

Custom oligonucleotides Bases

Item Number	Custom oligonucleotides DNA Bases
BA-DN001-001	DNA Base 10 nmol (15-60 bases)
BA-DN001-004	DNA Base 40 nmol (10-100 bases)
BA-DN001-020	DNA Base 200 nmol (5-140 bases)
BA-DN001-100	DNA Base 1000 nmol (5-140 bases)
BA-DN001-M02	DNA Base 2,5 µmol (5-140 bases)
BA-DN001-M05	DNA Base 5 µmol (5-140 bases)
BA-DN001-M10	DNA Base 10 µmol (5-140 bases)
BA-DN010-004	DNA Base + Phosphorothioate Link 40 nmol
BA-DN010-020	DNA Base + Phosphorothioate Link 200 nmol
BA-DN010-100	DNA Base + Phosphorothioate Link 1000 nmol
BA-DN010-M02	DNA Base + Phosphorothioate Link 2,5 µmol
BA-DN010-M05	DNA Base + Phosphorothioate Link 5 µmol
BA-DN010-M10	DNA Base + Phosphorothioate Link 10 µmol
BA-DN020-004	DNA Base + Methylphosphonate Link 40 nmol
BA-DN020-020	DNA Base + Methylphosphonate Link 200 nmol
BA-DN020-100	DNA Base + Methylphosphonate Link 1000 nmol
Custom oligonucleotides RNA Bases	
BA-RN001-004	RNA Base 40 nmol (15-30 bases)
BA-RN001-020	RNA Base 200 nmol (8-80 bases)
BA-RN001-100	RNA Base 1000 nmol (8-80 bases)
BA-RN001-M02	RNA Base 2,5 µmol (8-80 bases)
BA-RN001-M05	RNA Base 5 µmol (8-80 bases)
BA-RN001-M10	RNA Base 10 µmol (8-80 bases)
BA-RN010-004	RNA Base + Phosphorothioate Link 40 nmol
BA-RN010-020	RNA Base + Phosphorothioate Link 200 nmol
BA-RN010-100	RNA Base + Phosphorothioate Link 1000 nmol
BA-RN010-M02	RNA Base + Phosphorothioate Link 2,5 µmol
BA-RN010-M05	RNA Base + Phosphorothioate Link 5 µmol
BA-RN010-M10	RNA Base + Phosphorothioate Link 10 µmol
Custom oligonucleotides LNA® Bases	
BA-LN001-004	LNA Base 40 nmol
BA-LN001-020	LNA Base 200 nmol
BA-LN001-100	LNA Base 1000 nmol
BA-LN010-004	LNA Base + Phosphorothioate Link 40 nmol
BA-LN010-020	LNA Base + Phosphorothioate Link 200 nmol
BA-LN010-100	LNA Base + Phosphorothioate Link 1000 nmol
Custom oligonucleotides 2' OMethyl RNA Bases	
BA-2M001-004	2' O-Me RNA Base 40 nmol (15-30 bases)
BA-2M001-020	2' O-Me RNA Base 200 nmol (8-80 bases)
BA-2M001-100	2' O-Me RNA Base 1000 nmol (8-80 bases)
BA-2M001-M02	2' O-Me RNA Base 2,5 µmol (8-80 bases)
BA-2M001-M05	2' O-Me RNA Base 5 µmol (8-80 bases)
BA-2M001-M10	2' O-Me RNA Base 10 µmol (8-80 bases)
BA-2M010-004	2' O-Me RNA Base + Phosphorothioate Link 40 nmol
BA-2M010-020	2' O-Me RNA Base + Phosphorothioate Link 200 nmol
BA-2M010-100	2' O-Me RNA Base + Phosphorothioate Link 1000 nmol
BA-2M010-M02	2' O-Me RNA Base + Phosphorothioate Link 2,5 µmol
BA-2M010-M05	2' O-Me RNA Base + Phosphorothioate Link 5 µmol
BA-2M010-M10	2' O-Me RNA Base + Phosphorothioate Link 10 µmol