

# flashPAGE™ Pre-cast Gels (Type A)

Store at 4°C.

**Catalog # (P/N):** AM10010

**Product Description:** Ambion® flashPAGE Pre-cast Gels (Type A) are a proprietary formulation of polyacrylamide gel columns in storage buffer. They are designed for use with the flashPAGE Fractionator (P/N AM13100) and flashPAGE Buffer Kit Type A (P/N AM9015) for gel purification of nucleic acids <40 nt.

**Storage Conditions:** Store at 4°C. flashPAGE™ Pre-cast Gels (Type A) are guaranteed until the expiration date on the package, if stored properly. **Do not freeze flashPAGE Pre-cast Gels.**

## USER INFORMATION

### Applications:

#### Inspect the Pre-Cast Gel

IMPORTANT: Inspect each flashPAGE Pre-cast Gel before use to make sure that the bottom of the gel inside the cartridge is flush with the bottom of the column. If the gel has shifted during shipping, gently push the gel down so that the bottom of the gel aligns with the bottom of the column. We suggest using the wide end of a nuclease-free 10 µL pipette tip to push the gel, but any blunt nuclease-free implement can be used.

#### Running flashPAGE gels

Instructions for running flashPAGE Pre-cast Gels are provided on the flashPAGE Fractionator Protocol card which is supplied with the device and is available at the following web address:  
[www.ambion.com/catalog/CatNum.php?13100](http://www.ambion.com/catalog/CatNum.php?13100).

## QUALITY CONTROL

**Nonspecific Endonuclease Activity:** A sample is incubated for 14–16 hr with supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.

**Exonuclease Activity:** A sample is incubated for 14–16 hr with labeled double-stranded DNA, followed by PAGE analysis.

**RNase Activity:** A sample is incubated for 14–16 hr with labeled RNA, followed by PAGE analysis.

**Functional Testing:** Each lot of flashPAGE Pre-cast Gels is functionally tested using the Ambion flashPAGE Buffer Kit (Type A), the flashPAGE Fractionator, and Decade™ Markers (P/N AM7778). Each lot of flashPAGE Pre-cast Gels is randomly sampled and tested for RNase and DNase contamination.

## 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](http://www.ambion.com/techlib/msds). Alternatively, e-mail your request to [MSDS\\_Inquiry\\_CCRM@appliedbiosystems.com](mailto: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:** *For research use only. Not for use in diagnostic procedures.*

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](http://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.

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

Applied Biosystems disclaims all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall Applied Biosystems be liable, whether in contract, tort, warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

### Trademarks, Patents, and Licensing:

Applied Biosystems, AB (Design), and Ambion are registered trademarks, and flashPAGE and Decade are trademarks of Applied Biosystems 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. 4386450A