

**PRODUCT INFORMATION**

**0.5 M EDTA, pH 8.0,**

molecular biology grade

#\_ —

Lot: —

**Store at 4°C**

[www.thermoscientific.com/onebio](http://www.thermoscientific.com/onebio)

**Description**

0.5 M EDTA, pH 8.0, is 0.22 µm membrane-filtered molecular biology grade solution of EDTA in water.

**Application**

Suitable for biochemistry or molecular biology applications that require a chelator for divalent cations.

**CERTIFICATE OF ANALYSIS**

**Nicking Activity Assay**

Incubation of 1 µg of pUC19 DNA with 1 µl of 0.5 M EDTA, at 37°C for 17 hours and analysis on gel generated no detectable conversion of covalently closed circular DNA to nicked DNA.

**Labeled Oligonucleotide (LO) Assay**

No detectable degradation of single-stranded and double-stranded 5'-[<sup>33</sup>P]-labeled oligonucleotides was observed after incubation with 0.5 µl of EDTA at 37°C for 16 hours.

**Ribonuclease Assay**

No detectable RNA degradation was observed after incubation of 0.08 µg 2000 base RNA transcript with 0.5 µl of EDTA at 37°C for 4 hours.

**Quality authorized by:**

 Jurgita Zilinskiene

**PRODUCT USE LIMITATION**

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to [www.thermoscientific.com/onebio](http://www.thermoscientific.com/onebio) for Material Safety Data Sheet of the product.

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