



# **Product Information**

## RNase-Free Calf Thymus DNA, 1 mg/mL

Catalog Number: 31065

Unit Size: 1 mL

### Storage and Handling

Store at 4°C. Product is stable for at least 12 months from date of receipt when stored as recommended.

### **Product Description**

Many commercial preparations of DNA contain high levels of residual RNase activity, even after re-purification. Preparing DNA solutions in DEPC-treated water will not inactivate RNase already present in the DNA. Similarly, commercial RNase inhibitor additives are usually poorly effective, expensive, and interfere in downstream applications.

Biotium has developed a unique proprietary method that irreversibly inactivates RNases in purified DNA. The resulting DNA solution is free of excess protein or other chemicals that may interfere with downstream applications. Our RNase-free Calf Thymus DNA solution undergoes stringent quality control testing to confirm the product is 100% free of RNase activity.

RNase-Free DNA is useful in applications where RNase contamination needs to be avoided. It is suitable for preparing RNase-free dsDNA standards for use with DNA quantitation assays such as the AccuClear® Ultra High Sensitivity DNA Quantitation Solution (31027), the AccuBlue® Broad Range dsDNA Quantitation Solution (31009), or the AccuBlue® NextGen dsDNA Quantitation Solution (31061), if RNase-free reagents are desired. See the Related Products table and our website for more information about AccuClear® and AccuBlue® dsDNA Quantitation Kits.

Biotium also offers the AccuBlue® Broad Range RNA Quantitation Kit (31073) for sensitive and accurate quantitation of RNA samples in fluorescent plate reader or Qubit® instruments.

#### **Related Products**

Catalog number	Product
31073	AccuBlue® Broad Range RNA Quantitation Kit
31027	AccuClear® Ultra High Sensitivity dsDNA Quantitation Solution
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit with 7 DNA Standards
31009	AccuBlue® Broad Range dsDNA Quantitation Solution
31007	AccuBlue® Broad Range dsDNA Quantitation Kit with 9 DNA Standards
31061	AccuBlue® NextGen dsDNA Quantitation Solution
31060	AccuBlue® NextGen dsDNA Quantitation Kit
41024	Water, Ultrapure Molecular Biology Grade
41028	Agarose LE, Ultra-Pure Molecular Biology Grade
41029	GelRed® Agarose LE
41003	GelRed® Nucleic Acid Gel Stain, 10,000X in Water
41005	GelGreen® Nucleic Acid Gel Stain, 10,000X in Water
31041	Forget-Me-Not™ EvaGreen® qPCR Master Mix (2-Color Tracking)
31045	Forget-Me-Not™ EvaGreen® qPCR Master Mix (Low Rox)
31043	Forget-Me-Not™ Universal Probe Master Mix
CD202	DNAstorm™ Kit for Isolation of DNA from FFPE Tissues Samples
CD201	RNAstorm™ Kit for Isolation of RNA from FFPE Tissues Samples
CD504	RNAstorm™ RNA Isolation Kit

Please visit our website at <a href="www.biotium.com">www.biotium.com</a> for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, DNA and protein gel stains, DNA and RNA quantitation kits, fluorescent probes, and kits for cell biology research.

Qubit is a registered trademark of Thermo Fisher Scientific.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.