

# Reaction Set-up for Digestion of Multiple DNA Samples

This protocol is for the Reaction Set-up for Digestion of Multiple DNA Samples

1. Pipette 2 µl of each DNA sample\* into labeled tubes.
2. Prepare a master mix for n+1 samples.

Example of master mix (for 10 samples of plasmid DNA):

<b>Water, nuclease-free</b>	$(10 + 1) \times 15 \mu\text{l} = 165 \mu\text{l}$
<b>10X FastDigest® buffer or 10X FastDigest® Green buffer</b>	$(10 + 1) \times 2 \mu\text{l} = 22 \mu\text{l}$
<b>FastDigest® enzyme</b>	$(10 + 1) \times 1 \mu\text{l} = 11 \mu\text{l}$

3. Add 18 µl of master mix into tubes containing DNA

Note

\* The volume of DNA can be scaled up to 10 µl or down to 0.5 µl depending on the DNA concentration. The volume of water and master mix should be corrected to keep the indicated total reaction volume.

[thermoscientific.com](http://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  
**Customer Services:**  
 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

**Thermo**  
 S C I E N T I F I C

Part of Thermo Fisher Scientific