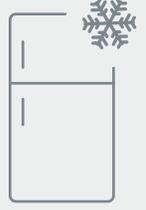


# Selecting the right reagent

Here's a quick guide to help you choose the right reagent for your genomic application.

REAGENT	NUCLEIC ACID INPUT	STORAGE TEMPERATURE	KEY DIFFERENTIATOR	PART NUMBERS & VOLUMES
 <b>AMPURE XP REAGENT</b>	 Works with DNA	 4°C	Known as the gold standard and suggested in over 200 library preparation kits, including those from the industry's most trusted sequencing companies.	A63880, 5 mL A63881, 60 mL A63882, 450 mL
 <b>SPRISELECT REAGENT</b>	 Works with fragmented DNA	 Room Temperature	Quality controlled for size selection, and manufacturing practices to limit lot-to-lot variation.	B23317, 5 mL B23318, 60 mL B23319, 450 mL
 <b>RNACLEAN XP REAGENT</b>	 Works with RNA and cDNA	 4°C	The only cleanup kit that has non detectable levels of RNase.	A63987, 40 mL A66514, 450 mL

Not intended or validated for use in the diagnosis of disease or other conditions.