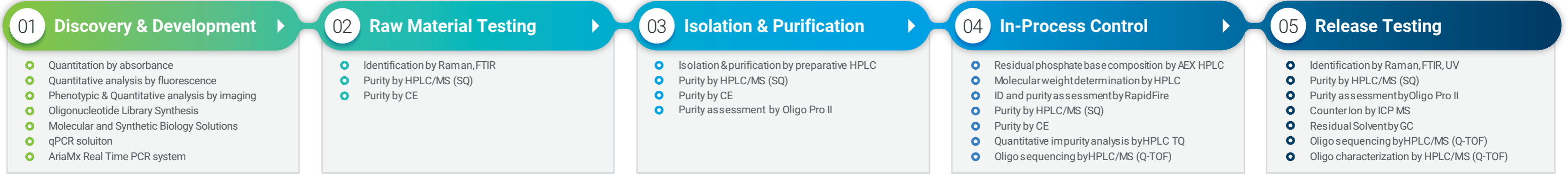


## Oligo Bio-workflow



## 응용자료 가이드: Method and System Overview

응용자료 이름 및 분석장비를 클릭하여 자세한 사항을 확인하세요

Workflow	Analytical Tools	Application Name	Bio LCs/Conventional LCs		LCMS	Automation	GC	CE	ICP	Molecular Spectroscopy	Microplate Reader	Cell Imaging System	PCR	Oligonucleotide Synthesis	Gene Cloning	Consumable		
			UHPLC	HPLC	2D-LC	Prep LC	SQ, TQ, QTOF	RapidFire	GC, HSS	CE, Oligo Pro II	ICP-MS	UV, FTIR, Raman	Epoch	Synergy	Cytation	qPCR	Sure Print	Gene Cloning
Discovery & Development	Quantitation by absorbance	Analytical Performance of Nucleic Acid Microvolume Quantification Using the Epoch Spectrophotometer System															-	
	Quantitative analysis by fluorescence	Fluorometric Quantitation of dsDNA using PicoGreen															-	
	Phenotypic & Quantitative analysis by imaging	Quantification of RNA using Microplate-based UV and Fluorescence Methods															-	
	Oligonucleotide Library Synthesis	Multiplexed Detection of Cytokine Cancer Biomarkers using Fluorescence RNA In Situ Hybridization and Cellular Imaging															-	
	Molecular and Synthetic Biology Solutions	Sureprint Oligonucleotide Libraries															-	
	qPCR solution	cloning, mutagenesis, viral expression system and competent cells by Agilent															-	
	AriaMx Real Time PCR system	Agilent solution for qPCR															-	
Raw Material Testing	Identification by Raman, FTIR	Realtime quantification of target gene															-	
		Identification of Oligonucleotide Starting Materials Directly Through Containers															-	
		Achieving Faster Ramp Rates for Thermal Melt Experiments															-	
		Best Practice for Nucleic Acid Thermal Stability Measurements using 3500 UV-Vis															-	
	Purity by HPLC/MS (SQ)	Analyzing Raw Material for Oligonucleotide Synthesis															ZORBAX RRHD Eclipse Plus C18	
Isolation & Purification	Purity by HPLC/MS (SQ)	Comparability Studies for the Analysis of Nucleotides on Four Different LC Systems															Poroshell 120 HILIC-Z, PEEK-lined	
		Polar-Modified Stationary Phases are Ideal for the Analysis of Nucleotides															Polaris C18-A	
	Purity by CE	Oligonucleotides—analysis of antisense therapeutics															Capillary (PVA)	
		Analysis of Oligonucleotides by Capillary Gel Electrophoresis															Capillaries (PVA μSIL WAX, μSIL DNA)	
	Isolation & purification by preparative HPLC	Purification of single-stranded RNA oligonucleotides using HPLC															PLRP-S	
		Generating Pure Oligonucleotides with PLRP-S															PLRP-S	
		Purification your way: siRNA to mRNA															PLRP-S, PL-SAX	
	Purity by HPLC/MS (SQ)	High performance DNA oligonucleotide purification using Agilent TOP-DNA															TOP-DNA	
		Evaluation of different ion-pairing reagents of oligonucleotides																AdvanceBio oligonucleotide
		Fast impurity profiling of synthetic oligonucleotides with LCQTOF																ZORBAX Eclipse Plus C18
In-Process Control	Purity by HPLC/MS (SQ)	High resolution separation of oligonucleotides using IEX column															PL-SAX 1000Å	
		Oligonucleotides—analysis of antisense therapeutics															Capillary (PVA)	
	Purity by CE	Analysis of Oligonucleotides by Capillary Gel Electrophoresis															Capillaries (PVA μSIL WAX, μSIL DNA)	
	Purity assessment by Oligo Pro II	Accurate Assessment of Oligonucleotide Purity															-	
	Residual phosphate base composition by HPLC	Comparability Studies for the Analysis of Nucleotides on Four Different LC Systems															Poroshell 120 HILIC-Z, PEEK-lined	
	Molecular weight determination by HPLC	Determination of Molecular Weight and Size Using the 1260 MDS with LSD															Bio SEC-3	
	ID and purity assessment by RapidFire	High-throughput Mass Spectrometry Analysis of Synthetic Oligonucleotides															AdvanceBio oligonucleotide	
		Direct Analysis of In-Process Oligonucleotides Without Manual Purification																PLRP-S, AdvanceBio Oligonucleotide
	Purity by HPLC/MS (SQ)	Evaluation of different ion-pairing reagents of oligonucleotides																AdvanceBio oligonucleotide
		Fast impurity profiling of synthetic oligonucleotides with LCQTOF																ZORBAX Eclipse Plus C18
Release Testing	Purity by HPLC/MS (SQ)	High resolution separation of oligonucleotides using IEX column															PL-SAX 1000Å	
		Fast and High-Resolution Reversed-Phase Separation of Synthetic Oligonucleotides															AdvanceBio oligonucleotide	
	Purity by CE	Oligonucleotides—analysis of antisense therapeutics															Capillary (PVA)	
		Analysis of Oligonucleotides by Capillary Gel Electrophoresis															Capillaries (PVA μSIL WAX, μSIL DNA)	
	Quantitative impurity analysis by HPLC-TQ	Quantification Using LCQTOF															-	
	Oligo sequencing by HPLC/MS (Q-TOF)	Analysis of mRNA Poly-A sequence variants by High-Resolution LC/MS															Poroshell 120 HPH-C18, PLRP-S	
	Identification by Raman, FTIR, UV	Measuring raw materials without sampling—directly through their containers															-	
Purity by HPLC/MS (SQ)	Fast impurity profiling of synthetic oligonucleotides with LCQTOF															ZORBAX Eclipse Plus C18		
	High resolution separation of oligonucleotides using IEX column															PL-SAX 1000Å		
	Evaluation of different ion-pairing reagents of oligonucleotides																AdvanceBio oligonucleotide	
	Performance Characteristics of the Agilent Oligo Pro II System																-	
Purity assessment by Oligo Pro II	Validating performance of an Agilent ICP-MS for USP <232>/<233> & ICH Q3D/Q2(R1)																-	
Counter ion by ICP-MS	Analysis of USP Method <467> Residual Solvents on the Agilent 8890 GC System																J&W DB-Select 624 UI, J&W HP-INNOWax	
Residual solvent by GC	Analysis of mRNA Poly-A sequence variants by High-Resolution LC/MS																Poroshell 120 HPH-C18, PLRP-S	
Oligo sequencing by HPLC/MS (Q-TOF)	Rapid Analysis of mRNA 5' Capping with High Resolution LC/MS																AdvanceBio Oligonucleotide	
Oligo Characterization by HPLC/MS (Q-TOF)	Mass Spectrometric Characterization of Antibody-siRNA Conjugates using the LC/Q-TOF																PLRP-S, AdvanceBio SEC	

## Agilent solutions

제품 사진 및 이름을 클릭하여 더 상세한 내용 확인하세요  
제품과 관련된 상담 및 견적은 영업지원 또는 제품 페이지에서 24시간 언제나 가능합니다.

Bio LC	Conventional LC	Quadrupole Time of Flight LC/MS	High Throughput LC/MS	CE & CE/MS Systems	Automated Electrophoresis	Cell Imaging Multimode Readers	Atomic Spectroscopy	GC System	Columns and supplies
애질런트 바이오 LC 시스템에 대해 더 알아보기	애질런트 LC 시스템에 대해 더 알아보기	6545XT AdvanceBio LC/Q-TOF	RapidFire 400	7100 CE	Oligo Pro II	Cytation 5	7850 ICP-MS	8890 GC	Oligonucleotide Column

더 많은 Agilent Oligonucleotide 워크플로우 응용 및 관련 자료 보기