



Genomics Reagent Selection Guide

- = Protocol detailed in product's Instructions For Use or Manual
- = Supplemental data available from Beckman Coulter Life Sciences
- △ = Customer-published data available

Application:		DNA Extraction				RNA Extraction				DNA & RNA Extraction		Cleanup and Size Selection					
		GenFind V3 kit	DNAdvance kit	FormaPure XL DNA kit	CosMCPrep kit	RNAAdvance Blood kit	RNAAdvance Tissue kit	RNAAdvance Cell V2 kit	RNAAdvance Viral kit	FormaPure XL RNA kit	FormaPure XL Total kit	Mixed kit components	AMPure XP reagent	SPRIselect reagent	RNAClean XP reagent	CleanSEQ reagent	
Input Sample Type																	
Biofluids	Blood, fresh or frozen whole	•				△											
	EDTA-preserved blood	•															
	Citrate-preserved blood	•															
	Heparin-preserved blood	•															
	PAXgene™-preserved blood	○				•											
	RNAgard®-preserved blood					△											
	Blood spots, dried	○															
	Archived neonatal blood spots	△															
	Filter paper	○															
	FTA cards	○															
	Buccal swabs		○														
	Cells isolated from blood samples	△															
	Circulating tumor cells (CTCs)							△									
	Hematopoietic progenitor cells (HPCs)	△															
	Hematopoietic stem cells (HSCs)	△															
	PBMCs	△															
	Mouthwash	○															
	Naso- or Oropharyngeal Swabs								•								
	Saliva	○	○						•								
	Serum	•															



© 2025 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

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

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
2024-GBL-EN-105996-V2

