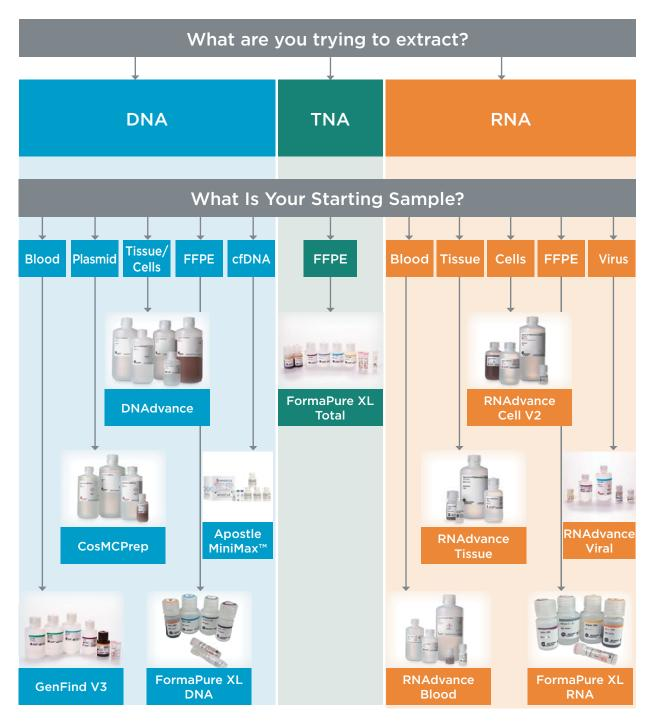


DON'T LOSE CRITICAL DATA EVALUATE BECKMAN COULTER'S EXTRACTION KITS

No matter how advanced the technology becomes, sequencing results are always impacted by sample quality. It's important to optimize every, equally-important step of your sample prep workflow starting with nucleic acid extraction.

Do you want to get more data out of your extraction? Are you looking to evaluate better, scalable reagent solutions? Review the chart below to determine which reagent kit can best support your workflow based on your extraction needs and starting samples.



© 2022 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. All other trademarks are the property of their respective owners.