PRODUCT INFORMATION SHEET

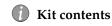
Pub. No. MAN0000840

Rev. A.0



Contents

Catalog Number Size 110488058 50 μq



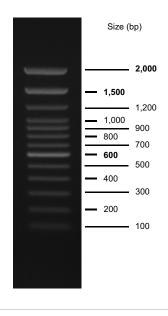


- Storage
- Product is shipped at ambient temperature.
- Store at room temperature or at 4°C for up to 6 months, or at -20°C for long term storage.



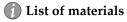
Product description

- The Invitrogen[™] TrackIt[™] 100 bp DNA Ladder is designed for sizing and quantification of double stranded DNA on 1% to 2% agarose gels.
- The 100 bp DNA Ladder consists of 13 individual chromatography-purified DNA fragments ranging in size from 100 bp to 2,000 bp.
- Three reference bands at 2,000 bp, 1,500 bp, and 600 bp are included for easy orientation.
- The ladder is supplied with 6X TrackIt[™] Cyan/ Orange Loading Buffer for sample DNA.





Required materials





Online resources

- Visit our product pages for additional information and protocols.
- Go online to view related DNA ladders and markers.
- For support, visit thermofisher.com/support.



Important guidelines

- Do not heat the TrackIt[™] 100 bp DNA Ladder before loading.
- Load the same volume of DNA sample and DNA ladder.
- For quantification, adjust the concentration of the sample to equalize it approximately with the amount of DNA in the nearest band of the ladder.
- For DNA bands visualization with GelRed[™] use gel staining after electrophoresis to avoid aberrant DNA migration.



Guidelines for agarose gel preparation

 Determine the required agarose concentration for your gel based on the size of DNA fragments to be separated.

Fragment size	Recommended agarose gel %				
	1X TAE	1X TBE			
800-10,000	0.8	0.7			
400-8,000	1.0	0.85			
300-7,000	1.2	1.0			

- Prepare agarose in a flask with 2-4 times the volume of the agarose solution.
- Exercise caution when handling microwaved agarose. The solution may become superheated and foam over when agitated.
- Refer to the product insert for UltraPure[™] Agarose for detailed instructions on agarose preparation.
- Guidelines for staining gels
- Troubleshooting
- Limited product warranty and disclaimer details

Prepare DNA ladders and samples for electrophoresis

Step		Action			
1	TILLILI.	Cast agarose gel	a. Prepare agarose solution (w/v) for the gel percentage appropriate for separating your DNA fragments.b. Microwave agarose solution.c. Cast agarose gel.		
2		Prepare DNA ladder	a. Thaw, mix and briefly centrifuge DNA ladder before use. b. Mix gently. c. Load the gel with 1 μL of DNA ladder per 1 mm of well width.		
3		Prepare samples	 a. Dilute your sample with 6X TrackIt™ Cyan/Orange Loading Buffer (Cat. no. 10482028): mix 1 volume of loading dye with 5 volumes of the DNA sample. b. Mix gently. c. Load DNA ladder on gel. 		
4		a. Add appropriate amount of UltraPure TAE or UltraPure TBE buffer to chamber.			
	Perform electrophoresis	b. Set appropriate voltage and perform electrophoresis of samples.			
		DNA size	Voltage	Buffer	
		<1 kb	5–10 V/cm	TBE	
		1-5 kb	4–10 V/cm	TAE or TBE	
		>5 kb	1–3 V/cm	TAE	
5		Stain agarose gel	 a. Incubate gel in staining buffer for 30 minutes. b. Visualize DNA ladder and samples. • Use UV transilluminator to detect DNA bands stained with ethidium bromide. • Use blue light transilluminator to detect DNA bands stained with SYBR™ stains. 		