

Recommendations for Accurate Gel Quantification

This protocol is for the Recommendations for Accurate Gel Quantification

- Always use the same DNA loading dye (supplied with the DNA ladder/marker) for both the sample DNA and the ladder/marker DNA.
- Always compare the sample band with the ladder band of the closest size.
- If possible, adjust the concentration of the sample to approximately equalize it with the amount of DNA in the nearest band.
- dNTPs, oligonucleotides, genomic DNA, RNA, NTPs or buffer components can interfere with spectrophotometrical measurements and lead to inaccurate quantification of sample DNA. In these cases, it is best to rely on gel quantification data.
- For the most accurate quantification, use video-densitometry analysis.

Reference

1. Sambrook, J., et al., *Molecular Cloning. A Laboratory Manual*, Cold Spring Harbor Laboratory, Cold Spring Harbor, N.Y., 12.89, 5.42, 2001.

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