Fast Digestion of DNA

This protocol is for the Fast Digestion of DNA

1. Prepare the reaction mixture at room temperature in the order indicated:

	Volume		
Component	Plasmid DNA	Unpurified PCR product	Genomic DNA
Water, nuclease-free*	15 µl	17 µl	30 µl
10X FastDigest® buffer or 10X FastDigest® Green buffer	2 µl	2 µl**	5 µl
DNA*	2 μl (up to 1 μg)	10 µl (~0.2 µg)	10 µl (5 µg)
FastDigest® enzyme	1 µl	1 µl	5 µl
Total volume	20 µl	30 µl	50 µ l

- 2. Mix gently and spin down.
- 3. Incubate at 37°C in a heat block or water thermostat for 5 min.***
- 4. Inactivate the enzyme (optional).***

Note

- * The volume of water should be corrected to keep the indicated total reaction volume. The volume of DNA can be scaled up to 10 μ l or down to 0.5 μ l depending on the DNA concentration.
- ** Only 2 μl of 10X FastDigest® buffer is required for unpurified PCR product in a 30 μl reaction volume.
- *** See the Certificate of Analysis for enzyme and substrate specific incubation time and enzyme inactivation conditions.

thermoscientific.com

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

North America Technical Services: techservice.genomics@thermofisher.com CustomerService.genomics@ thermofisher.com Tel 800 235 9880 Fax 800 292 6088 Europe and Asia Technical Services: techservice.emea.genomics@ thermofisher.com Customer Services: customer service.emea.genomics@ thermofisher.com

