Scaling up DNA Digestion Reaction

This protocol is for the Scaling up DNA Digestion Reaction

DNA	1 µg	2 µg	3 µg	4 µg	5 µg
FastDigest® enzyme	1 µl	2 µl	3 µl	4 µl	5 µl
10X FastDigest® buffer or 10X FastDigest® Green buffer	2 µl	2 µl	3 µl	4 µl	5 µl
Total volume	20 µl	20 µl	30 µl	40 µl	50 µl

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: customerservice.emea.genomics@ thermofisher.com

