

Item Number	RealTime qPCR oligonucleotides NASBA kits
NB-CC02-048	U1A cDNA Control
NB-CI01-048	Mix EasyQ® Control U1A (48 tests)
NB-CMV01-048	Mix EasyQ® Cytomegalovirus (48 tests)
NB-CT01-048	Mix EasyQ® Chlamydia Trachomatis(48 tests)
NB-DE01-048	Mix EasyQ® Dengue (48 tests)
NB-MT01-048	Mix EasyQ® Mycobacterium tuberculosis (48 tests)
NB-PAI01-048	Mix EasyQ® Parainfluenza 1 (48 tests)
NB-PAI02-048	Mix EasyQ® Parainfluenza 2 (48 tests)
NB-PAI03-048	Mix EasyQ® Parainfluenza 3 (48 tests)
NB-PAI04-048	Mix EasyQ® Parainfluenza 4 (48 tests)
NB-UC02-048	Mix EasyQ® Control U1A (48 tests)