

# Removal of Template DNA after *in vitro* Transcription

This protocol is for the Removal of Template DNA after *in vitro* Transcription

1. Add 2 u of DNase I, RNase-free per 1 µg of template DNA directly to a transcription reaction mixture.  
In some cases, the amount of enzyme should be determined empirically.
2. Incubate at 37°C for 15 minutes.
3. Inactivate DNase I by phenol/chloroform extraction

[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**  
SCIENTIFIC

Part of Thermo Fisher Scientific