

**MapMarker®400**  
**MW-0190-80ROX • MW-0190-80TMR • MW-0190-80FAM**

MapMarkers® are sets of high quality fluorescently labelled DNA fragments designed to provide consistent intensities and migration patterns for DNA sizing standards. They are compatible with most of fluorescent based separation instruments systems.

**Labelling**

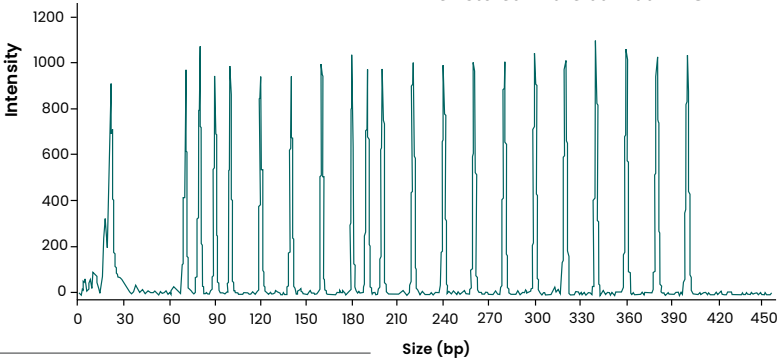
Either labelled with the x rhodamine (ROX), tetramethylrhodamine (TMR) or fluorescein (FAM) at 8 fmol/band/µl.

**Package content**

Reagent	References	Description
<b>MapMaker®400</b> 800 lanes, 400 µl 8 fmol/band/µl	MW-0190-80ROX MW-0190-80TMR MW-0190-80FAM	Labelled with: x rhodamine tetramethylrhodamine fluorescein
<b>Tracking Dye</b>	-	Tracking Dye contains 25 mg/mL blue dextran (average Molecular Weight of 2 000 000) dissolved in 25 mM EDTA

**Description**

MapMarker®400 contains 20 single stranded DNA fragments, ranging in size from 70 to 400 base pairs and labelled with a single fluorophore. The exact fragment sizes are: 70, 80, 90, 100, 120, 140, 160, 180, 190, 200, 220, 240, 260, 280, 300, 320, 340, 360, 380 and 400 base pairs.



MapMarker® is a registered trademark of BioVentures, Inc.  
Cy® and CyDye® are registered trademarks of Cytiva.

MapMarkers® are evaluated for resolution, intensity and background on capillary electrophoresis instrument using laser excitation. MapMarkers® are free of extraneous peaks over their entire sizing range and are free of “pull up” signal in other channels.

MapMarkers® may be used under either denaturing or non denaturing conditions.

MapMarkers® are compatible with many fluorescent based instrument systems, as well as a wide variety of fluorescent dyes including TAMRA, ROX, FAM, Cy5® equivalent and more.

**Recommendation for use**

The tracking dye could be utilized on any gel-based instrument. It is however not necessary when using a capillary or electrophoresis instrument.

**Sample Preparation**

For each sample combine 0.5 µL of MapMarker®, 0.5µL of Tracking Dye, 2.5 µL of dionized formamide and 1-2 µL of sample. A Master Mix can be prepared by adjusting the above component volumes to the number of samples to analyze. After the sample is added to the MapMarker/ Tracking Dye/Formamide mix, the solution should be briefly vortexed followed by brief centrifugation and denatured at 95°C for 5 minutes.

The solution is then cooled to 4°C.  
Load from 1 µL to 3 µL of this mixture per lane. The volume depends on the well capacity.

**Storage**

MapMarkers® are stable for a minimum of 18 months when stored in the dark at 4 °C.