

Item Number	DoubleDye Probes 3' TAMRA
PB-DD101-001	DD Probe 5' FAM + 3' TAMRA 10 nmol
PB-DD101-004	DD Probe 5' FAM + 3' TAMRA 40 nmol
PB-DD101-020	DD Probe 5' FAM + 3' TAMRA 200 nmol
PB-DD101-100	DD Probe 5' FAM + 3' TAMRA 1000 nmol
PB-DD110-004	DD Probe 5' HEX + 3' TAMRA 40 nmol
PB-DD110-020	DD Probe 5' HEX + 3' TAMRA 200 nmol
PB-DD110-100	DD Probe 5' HEX + 3' TAMRA 1000 nmol
PB-DD112-004	DD Probe 5' TET + 3' TAMRA 40 nmol
PB-DD112-020	DD Probe 5' TET + 3' TAMRA 200 nmol
PB-DD112-100	DD Probe 5' TET + 3' TAMRA 1000 nmol
PB-DD140-004	DD Probe 5' Yakima Yellow® + 3' TAMRA 40 nmol
PB-DD140-020	DD Probe 5' Yakima Yellow® + 3' TAMRA 200 nmol
PB-DD140-100	DD Probe 5' Yakima Yellow® + 3' TAMRA 1000 nmol