

Item Number	DoubleDye Probes 3' BHQ™
PB-DD600-004	DD Probe 5' TAMRA + 3' BHQ-2™ 40 nmol
PB-DD600-020	DD Probe 5' TAMRA + 3' BHQ-2™ 200 nmol
PB-DD600-100	DD Probe 5' TAMRA + 3' BHQ-2™ 1000 nmol
PB-DD601-004	DD Probe 5' FAM + 3' BHQ-1™ 40 nmol
PB-DD601-020	DD Probe 5' FAM + 3' BHQ-1™ 200 nmol
PB-DD601-100	DD Probe 5' FAM + 3' BHQ-1™ 1000 nmol
PB-DD601-M05	DD Probe 5' FAM + 3' BHQ-1™ 5 µmol
PB-DD610-004	DD Probe 5' HEX + 3' BHQ-1™ 40 nmol
PB-DD610-020	DD Probe 5' HEX + 3' BHQ-1™ 200 nmol
PB-DD610-100	DD Probe 5' HEX + 3' BHQ-1™ 1000 nmol
PB-DD612-004	DD Probe 5' TET + 3' BHQ-1™ 40 nmol
PB-DD612-020	DD Probe 5' TET + 3' BHQ-1™ 200 nmol
PB-DD612-100	DD Probe 5' TET + 3' BHQ-1™ 1000 nmol
PB-DD614-004	DD Probe 5' JOE + 3' BHQ-1™ 40 nmol
PB-DD614-020	DD Probe 5' JOE + 3' BHQ-1™ 200 nmol
PB-DD614-100	DD Probe 5' JOE + 3' BHQ-1™ 1000 nmol
PB-DD616-004	DD Probe 5' ROX + 3' BHQ-2™ 40 nmol
PB-DD616-020	DD Probe 5' ROX + 3' BHQ-2™ 200 nmol
PB-DD616-100	DD Probe 5' ROX + 3' BHQ-2™ 1000 nmol
PB-DD620-004	DD Probe 5' Rhodamine-6G + 3' BHQ-1™ 40 nmol
PB-DD620-020	DD Probe 5' Rhodamine-6G + 3' BHQ-1™ 200 nmol
PB-DD620-100	DD Probe 5' Rhodamine-6G + 3' BHQ-1™ 1000 nmol
PB-DD630-004	DD Probe 5' Texas Red® + 3' BHQ-2™ 40 nmol
PB-DD630-020	DD Probe 5' Texas Red® + 3' BHQ-2™ 200 nmol
PB-DD630-100	DD Probe 5' Texas Red® + 3' BHQ-2™ 1000 nmol
PB-DD640-004	DD Probe 5' Yakima Yellow® + 3' BHQ-1™ 40 nmol
PB-DD640-020	DD Probe 5' Yakima Yellow® + 3' BHQ-1™ 200 nmol
PB-DD640-100	DD Probe 5' Yakima Yellow® + 3' BHQ-1™ 1000 nmol
PB-DD645-004	DD Probe 5' Dragonfly Orange™ + 3' BHQ-2™ 40 nmol
PB-DD645-020	DD Probe 5' Dragonfly Orange™ + 3' BHQ-2™ 200 nmol
PB-DD645-100	DD Probe 5' Dragonfly Orange™ + 3' BHQ-2™ 1000 nmol
PB-DD650-004	DD Probe 5' Cy® 3 + 3' BHQ-2™ 40 nmol
PB-DD650-020	DD Probe 5' Cy® 3 + 3' BHQ-2™ 200 nmol
PB-DD650-100	DD Probe 5' Cy® 3 + 3' BHQ-2™ 1000 nmol
PB-DD655-004	DD Probe 5' Cy5® + 3' BHQ-2™ 40 nmol
PB-DD655-020	DD Probe 5' Cy5® + 3' BHQ-2™ 200 nmol
PB-DD655-100	DD Probe 5' Cy5® + 3' BHQ-2™ 1000 nmol
PB-DD680-004	DD Probe 5' DYOMICS + 3' BHQ™ 40 nmol
PB-DD680-020	DD Probe 5' DYOMICS + 3' BHQ™ 200 nmol
PB-DD680-100	DD Probe 5' DYOMICS + 3' BHQ™ 1000 nmol
PB-DD690-004	DD Probe 5' ATTO + 3' BHQ™ 40 nmol
PB-DD690-020	DD Probe 5' ATTO + 3' BHQ™ 200 nmol
PB-DD690-100	DD Probe 5' ATTO + 3' BHQ™ 1000 nmol