

Item Number	Modifications for custom oligonucleotides Spacers
MD-SP003-03004	3' Spacer C3 40 nmol
MD-SP003-03020	3' Spacer C3 200 nmol
MD-SP003-03100	3' Spacer C3 1000 nmol
MD-SP003-IN004	Spacer C3 40 nmol
MD-SP003-IN020	Spacer C3 200 nmol
MD-SP003-IN100	Spacer C3 1000 nmol
MD-SP008-IN004	Spacer C8 (Octanediol) 40 nmol
MD-SP008-IN020	Spacer C8 (Octanediol) 200 nmol
MD-SP008-IN100	Spacer C8 (Octanediol) 1000 nmol
MD-SP009-IN004	Spacer C9 40 nmol
MD-SP009-IN020	Spacer C9 200 nmol
MD-SP009-IN100	Spacer C9 1000 nmol
MD-SP012-IN004	Spacer C12 40 nmol
MD-SP012-IN020	Spacer C12 200 nmol
MD-SP012-IN100	Spacer C12 1000 nmol
MD-SP018-IN004	Spacer C18 /HEG 40 nmol
MD-SP018-IN020	Spacer C18 /HEG 200 nmol
MD-SP018-IN100	Spacer C18 /HEG 1000 nmol
MD-SP025-IN004	THF (Tetrahydrofuran) / dSpacer (Abasic site) 40 nmol
MD-SP025-IN020	THF (Tetrahydrofuran) / dSpacer (Abasic site) 200 nmol
MD-SP025-IN100	THF (Tetrahydrofuran) / dSpacer (Abasic site) 1000 nmol