touch Off</td><td>–</td><td>Collection</td><td>–</td><td>–</td><td>–</td></tr><tr><td colspan="7">a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.</td></tr><tr><td colspan="7">b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.</td></tr><tr><td colspan="7">c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.</td></tr><tr><td rowspan="2">2</td><td rowspan="2">Perform the purification run</td><td>a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.</td></tr><tr><td>b. Start the run.</td></tr></tbody></table></td></tr></tbody></table>	Step	Description	Volume (μL)	Position	Incubation (sec)	Vacuum (%)	Time (sec)	Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the 6100 waste station.							1	Load Samples ^a	650	Waste	0	80	300	2	Add BloodPrep DNA Purification Solution ^a	650	Waste	0	80	400	3	Add BloodPrep DNA Wash Solution	650	Waste	0	80	60	4	Add BloodPrep DNA Wash Solution	600	Waste	0	80	60	5	Add BloodPrep DNA Wash Solution	300	Waste	0	80	60	6	Pre-Elution Vacuum	–	Waste	0	100	120	7	Touch Off	–	Waste	–	–	–	8	BloodPrep DNA Elution Solution 1 ^b	100	Collection	180	60	120	9	BloodPrep DNA Elution Solution 2 ^c	100	Collection	0	60	120	10	Touch Off	–	Collection	–	–	–	a.If the well evacuates slowly, increase the vacuum setting to 100% and repeat the step.							b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. You may lower the total elution volume to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.							c.It is very important that you use the elution solutions in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.								Set up the 6100 PrepStation for tissue culture or buccal swab samples	Set up the DNA purification parameters for the blood sample on the protocol tab of the 6100 PrepStation software:	<table><thead><tr><th>Step</th><th>Description</th><th>Volume (μL)</th><th>Position</th><th>Incubation (sec)</th><th>Vacuum (%)</th><th>Time (sec)</th></tr></thead><tbody><tr><td colspan="7">Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..</td></tr><tr><td>1</td><td>Load Samples^a</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>2</td><td>Add BloodPrep DNA Purification Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>3</td><td>Add BloodPrep DNA Wash Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>4</td><td>Add BloodPrep DNA Wash Solution</td><td>600</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>5</td><td>Add BloodPrep DNA Wash Solution</td><td>300</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>9</td><td>Pre-Elution Vacuum</td><td>–</td><td>Waste</td><td>0</td><td>100</td><td>120</td></tr><tr><td>10</td><td>Touch Off</td><td>–</td><td>Waste</td><td>–</td><td>–</td><td>–</td></tr><tr><td>11</td><td>BloodPrep DNA Elution Solution 1^b</td><td>100</td><td>Collection</td><td>180</td><td>60</td><td>120</td></tr><tr><td>12</td><td>BloodPrep DNA Elution Solution 2^c</td><td>100</td><td>Collection</td><td>0</td><td>60</td><td>120</td></tr><tr><td>13</td><td>Touch Off</td><td>–</td><td>Collection</td><td>–</td><td>–</td><td>–</td></tr><tr><td colspan="7">a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.</td></tr><tr><td colspan="7">b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.</td></tr><tr><td colspan="7">c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.</td></tr><tr><td rowspan="2">2</td><td rowspan="2">Perform the purification run</td><td>a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.</td></tr><tr><td>b. Start the run.</td></tr></tbody></table>	Step	Description	Volume (μL)	Position	Incubation (sec)	Vacuum (%)	Time (sec)	Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..							1	Load Samples ^a	650	Waste	0	80	300	2	Add BloodPrep DNA Purification Solution	650	Waste	0	80	300	3	Add BloodPrep DNA Wash Solution	650	Waste	0	80	60	4	Add BloodPrep DNA Wash Solution	600	Waste	0	80	60	5	Add BloodPrep DNA Wash Solution	300	Waste	0	80	60	9	Pre-Elution Vacuum	–	Waste	0	100	120	10	Touch Off	–	Waste	–	–	–	11	BloodPrep DNA Elution Solution 1 ^b	100	Collection	180	60	120	12	BloodPrep DNA Elution Solution 2 ^c	100	Collection	0	60	120	13	Touch Off	–	Collection	–	–	–	a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.							b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.							c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.							2	Perform the purification run	a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.	b. Start the run.
		Step	Description	Volume (μL)	Position	Incubation (sec)	Vacuum (%)	Time (sec)																																																																																																																																																																																																																				
		Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the 6100 waste station.																																																																																																																																																																																																																										
		1	Load Samples ^a	650	Waste	0	80	300																																																																																																																																																																																																																				
		2	Add BloodPrep DNA Purification Solution ^a	650	Waste	0	80	400																																																																																																																																																																																																																				
		3	Add BloodPrep DNA Wash Solution	650	Waste	0	80	60																																																																																																																																																																																																																				
		4	Add BloodPrep DNA Wash Solution	600	Waste	0	80	60																																																																																																																																																																																																																				
		5	Add BloodPrep DNA Wash Solution	300	Waste	0	80	60																																																																																																																																																																																																																				
		6	Pre-Elution Vacuum	–	Waste	0	100	120																																																																																																																																																																																																																				
7	Touch Off	–	Waste	–	–	–																																																																																																																																																																																																																						
8	BloodPrep DNA Elution Solution 1 ^b	100	Collection	180	60	120																																																																																																																																																																																																																						
9	BloodPrep DNA Elution Solution 2 ^c	100	Collection	0	60	120																																																																																																																																																																																																																						
10	Touch Off	–	Collection	–	–	–																																																																																																																																																																																																																						
a.If the well evacuates slowly, increase the vacuum setting to 100% and repeat the step.																																																																																																																																																																																																																												
b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. You may lower the total elution volume to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.																																																																																																																																																																																																																												
c.It is very important that you use the elution solutions in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.																																																																																																																																																																																																																												
	Set up the 6100 PrepStation for tissue culture or buccal swab samples	Set up the DNA purification parameters for the blood sample on the protocol tab of the 6100 PrepStation software:																																																																																																																																																																																																																										
		<table><thead><tr><th>Step</th><th>Description</th><th>Volume (μL)</th><th>Position</th><th>Incubation (sec)</th><th>Vacuum (%)</th><th>Time (sec)</th></tr></thead><tbody><tr><td colspan="7">Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..</td></tr><tr><td>1</td><td>Load Samples^a</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>2</td><td>Add BloodPrep DNA Purification Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>300</td></tr><tr><td>3</td><td>Add BloodPrep DNA Wash Solution</td><td>650</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>4</td><td>Add BloodPrep DNA Wash Solution</td><td>600</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>5</td><td>Add BloodPrep DNA Wash Solution</td><td>300</td><td>Waste</td><td>0</td><td>80</td><td>60</td></tr><tr><td>9</td><td>Pre-Elution Vacuum</td><td>–</td><td>Waste</td><td>0</td><td>100</td><td>120</td></tr><tr><td>10</td><td>Touch Off</td><td>–</td><td>Waste</td><td>–</td><td>–</td><td>–</td></tr><tr><td>11</td><td>BloodPrep DNA Elution Solution 1^b</td><td>100</td><td>Collection</td><td>180</td><td>60</td><td>120</td></tr><tr><td>12</td><td>BloodPrep DNA Elution Solution 2^c</td><td>100</td><td>Collection</td><td>0</td><td>60</td><td>120</td></tr><tr><td>13</td><td>Touch Off</td><td>–</td><td>Collection</td><td>–</td><td>–</td><td>–</td></tr><tr><td colspan="7">a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.</td></tr><tr><td colspan="7">b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.</td></tr><tr><td colspan="7">c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.</td></tr><tr><td rowspan="2">2</td><td rowspan="2">Perform the purification run</td><td>a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.</td></tr><tr><td>b. Start the run.</td></tr></tbody></table>	Step	Description	Volume (μL)	Position	Incubation (sec)	Vacuum (%)	Time (sec)	Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..							1	Load Samples ^a	650	Waste	0	80	300	2	Add BloodPrep DNA Purification Solution	650	Waste	0	80	300	3	Add BloodPrep DNA Wash Solution	650	Waste	0	80	60	4	Add BloodPrep DNA Wash Solution	600	Waste	0	80	60	5	Add BloodPrep DNA Wash Solution	300	Waste	0	80	60	9	Pre-Elution Vacuum	–	Waste	0	100	120	10	Touch Off	–	Waste	–	–	–	11	BloodPrep DNA Elution Solution 1 ^b	100	Collection	180	60	120	12	BloodPrep DNA Elution Solution 2 ^c	100	Collection	0	60	120	13	Touch Off	–	Collection	–	–	–	a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.							b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.							c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.							2	Perform the purification run	a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.	b. Start the run.																																																																																																													
		Step	Description	Volume (μL)	Position	Incubation (sec)	Vacuum (%)	Time (sec)																																																																																																																																																																																																																				
		Note: Tape over empty wells of the purification tray with adhesive tape or an adhesive tray cover, or pre-wet all empty wells with 50 μL of BloodPrep DNA Purification Solution to ensure even vacuum. Position the carriage with the purification tray over the waste station..																																																																																																																																																																																																																										
		1	Load Samples ^a	650	Waste	0	80	300																																																																																																																																																																																																																				
		2	Add BloodPrep DNA Purification Solution	650	Waste	0	80	300																																																																																																																																																																																																																				
		3	Add BloodPrep DNA Wash Solution	650	Waste	0	80	60																																																																																																																																																																																																																				
		4	Add BloodPrep DNA Wash Solution	600	Waste	0	80	60																																																																																																																																																																																																																				
		5	Add BloodPrep DNA Wash Solution	300	Waste	0	80	60																																																																																																																																																																																																																				
		9	Pre-Elution Vacuum	–	Waste	0	100	120																																																																																																																																																																																																																				
10	Touch Off	–	Waste	–	–	–																																																																																																																																																																																																																						
11	BloodPrep DNA Elution Solution 1 ^b	100	Collection	180	60	120																																																																																																																																																																																																																						
12	BloodPrep DNA Elution Solution 2 ^c	100	Collection	0	60	120																																																																																																																																																																																																																						
13	Touch Off	–	Collection	–	–	–																																																																																																																																																																																																																						
a.For lysates with volumes in excess of 650 μL, use the Quick Run feature to pull 650 μL aliquots of lysate across the purification tray membrane, operating vacuum at 80% for 180 secs. Repeat until one aliquot remains to be added and then proceed to step 1 of the purification protocol. If the well starts to evacuate slowly after a number of additions, increase the vacuum setting to 100% and repeat the step.																																																																																																																																																																																																																												
b.Elution volume is between 75 μL and 200 μL in a standard PCR microplate. Elution volumes below 75 μL result in lowered yields of DNA. The total elution volume may be lowered to 75 μL, but the volumes of Elution Solution 1 and 2 must remain equal.																																																																																																																																																																																																																												
c.It is very important that the elution solutions are used in the correct order. Elution Solution 1 must be incubated with the DNA on the membrane for 3 minutes to ensure maximum yield. Then an equal volume of Elution Solution 2 must follow to give the correct pH for DNA storage.																																																																																																																																																																																																																												
2	Perform the purification run	a. Pipette the lysate samples into the purification tray. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 μL.																																																																																																																																																																																																																										
		b. Start the run.																																																																																																																																																																																																																										

DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs on the ABI PRISM™ 6100 Nucleic Acid PrepStation

© Copyright 2003, Applied Biosystems. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document. This document is believed to be complete and accurate at the time of publication. In no event shall Applied Biosystems be liable for incidental, special, multiple, or consequential damages in connection with or arising from the use of this document.

Applied Biosystems is a registered trademark of Applied Biosystems Corporation or its subsidiaries in the U.S. and/or certain other countries.

AB (Design), ABI PRISM, Applied Biosystems, and BloodPrep are trademarks of Applied Biosystems Corporation or its subsidiaries in the U.S. and/or certain other countries.

Printed in the USA. 05/2003

PN 4345297, Rev. A