

## Custom oligonucleotides Purifications

Item Number	Purifications for custom oligonucleotides Single purifications
PU-GOLD0-001	RP-Cartridge - GoldTM 10 nmol
PU-GOLD0-004	RP-Cartridge - GoldTM 40 nmol
PU-GOLD0-020	RP-Cartridge - GoldTM 200 nmol
PU-GOLD0-100	RP-Cartridge - GoldTM 1000 nmol
PU-HPLIE-004	Ion-Exchange HPLC (IEX-HPLC) 40 nmol
PU-HPLIE-020	Ion-Exchange HPLC (IEX-HPLC) 200 nmol
PU-HPLIE-100	Ion-Exchange HPLC (IEX-HPLC) 1000 nmol
PU-HPLIE-M02	Ion-Exchange HPLC (IEX-HPLC) 2,5 µmol
PU-HPLIE-M05	Ion-Exchange HPLC (IEX-HPLC) 5 µmol
PU-HPLIE-M10	Ion-Exchange HPLC (IEX-HPLC) 10 µmol
PU-HPLRP-004	Reverse Phase HPLC (RP-HPLC) 40 nmol
PU-HPLRP-020	Reverse Phase HPLC (RP-HPLC) 200 nmol
PU-HPLRP-100	Reverse Phase HPLC (RP-HPLC) 1000 nmol
PU-HPLRP-M02	Reverse Phase HPLC (RP-HPLC) 2,5 µmol
PU-HPLRP-M05	Reverse Phase HPLC (RP-HPLC) 5 µmol
PU-HPLRP-M10	Reverse Phase HPLC (RP-HPLC) 10 µmol
PU-PAGE0-004	PolyAcrylamide Gel Electrophoresis (PAGE) 40 nmol
PU-PAGE0-020	PolyAcrylamide Gel Electrophoresis (PAGE) 200 nmol
PU-PAGE0-100	PolyAcrylamide Gel Electrophoresis (PAGE) 1000 nmol
PU-PAGE0-M02	PolyAcrylamide Gel Electrophoresis (PAGE) 2,5 µmol
PU-PAGE0-M05	PolyAcrylamide Gel Electrophoresis (PAGE) 5 µmol
PU-PAGE0-M10	PolyAcrylamide Gel Electrophoresis (PAGE) 10 µmol
PU-SG025-004	Sephadex G-25 40 nmol
PU-SG025-020	Sephadex G-25 200 nmol
PU-SG025-100	Sephadex G-25 1000 nmol
	Purifications for custom oligonucleotides Dual Purifications
PU-DRPIE-004	Dual HPLC (RP+IEX) 40 nmol
PU-DRPIE-020	Dual HPLC (RP+IEX) 200 nmol
PU-DRPIE-100	Dual HPLC (RP+IEX) 1000 nmol
PU-DRPIE-M02	Dual HPLC (RP+IEX) 2,5 µmol
PU-DRPIE-M05	Dual HPLC (RP+IEX) 5 µmol
PU-DRPIE-M10	Dual HPLC (RP+IEX) 10 µmol
PU-DRPLC-004	Dual HPLC (RP+RP) 40 nmol
PU-DRPLC-020	Dual HPLC (RP+RP) 200 nmol
PU-DRPLC-100	Dual HPLC (RP+RP) 1000 nmol
PU-DRPLC-M02	Dual HPLC (RP+RP) 2,5 µmol
PU-DRPLC-M05	Dual HPLC (RP+RP) 5 µmol
PU-DRPLC-M10	Dual HPLC (RP+RP) 10 µmol
PU-UPGD0-020	UltraPureGoldTM 200 nmol
PU-UPGD0-100	UltraPureGoldTM 1000 nmol