

10X TAE Ultra Pure MB Grade

Store at room temperature.

Catalog #:	AM9869
Volume:	1 L
Appearance:	Colorless liquid
Product Description:	A pH-adjusted blend of ultra pure grade Tris-Acetate buffer and EDTA which, when diluted to a 1X working concentration, contains 40 mM Tris-Acetate and 1 mM Na ₂ EDTA.
Molecular Weight:	Tris component 121.14 EDTA component 372.3
Storage Conditions:	Store at room temperature.

USER INFORMATION

General Information:	Tris-Acetate-EDTA (TAE) buffers are used primarily in gel electrophoresis of nucleic acids. Each solution of 10X TAE is ready to use upon dilution to the desired concentration. All components are ultra pure molecular biology grade.
-----------------------------	--

QUALITY CONTROL

Nonspecific Endonuclease Activity:	Meets or exceeds specification when a sample is incubated for 14–16 hr with 300 ng supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.
Exonuclease Activity:	Meets or exceeds specification when a sample is incubated for 14–16 hr with 40 ng labeled <i>Sau3A</i> fragments of pUC19 and analyzed by PAGE.
RNase Activity:	Meets or exceeds specification when a sample is incubated for 14–16 hr with 25 ng labeled RNA and analyzed by PAGE.
Functional Testing:	Molecular weight markers and plasmid DNA are run on a 1% Agarose–LE gel using 1X TAE. All bands are sharp and intact.

OTHER INFORMATION

Material Safety Data Sheets:	Material Safety Data Sheets (MSDSs) can be printed or downloaded from product-specific links on our website at the following address: www.ambion.com/techlib/msds . Alternatively, e-mail your request to MSDS_Inquiry_CCRM@appliedbiosystems.com . Specify the catalog or part number(s) of the product(s), and we will e-mail the associated MSDSs unless you specify a preference for fax delivery. For customers without access to the internet or fax, our technical service department can fulfill MSDS requests placed by telephone or postal mail. (Requests for postal delivery require 1–2 weeks for processing.)
-------------------------------------	---

Warranty and Liability:	<p><i>For research use only. Not for use in diagnostic procedures.</i></p> <p>Applied Biosystems is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at www.ambion.com/info/warranty in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Applied Biosystems if you have any questions about our warranties or would like information about post-warranty support.</p> <p>Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.</p>
--------------------------------	--

Trademarks, Patents, and Licensing:	Applied Biosystems, AB (Design), and Ambion are registered trademarks of Applied Biosystems Corporation or its subsidiaries in the US and/or certain other countries. All other trademarks are the sole property of their respective owners.
--	--

© 2008 Ambion, Inc. All rights reserved. 4386656A