

# Verified Agilent Bravo NGS SureSelect Automation Protocols

Agilent NGS Reagents	Compatible Panels	Option*	Reagent User Guide
SureSelect XT HS2 DNA	SureSelect Human All Exon including V8, pre-designed probes plus custom, and SureSelect Custom probes	A	<a href="#">G9985-90020</a>
		B	<a href="#">G9985-90010</a>
SureSelect XT HS2 RNA (Targeted)		A	<a href="#">G9993-90020</a>
		B	<a href="#">G9993-90010</a>
SureSelect <sup>XT</sup>		B	<a href="#">G7550-90000</a>
SureSelect <sup>XT2</sup>	SureSelect Human All Exon, pre-designed probes plus custom, and SureSelect Custom probes	B	<a href="#">G9450-90000</a>
SureSelect <sup>QXT</sup>	N/A – whole genome library prep	A	<a href="#">G9682-90020</a>
		B	<a href="#">G9682-90010</a>
SureSelect <sup>XT</sup> mRNA Target Enrichment	Automated poly-A selection + target enrichment (SureSelect Human All Exon, pre-designed probes plus custom, SureSelect Custom probes)	B	<a href="#">G9691-90020</a>
SureSelect XT HS2 mRNA	N/A – Automated poly-A selection	B	<a href="#">G9995-90010</a>
SureSelect <sup>XT</sup> Strand-specific RNA Library Prep		B	<a href="#">G9691-90030</a>
SureSelect <sup>XT</sup> Low Input	SureSelect Human All Exon including V8, pre-designed probes plus custom, and SureSelect Custom probes	A	<a href="#">G9703-90020</a>
SureSelect XT HS2 DNA	DNA Library Preparation, Pre-Capture Pooling (optional), and Target Enrichment	B+	<a href="#">G9985-90015</a>
SureSelect Max DNA-Enzymatic Fragmentation	SureSelect Human All Exon including V8, SureSelect Cancer CGP, pre-designed probes plus custom, and SureSelect custom probes	B	<a href="#">G9660-90010</a>
SureSelect Max DNA-Sheared DNA	SureSelect Human All Exon including V8, SureSelect Cancer CGP, pre-designed probes plus custom, and SureSelect custom probes	B	<a href="#">G9663-90010</a>

\*Bravo NGS (Option A) Bravo NGS Workstations (Options B) Bravo NGS Workstation with ODTC (Option B+)