



# GENOMIC PRODUCTS PORTFOLIO OVERVIEW

## Advancing science is in our DNA

From the Human Genome Project to today's quest for precision medicine, we've long invested in providing genomic products that reduce complexity and improve productivity.

Our reagent portfolio is powered by Solid Phase Reversible Immobilization (SPRI) technology, widely known for use in our Agencourt AMPure XP, which uses paramagnetic beads to selectively bind nucleic acids by type and size. SPRI enables our chemistries to deliver high-performance isolation, purification and clean-up protocols supporting applications such as qPCR, ddPCR, Sanger sequencing, next-generation sequencing (NGS) and microarrays.

Our chemistries can be used interchangeably between manual methods and automated methods on Biomek liquid handlers, combining optimum performance with unprecedented flexibility.

### SOLUTIONS FOR:

- DNA Isolation
- RNA Isolation
- Total Nucleic Acid Isolation
- Purification & Clean-up

### SUPPORTED APPLICATIONS:

- Next-Generation Sequencing
- PCR/qPCR/ddPCR
- Sanger Sequencing
- Enzymatic post-reaction clean-ups





# Selection Guide

STARTING MATERIAL		DNA ISOLATION				RNA ISOLATION				PURIFICATION & CLEAN-UP			
		CosMC Prep	DNAdvance	FormaPure DNA <sup>1</sup>	GenFind v2	FormaPure	RNAdvance Blood	RNAdvance Cell v2	RNAdvance Tissue	AMPure XP	CleanSeq	RNAClean XP	SPRIselect
MAMMALIAN SAMPLES	Buccal Swabs		■										
	FFPE Tissue			■		■							
	FTA Cards				■								
	Oragene Saliva Collection Tubes				■								
	Plasma				■								
	Serum				■								
	Tissue		■						■				
	Whole Blood				■		■						
CELL CULTURES	293 T				■			■					
	Adherent				■			■					
	CHO				■			■					
	HeLa				■			■					
	Hepatocytes <50k Cells				■			■					
	Hepatocytes ≥50k Cells				■			■					
	Non-Adherent				■			■					
BACTERIA	<i>B. subtilis</i> (Gram +)				■			■	■				
	<i>E. coli</i> (Gram -)				■			■	■				
PLASMIDS	BACs, Cosmids, Fosmids	■											
OTHER MATERIALS	Avian Blood				■								
	Insects (including Drosophila)		■					■	■				
	Zebra Fish		■					■					
NUCLEIC ACID SAMPLES	cDNA Synthesis										■	■	
	Dye Terminator Removal									■	■	■	
	In vitro Transcription										■	■	
	NGS Library Construction Steps								■		■	■	
	PCR Products								■				

**DON'T SEE WHAT YOU'RE LOOKING FOR?** Please contact your local genomics specialist with questions. Our products can be optimized for individual needs, and new protocols are developed regularly.



1. FormaPure DNA can be optimized for Total Nucleic Acid Isolation

© 2017 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. Genomic reagents and kits are not intended or validated for use in the diagnosis of disease or other conditions.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at [beckman.com](http://beckman.com)

AAG-2765FLY06.17