

# Oligonucleotide Quality Control

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

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

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

<sup>1</sup>Always provided up to 60 bases long Oligonucleotides.

<sup>2</sup>Randomized high throughput control

<sup>3</sup>Except for SePOP desalted oligonucleotides.

<sup>4</sup>Optional.

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