

Oligonucleotide Quality Control

According to the oligo specifications, we perform quality controls for free as reported in the table below.

Oligo Type		MS ¹	UHPLC
Custom Oligonucleotides	Unmodified	✓ ²	
	Modified	✓ ³	
	UltraPureGold™	✓	✓
qPCR Probes	Double Dye Probes	✓	✓
	Molecular Beacons	✓	✓
	MGB Probes	✓	✓
RNAi Oligonucleotides	siRNA Duplexes	✓	
NGS Oligonucleotides		✓	✓
Complex Oligonucleotides		✓	✓ ⁴
Calibration Oligos		✓	✓
Universal Primers		✓	✓

MS: Mass Spectrometry; UHPLC: Ultra Performance Liquid Chromatography.

¹Always provided up to 60 bases long Oligonucleotides.

²Randomized high throughput control

³Except for SePOP desalted oligonucleotides.

⁴Optional.

For technical reasons this general rule may be adapted to provide you with the most suitable accurate oligonucleotide.