

Phenol/Chloroform Extraction and Ethanol Precipitation

This protocol is for the Phenol/Chloroform Extraction and Ethanol Precipitation

1. Mix your sample with 1 volume of Tris-saturated phenol and 1 volume of chloroform. Centrifuge at 10,000 rpm for 5 min at room temperature.
2. Transfer the upper aqueous phase to a fresh tube. Add an equal volume of chloroform and mix. Centrifuge at 10,000 rpm for 5 min at room temperature. Repeat.
3. Transfer the upper aqueous phase to a fresh tube. Add 1/10 the volume of 3 M Sodium Acetate Solution or 2 M sodium chloride.
4. Add 2.5 volumes of ethanol or an equal volume of isopropanol to precipitate DNA.
5. Incubate the mixture for 30 min at -20°C.
6. Centrifuge for 10 min at 10,000 rpm. Then discard the supernatant and rinse the pellet with 70% cold ethanol.
7. Air-dry the pellet. Dissolve in Water, nuclease-free or TE buffer.

Note

Use Glycogen to maximize the yield of DNA during precipitation.

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