Double and Multiple Digestion of DNA

This protocol is for the Double and Multiple Digestion of DNA

FastDigest® enzymes allow simultaneous digestion of DNA with two or more enzymes in one digestion reaction.

- Use 1 µl of each enzyme and scale up the reaction conditions appropriately.
- The combined volume of all added enzymes should not exceed ¹/10 of the total reaction volume.
- If enzymes require different incubation temperature, perform sequential DNA cleavage: complete the first digestion reaction at the lower temperature, add the second enzyme and increase the digestion temperature for the second enzyme cleavage.

© 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 CustomerServices: 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

