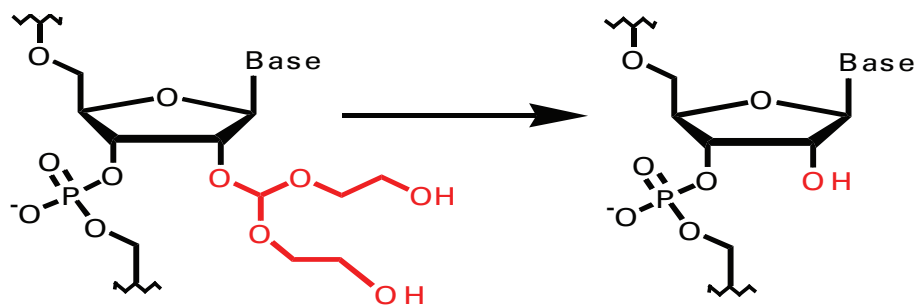


Thermo Scientific Deprotection 2'- ACE Protected RNA

This protocol is for the Thermo Scientific Deprotection 2'- ACE Protected RNA.

1. Centrifuge tubes briefly.
2. Add 400 μ L of 2'- Deprotection buffer to each tube of RNA.
3. Completely dissolve RNA pellet by pipetting up and down.
4. Vortex for 10 seconds and centrifuge for 10 seconds.
5. Incubate at 60 ° C for 30 minutes.
 - Incubate at 60 ° C for 2 hours for oligos with biotin modifications or homopolymer stretches of rA longer than 10 bases.
6. Lyophilize or SpeedVac to dryness before use.
7. The dry pellet can be stored at -20 ° C until use or resuspended in an appropriately buffered RNase-free solution.
8. If appropriate for your application you may perform desalting of the RNA oligo by Ethanol Precipitation (<http://www.thermoscientific.com/uploadedFiles/Home/Resources/Protocols/ethanol-precipitation-of-rna-oligonucleotide-protocol.pdf>) or desalting columns (<http://www.thermoscientific.com/uploadedFiles/Home/Resources/Protocols/reverse-phase-protocol.pdf>).

2'-Deprotection buffer = 100 mM acetic acid, adjusted to pH 3.8 with TEMED.



thermoscientific.com/onebio

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Europe

Customer Service
cs.molbio.eu@thermofisher.com

Technical Support
ts.molbio.eu@thermofisher.com

Tel 00800 222 00 888
Fax 00800 222 00 889

United States

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 235 9880
Fax 800 292 6088

Canada

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 340 9026
Fax 800 472 8322

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific