

Item Number	Custom oligonucleotides Special quote
	<b>Special quote HTP oligonucleotides</b>
SQ-HTPOL-XXXX	Special Quote - HTP Oligos
	<b>Special quote RNA oligonucleotides</b>
SQ-RNAOL-XXXX	Special Quote - RNA Oligos
	<b>Special quote siRNA</b>
SQ-SIRNA-XXXX	Special Quote - siRNA Duplexes
	<b>Special quote Large Scale oligonucleotides</b>
SQ-LASOL-XXXX	Special Quote - Large Scale Oligos
	<b>Special quote Medium Scale oligonucleotides</b>
SQ-MESOL-XXXX	Special Quote - Medium Scale Oligos
	<b>Special quote modified oligonucleotides</b>
SQ-MODOL-XXXX	Special Quote - Modified Oligos
	<b>Special quote PNA oligonucleotides</b>
SQ-PNAOL-XXXX	Special Quote - PNA Oligos
	<b>Special quote RealTime qPCR probes</b>
SQ-PROBE-XXXX	Special Quote - qPCR Oligos