

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
• Fax: +1 510-791-9572 • Toll-free: +1 800-452-5530 • service@anaspec.com
• www.anaspec.com

TECHNICAL DATA SHEET

BioTracker® 6X Tracking Dye EP-5010-10

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

Description

BioTracker® 6X is a Bromophenol Blue loading dye for agarose and non-denaturing polyacrylamide gel electrophoresis. BioTracker® is composed of 60 % sterile DI H₂O, 30 % glycerol, 10 % 10X TE, 0.25% BPB dye, and 0.75% SDS. Tracking dye contains 25 mg/mL blue dextran of an average molecular weight of 2,000,000 dissolved in 25mM EDTA.

Package contents

Reagent	Volume	Reference
BioTracker® 6X Tracking Dye	10 mL	EP-5010-10

Shipping conditions

Shipped at 4°C.

Storage conditions

Storage at 4°C is recommended.

Sample preparation

For each sample combine 0.5µL MapMarker®, 0.5 µL of Tracking Dye, 2.5µL deionized formamide and 1-2µL of Sample. A marker-dye master mix can be prepared by combining the above components in suitable quantities for the number of samples being analyzed.

After combining samples with marker/dye/formamide mix, the mixture should be briefly vortexed followed by brief centrifugation and denatured at 95°C for 5 min. followed by cooling to 4°C. Load from 1-3µL of this material per lane depending on well capacity.

Related Products

Reagent	Size	Reference
Mupid®-One Electrophoresis System	1 Apparatus	MU-0041
SmartLadder - 200 to 10000bp	1000 lanes	MW-1700-10
Molecular Biology Grade Agarose	500 g	EP-0010-05
MapMarker® 400 labelled with FAM	800 lanes	MW-0190-80FAM
MapMarker® 400 labelled with DY-632	800 lanes	MW-0190-80ORANGE
MapMarker® 400 labelled with ROX	800 lanes	MW-0190-80ROX
MapMarker® 400 labelled with TAMRA	800 lanes	MW-0190-80TMR
MapMarker® 500 labelled with DY-632	800 lanes	MW-0191-80 ORANGE
MapMarker® 1000 labelled with FAM	800 lanes	MW-0195-80FAM
MapMarker® 1000 labelled with ROX	800 lanes	MW-0195-80ROX
MapMarker® 1000 labelled with TAMRA	800 lanes	MW-0195-80TMR
MapMarker® 1000 labelled with DY-632	800 lanes	MW-0195-80 ORANGE

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
• Fax: +1 510-791-9572 • Toll-free: +1 800-452-5530 • service@anaspec.com
• www.anaspec.com

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65 • Toll-free: + 800 666 00 123

For USA:

E-mail: service@anaspec.com

Tel.: +1-510-791-9560 • Toll-free: +1-800-452-5530