EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00
• Fax: +32 4 372 75 00 • Toll-free: +800 666 00 123 • info@eurogentec.com
www.eurogentec.com



NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560 • Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com www.anaspec.com

MapMarker®500 MW-0191-800RANGE

Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

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

MapMarker® 500 is labelled with DY-632.

Package content

Reagent	References	Description
MapMaker®500 800 lanes, 400 μl	MW-0191- 80Orange	Labelled with: DY-632
Tracking Dye	-	Tracking Dye contains 25 mg/mL blue dextran (average Molecular Weight of 2 000 000) dissolved in 25 mM EDTA

Description

MapMarker®500 contains 16 single stranded DNA fragments, ranging in size from 35 to 500 base pairs and labelled with a single fluorophore. The exact fragment sizes are: 35,50,75,100,139,150,160,200,250, 300,340, 350,400,450,490 and 500 base pairs.

MapMarkers® are evaluated for resolution, intensity and background on denaturing polyacrylamide sequencing gels using laser excitation. MapMarkers® are free of extraneous peaks over their entire sizing range and are free of "pull up" signal in other channels.

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®, 0.5 μL 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 $1 \mu L$ to $3 \mu L$ of this mixture per lane. The volume depends on the well capacity.

Storage

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

For further information please contact our Customer Help Desk:

For Europe:

e-mail: info@eurogentec.com

Tel.: +32 4 372 76 65

USA:

e-mail: info.anaspec@eurogentec.com

Tel: +1 510 791 9560