

Item Number	Molecular Beacons 3' Dabcyl
PB-MB200-004	M Beacon 5' TAMRA + 3' DABCYL 40 nmol
PB-MB200-020	M Beacon 5' TAMRA + 3' DABCYL 200 nmol
PB-MB200-100	M Beacon 5' TAMRA + 3' DABCYL 1000 nmol
PB-MB201-004	M Beacon 5' FAM + 3' DABCYL 40 nmol
PB-MB201-020	M Beacon 5' FAM + 3' DABCYL 200 nmol
PB-MB201-100	M Beacon 5' FAM + 3' DABCYL 1000 nmol
PB-MB210-004	M Beacon 5' HEX + 3' DABCYL 40 nmol
PB-MB210-020	M Beacon 5' HEX + 3' DABCYL 200 nmol
PB-MB210-100	M Beacon 5' HEX + 3' DABCYL 1000 nmol
PB-MB212-004	M Beacon 5' TET + 3' DABCYL 40 nmol
PB-MB212-020	M Beacon 5' TET + 3' DABCYL 200 nmol
PB-MB212-100	M Beacon 5' TET + 3' DABCYL 1000 nmol
PB-MB214-004	M Beacon 5' JOE + 3' DABCYL 40 nmol
PB-MB214-020	M Beacon 5' JOE + 3' DABCYL 200 nmol
PB-MB214-100	M Beacon 5' JOE + 3' DABCYL 1000 nmol
PB-MB216-004	M Beacon 5' ROX + 3' DABCYL 40 nmol
PB-MB216-020	M Beacon 5' ROX + 3' DABCYL 200 nmol
PB-MB216-100	M Beacon 5' ROX + 3' DABCYL 1000 nmol
PB-MB220-004	M Beacon 5' Rhodamine-6G + 3' DABCYL 40 nmol
PB-MB220-020	M Beacon 5' Rhodamine-6G + 3' DABCYL 200 nmol
PB-MB220-100	M Beacon 5' Rhodamine-6G + 3' DABCYL 1000 nmol
PB-MB230-004	M Beacon 5' Texas Red® + 3' DABCYL 40 nmol
PB-MB230-020	M Beacon 5' Texas Red® + 3' DABCYL 200 nmol
PB-MB230-100	M Beacon 5' Texas Red® + 3' DABCYL 1000 nmol
PB-MB240-004	M Beacon 5' Yakima Yellow® + 3' DABCYL 40 nmol
PB-MB240-020	M Beacon 5' Yakima Yellow® + 3' DABCYL 200 nmol
PB-MB240-100	M Beacon 5' Yakima Yellow® + 3' DABCYL 1000 nmol
PB-MB245-004	M Beacon 5' Dragonfly Orange™ + 3' DABCYL 40 nmol
PB-MB245-020	M Beacon 5' Dragonfly Orange™ + 3' DABCYL 200 nmol
PB-MB245-100	M Beacon 5' Dragonfly Orange™ + 3' DABCYL 1000 nmol
PB-MB250-004	M Beacon 5' Cy® 3 + 3' DABCYL 40 nmol
PB-MB250-020	M Beacon 5' Cy® 3 + 3' DABCYL 200 nmol
PB-MB250-100	M Beacon 5' Cy® 3 + 3' DABCYL 1000 nmol
PB-MB255-004	M Beacon 5' Cy5® + 3' DABCYL 40 nmol
PB-MB255-020	M Beacon 5' Cy5® + 3' DABCYL 200 nmol
PB-MB255-100	M Beacon 5' Cy5® + 3' DABCYL 1000 nmol
PB-MB260-004	M Beacon 5' Alexa® + 3' DABCYL 40 nmol
PB-MB260-020	M Beacon 5' Alexa® + 3' DABCYL 200 nmol
PB-MB260-100	M Beacon 5' Alexa® + 3' DABCYL 1000 nmol
PB-MB270-004	M Beacon 5' Cascade Blue® + 3' DABCYL 40 nmol
PB-MB270-020	M Beacon 5' Cascade Blue® + 3' DABCYL 200 nmol
PB-MB270-100	M Beacon 5' Cascade Blue® + 3' DABCYL 1000 nmol
PB-MB280-004	M Beacon 5' DYOMICS + 3' DABCYL 40 nmol
PB-MB280-020	M Beacon 5' DYOMICS + 3' DABCYL 200 nmol
PB-MB280-100	M Beacon 5' DYOMICS + 3' DABCYL 1000 nmol
PB-MB285-004	M Beacon 5' BODIPY® + 3' DABCYL 40 nmol
PB-MB285-020	M Beacon 5' BODIPY® + 3' DABCYL 200 nmol
PB-MB285-100	M Beacon 5' BODIPY® + 3' DABCYL 1000 nmol
PB-MB290-004	M Beacon 5' ATTO + 3' DABCYL 40 nmol

PB-MB290-020	M Beacon 5' ATTO + 3' DABCYL 200 nmol
PB-MB290-100	M Beacon 5' ATTO + 3' DABCYL 1000 nmol