E-Gel[™] 1 Kb Plus Express DNA Ladder

PRODUCT INFORMATION SHEET

Pub. No. MAN0017358

P	Contents	Catalog No. 10488091	Amount 100 applications	i Kit contents		
	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[™] E-Gel[™] 1 Kb Plus Express DNA Ladder is designed for sizing and quantification of double stranded DNA on 0.8–1% E-Gel[™] agarose gels.
- The E-Gel[™] 1 Kb Plus Express DNA Ladder consists of 9 individual chromatography-purified DNA fragments ranging in size from 100 bp to 5,000 bp.
- Reference bands at 500 bp and 1,500 bp are included for easy orientation.

Online

resources

• The ladder is supplied with 1X E-Gel[™] Sample Loading Buffer for sample DNA.



Rev. A.0

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

Required materials

- E-Gel[™] CloneWell[™] II or other E-Gel[™] agarose gel (See Choosing the right DNA ladder for your E-Gel[™] agarose gel)
- TE Buffer (Cat. No. AM9858)
- Ultrapure[™] DNase/RNase-Free Distillated Water (Cat. No. 10977023)

Important guidelines

- Do not heat the E-Gel[™] 1 Kb Plus 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.
- Dilute sample DNA in TE buffer to avoid degradation of DNA sample.
- Choosing the right DNA ladder for your E-Gel™ agarose gel
- Troubleshooting
- Limited product warranty and disclaimer details



For Research Use Only. Not for use in diagnostic procedures.

Prepare and load DNA ladders and samples

This protocol provides a brief description of how to use the DNA ladder with E-Gel[™] agarose gels. For detailed instructions on using specific types of E-Gel[™] agarose gels, go to thermofisher.com or contact Technical Support.

Step			Action		
1		Prepare DNA ladder	 a. Thaw, mix and briefly centrifuge DNA ladder before use. b. Prepare DNA ladder. For E-Gel[™] CloneWell[™] II Agarose Gels, mix and use 25 µL of the ladder without dilution. For E-Gel[™] EX Agarose Gels, mix 2 µL of DNA ladder with 18 µL of water. For E-Gel[™] Agarose Gels, mix and use the ladder without dilution. For E-Gel[™] 48 Agarose Gels, mix 2 µL of DNA ladder with 13 µL of water. 		
2 Prepare sample		Prepare samples	 a. Dilute your sample 2- to 10-fold with TE Buffer (Cat. No. AM9858), 1X E-Gel[™] Sample Loading Buffer (Cat No. 10482055), or water. b. Mix gently. 		
3		Load samples and DNA ladders	 a. Load DNA ladders and DNA samples into the appropriate wells of the E-Gel[™] agarose gel. Add 25 µL for E-Gel[™] CloneWell[™] II Agarose Gels. Note: All wells of E-Gel[™] Clonewell[™] II Agarose Gels must be pre-filled with 50 µL of water. Add 20 µL for E-Gel[™] and E-Gel[™] EX Agarose Gels. Add 15 µL for E-Gel[™] 48 Agarose Gels. b. Add water to any empty wells, so that all wells contain an equal volume of liquid. c. Proceed to Perform electrophoresis of E-Gel[™] agarose gel. 		

Thermo Fisher

SCIENTIFIC

Perform electrophoresis of E-Gel[™] agarose gel

This protocol provides a brief description of how to perform electrophoresis with E-Gel[™] agarose gels. For detailed instructions on using specific types of E-Gel[™] agarose gels, go to thermofisher.com or contact Technical Support.

Step			Action			
		Perform electrophoresis	a. Choose the appropriate E-Gel [™] run protocol for your gel type based on the electrophoresis device being used.			
				Gel type	Program	Recommended run time
4				E-Gel™ Power Snap Electrophoresis Device (Cat. No. G8100)		
				E-Gel [™] Clonewell [™] II Agarose Gel (0.8%)	CloneWell 0.8%	12 min (40 min max)
				E-Gel™ EX Agarose Gel (1%, 2%)	E-Gel EX 1-2%	10 min (20 min max)
				E-Gel [™] Agarose Gel (0.8%, 1.2%, 2%)	E-Gel 0.8-2%	26 min (40 min max)
				E-Gel [™] Double Comb Agarose Gel (1%, 2%)	E-Gel Double Comb	13 min (20 min max)
				E-Gel [™] E-Base [™] Device		
				E-Gel [™] 48 Agarose Gel (1%, 2%)	EG	20 min
			b. Run the program to start electrophoresis.			
	Contraction of the second seco	Visualize agarose gel	Visualize DNA ladder and samples.			
5			 Use the E-Gel[™] Power Snap Camera (Cat. No. G8200), E-Gel[™] Imager (Cat. No. 466612), or other blue light imager to detect DNA bands stained with SYBR[™] stains. 			
				 UV transilluminator to detect DNA base 	ands stained with ethidium bromide.	

Thermo Fisher

SCIENTIFIC