



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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

Double-stranded RNase-free Calf Thymus DNA prepared using a proprietary method to irreversibly inactivate RNase activity. Useful in applications where RNase contamination needs to be avoided.



### Product attributes

### Product Description

RNase-free Calf Thymus is prepared using a proprietary method to irreversibly inactivate RNase activity and is guaranteed to be free of RNase activity. It is useful in applications where RNase contamination needs to be avoided.

- Calf thymus DNA in 1X Tris-EDTA
- 1 mg/mL concentration
- 1 mL volume
- Guaranteed RNase-Free
- Store at 4°C

During the isolation of purified DNA, RNase treatment is performed to eliminate contaminating RNA. Because RNase activity is robust and resistant to denaturation, 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 RNases 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.

This datasheet was generated on January 1, 2026 at 08:24:26 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/rnase-free-calf-thymus-dna-1-mgml/>