EUROGENTEC HEADQUARTERS

LIEGE SCIENCE PARK • 4102 Seraing • Belgium • Tel.: +32 4 372 74 00 Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

EUROGENTEC NORTH AMERICA, INC.

11111 Flintkote Avenue • San Diego CA 92121-1222 USA Tel.:+1 858 793 2661 • Fax:+1 858 793 2666 • info.usa@eurogentec.com www.eurogentec.com



SmartLadder SF MW-1800-04

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

Descriptions

The SmartLadder SF is a ready-to-use molecular weight marker, especially designed for easy quantification and size determination of short DNA fragments.

Package content

Reagent	Volume	Description
SmartLadder SF 400 lanes	2 x 1 ml	2 tubes of 200 lanes each, ready to use

Shipping conditions

Shipped at room temperature. For long term storage, freeze upon arrival.

Storage

The SmartLadder SF is stable for 1 month at RT or at 4 °C for 6 months. For long-term storage keep at -20 °C. Avoid multiple freeze-thaw cycles.

Size range

The SmartLadder SF produces a pattern of 10 regularly spaced bands ranging from 100 to 1000 bp. All bands have

a different intensity to allow a quick and easy identification. The size of each band is an exact multiple of 100 bp.

Quantification

Using a standard loading volume of 5 μ l, each band corresponds to an exact quantity of DNA, from 20 to 100 ng.

Band size	ng/band
1000	100
900	90
800	80
700	70
600	60
500	50
400	40
300	30
200	20
100	20

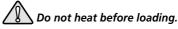
Loading Buffer composition

> Bromophenol blue	0.25 g/l
> Xylene cyanol	0.25 g/l
> Ficoll 400	25 g/l
> Sodium Azide	1 g/l
> Chloroform	1/1000

> TE (Tris 10mM, EDTA 1mM, pH 8)

Recommended Procedure

- 1. Vortex the ladder gently to ensure an homogenous solution.
- 2. Apply approximately 5 µl per 5 mm lane width.



Additional protocols:

T4 DNA Polymerase Labelling Protocol

- 1. Exonuclease Reaction (Degradation of DNA from both 3'-ends)
- > To a 1.5 ml microcentrifuge tube on ice, add the following:
 - 5X T4 DNA polymerase reaction buffer 4 μl (165 mM Tris acetate (pH 7.9), 330 mM sodium acetate, 50 mM magnesium acetate, 2.5 mM DTT, 500 μg/ml BSA)
 - SmartLadder
 T4 DNA polymerase
 40 units
 - Autoclaved water to 20 µl
- > Mix the tube thoroughly but not vigorously
- > Centrifuge briefly
- > Incubate 2 min at 37 °C (about 25 nucleotides/min are removed)
- > Cool reaction vial on ice
- 2. Resynthesis Reaction (Resynthesis of the degraded DNA strands)
- > Add into the reaction vial the following components:
 - Autoclaved water 8 μl
 - 5X T4 DNA polymerase reaction buffer 6 μl
 - dCTP (2 mM) 5 μl

5 µl

5 µl

- dGTP (2 mM)
- dTTP (2 mM)
- [α-³²P]dATP (3000 Ci/mmol; 10 mCi/ml) 1 μl

EUROGENTEC HEADQUARTERS

LIEGE SCIENCE PARK • 4102 Seraing • Belgium • Tel.: +32 4 372 74 00 Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

EUROGENTEC NORTH AMERICA, INC.

11111 Flintkote Avenue • San Diego CA 92121-1222 USA Tel.:+1 858 793 2661 • Fax:+1 858 793 2666 • info.usa@eurogentec.com www.eurogentec.com



- > Mix thoroughly
- > Centrifuge briefly
- > Incubate 2 min at 37 °C
- > Add 5 µl of 2 mM dATP
- > Incubate 2 min at 37 °C
- > Stop reaction by adding 2.5 µl of 0.5 M EDTA
- > Centrifuge for 10 s
- > The cpm incorporated is determined by adding 1 µl of reaction to 24 µl of 250 mM NaCl, 25 mM EDTA
- > Spot 5 μ l of dilution on a glass fiber filter
- > Place filter in 10 % (w/v) TCA + 1 % (w/v) pyrophosphate.
- > Wash filter 3 times with 5 % (w/v) TCA and then 2 times with ethanol
- > The filter is dried and then counted using an appropriate scintillant
- > Add 5 μ l 0.1 % (w/v) bromophenol blue, 0.1 mM EDTA, 50 % (v/v) glycerol to the sample
- > Load 1 x 10⁵ cpm in a lane.

5' DNA Terminus Labelling Protocol (Phosphate Exchange Reaction)

This reaction will yield specific activities of approximately 1-5 x 10⁵ cpm/pmol of ends.

T4 polynucleotide kinase 5'P–3'OH + [γ -³²P]ATP + ADP \longrightarrow 5'P³²–3'OH + ATP + ADP

- > Add the following components to a 0.5 ml microcentrifuge tube in the following order:
- Autoclaved water
 SmartLadder
 S μg
 SX exchange reaction buffer
 (250 mM imidazole (pH 6.4), 1.5 mM ADP, 60 mM MgCl₂, 75 mM, 2-mercaptoethanol)

– [γ- ³² P]ATP (10 μCi/μl)	3 µl
– T4 polynucleotide kinase (5 or 10 U/µl)	1 µl

- > Incubate the reaction mixture at 37 °C for 30 minutes. Increasing reaction times beyond 30 min will not increase labeling of the DNA.
- > Stop reaction by adding 1 μ l of 0.5 M EDTA.
- > Centrifuge for 10 s.
- > Determine radioactive incorporation as above.
- > Add 5 µl 0.1 % (w/v) bromophenol blue, 0.1 mM EDTA, 50 % (w/v) glycerol to the sample.
- > Load 1×10^5 cpm in a lane.

Related products

Reagent	Package size	Reference
Smart Ladder	1000 lanes	MW-1700-10
Agarose Molecular Biology Grade	100 g 500 g 1 Kg	EP-0010-01 EP-0010-05 EP-0010-10
Agarose AgaTabs	300 tablets	EP-0030-15
Mupid®-One Electrophoresis system	1	MU-0041

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com Tel: +32 4 372 76 65

For USA:

E-mail: info.usa@eurogentec.com Tel: +1 858 793 26 61