

Item Number	RNAi oligonucleotides
	<b>Custom siRNA Duplexes</b>
SR-HP001-001	siRNA Duplex HPLC 10 nmol
SR-HP001-004	siRNA Duplex HPLC 40 nmol
SR-HP001-020	siRNA Duplex HPLC 200 nmol
SR-HP001-100	siRNA Duplex HPLC 1000 nmol
SR-HP001-M02	siRNA Duplex HPLC 2,5 µmol
SR-HP001-M05	siRNA Duplex HPLC 5 µmol
SR-HP001-M10	siRNA Duplex HPLC 10 µmol
SR-NP001-001	siRNA Duplex SePOP 10 nmol
SR-NP001-004	siRNA Duplex SePOP 40 nmol
SR-NP001-020	siRNA Duplex SePOP 200 nmol
SR-NP001-100	siRNA Duplex SePOP 1000 nmol
SR-PA001-004	siRNA Duplex PAGE 40 nmol
SR-PA001-020	siRNA Duplex PAGE 200 nmol
SR-PA001-100	siRNA Duplex PAGE 1000 nmol
	<b>Control siRNA Duplexes</b>
SR-CL000-005	Control siRNA duplex negative control 5 nmol
SR-CL001-005	Control siRNA duplex LaminB1 (human) 5 nmol
SR-CL002-005	Control siRNA duplex Vimentin (human) 5 nmol
SR-CL003-005	Control siRNA duplex NuMA (human) 5 nmol
SR-CL004-005	Control siRNA duplex Beta-actin (human) 5 nmol
SR-CL005-005	Control siRNA duplex Eg-5 (human) 5 nmol
SR-CL006-005	Control siRNA duplex Cdk-1 (human) 5 nmol
SR-CL010-005	Control siRNA duplex pGL2 luciferase (firefly) 5 nmol
SR-CL011-005	Control siRNA duplex pGL3 luciferase (firefly) 5 nmol
SR-CL020-005	Control siRNA duplex GFP (jellyfish) 5 nmol
	<b>ClearMiR miRNA inhibitors</b>
MR-CLHSA-134A	Clear-miR hsa-miR34a 0.2 µmol
MR-CLHSA-146	Clear-MiR hsa-miR-146 0.2 µmol
MR-CLHSA-148b	Clear-MiR hsa-miR-148b 200 nmol
MR-CLHSA-152	Clear-MiR hsa-miR-152 200 nmol
MR-CLHSA-NEG	Clear-miR hsa-neg 0.2 µmol
MR-CLLET-7	Clear-MiR Plus hsa-let-7 200 nmol
MR-CLMMU-691	Clear-MiR mmu-miR-691 200 nmol
MR-CLMMU-709	Clear-MiR mmu-miR-709 200 nmol
MR-CPMMU-146a	Clear-MiR Plus mmu-miR-146a 5 nmol