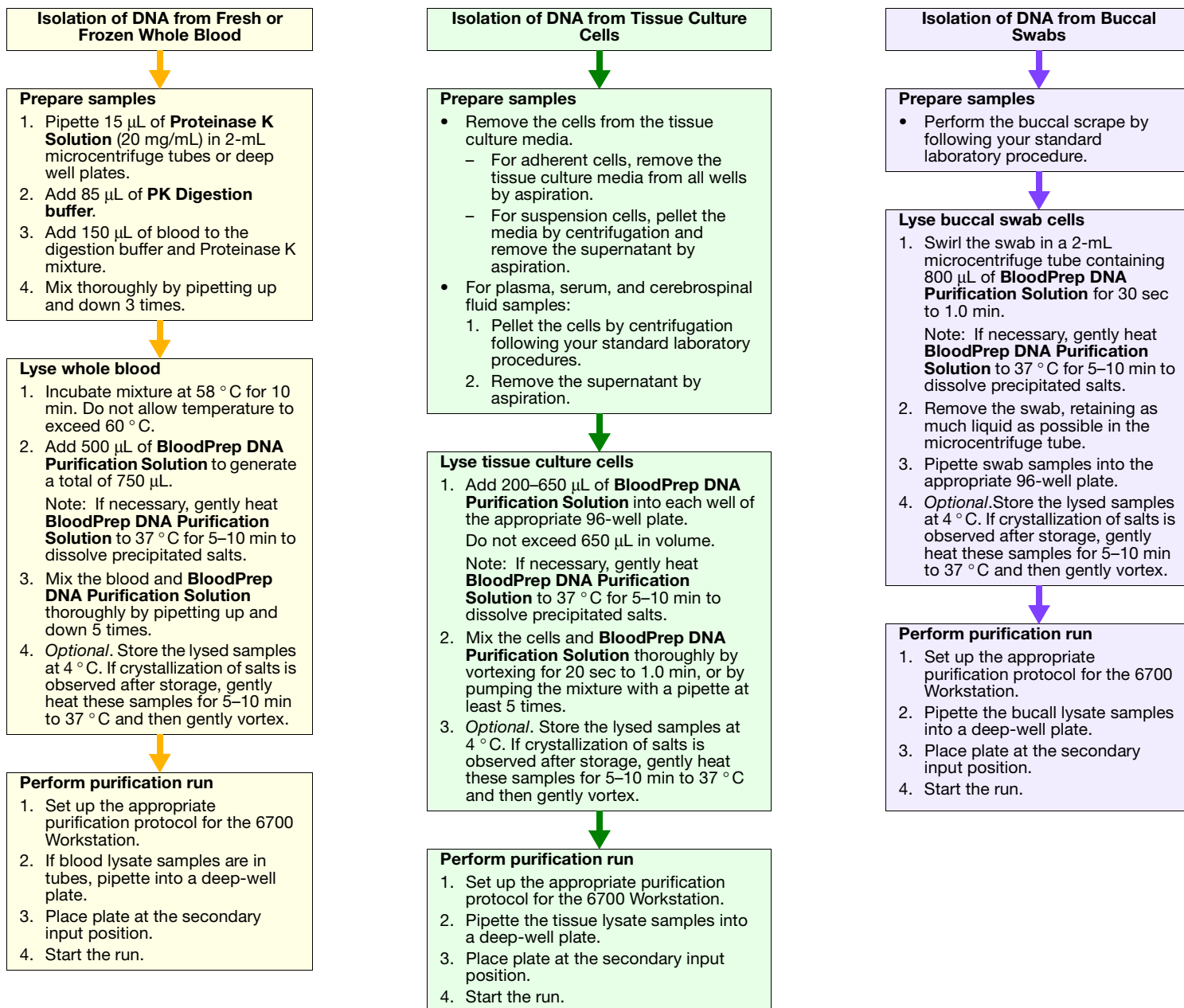


DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs on the ABI PRISM™ 6700 Nucleic Acid Workstation

⚠ 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

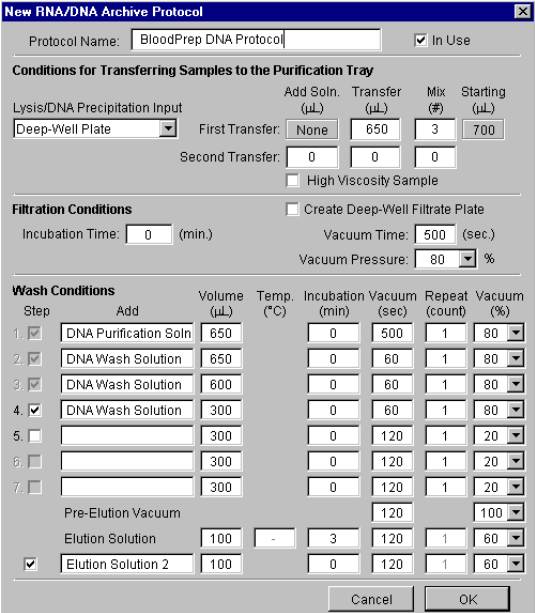
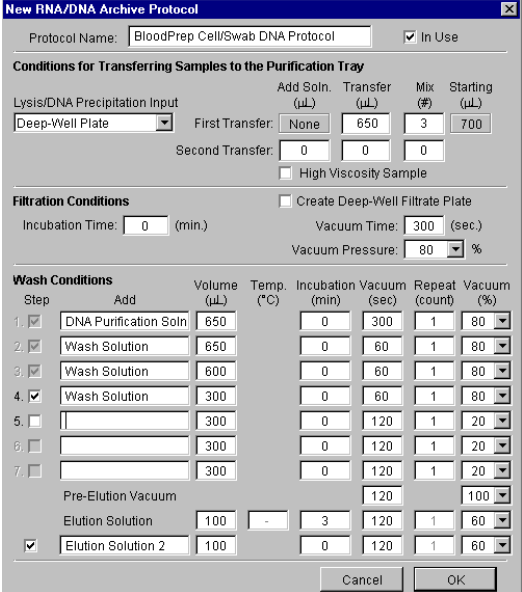
a. Major Laboratory Supplier

DNA Isolation from Fresh and Frozen Blood, Tissue Culture

Cells, and Buccal Swabs on the ABI PRISM™ 6700 Nucleic Acid Workstation *(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) in 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 appropriate 96-well plate.</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 6700 Workstation
1	Set up the 6700 Workstation	<p>Set up the appropriate DNA purification parameters for your sample on the protocol tab of the 6700 Workstation software.</p> <p>⚠ WARNING Read the MSDSs for BloodPrep DNA Purification Solution, BloodPrep DNA Wash Solution, and BloodPrep DNA Elution Solution 1 and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.</p> <p>Fresh or Frozen Whole Blood Parameters for the 6700 Workstation</p>  <p>Tissue Culture Cell or Buccal Swab Cell Parameters</p> 
2	Perform the purification run	<ol style="list-style-type: none"> Pipette the lysate samples into a deep-well plate. Place the plate at the secondary input position. Note: The maximum volume of lysate that the total DNA purification tray can accommodate is 600 µL. Start the run.

DNA Isolation from Fresh and Frozen Blood, Tissue Culture Cells, and Buccal Swabs on the ABI PRISM™ 6700 Nucleic Acid Workstation

© 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.

All other trademarks are the sole property of their respective owners.

Printed in the USA. 05/2003

PN 4345295, Rev. A