

# Adeno-X™ Maxi Purification Kit Protocol-at-a-Glance

(PT3680-2)

Please read the *User Manual* before using this Protocol-at-a-Glance. This abbreviated protocol is provided for your convenience, but is not intended for first-time users.

**Purification Protocol** (see other side for a graphic overview of the protocol)

1. Equilibrate the Filter Assembly with 5 ml 1X Equilibration Buffer.  
**Note:** The flow rate for this and all subsequent steps should be **3 ml/min (~1 drop/sec)**.
2. Pass the diluted and clarified lysate through the Purification Filter to allow the virus to bind.
3. Wash the filter with 20 ml 1X Wash Buffer.
4. Remove the filter from the Assembly.
5. To elute the adenovirus, attach the filter to a new syringe containing 3 ml 1X Elution Buffer. Push 1 ml of Elution Buffer through the filter into a sterile 15 ml conical tube.
6. Incubate the filter at room temperature for 5 min.
7. Push the remaining elution buffer through to elute the rest of the adenovirus. Use residual air in the syringe to push any remaining virus through the filter.
8. Determine the adenoviral titer.  
**Note:** We recommend using the Adeno-X Rapid Titer Kit (Cat. No. 632250) or the Adeno-X qPCR Titration Kit (Cat. No. 632252).
9. Use immediately, or aliquot and store the adenovirus at  $-70^{\circ}\text{C}$ .

**Note:**

For improved long-term stability, and proper tonicity for *in vivo* applications, we recommend a buffer exchange of the eluted adenovirus into 1X Formulation Buffer.

1X Formulation Buffer:

2.5% glycerol (w/v), 25 mM NaCl, and 20 mM Tris-HCl, pH 8.0 (GTS buffer; Hoganson, *et al.*, 2002)



Clontech

United States/Canada  
800.662.2566

Asia Pacific  
+1.650.919.7300

Europe  
+33.(0)1.3904.6880

Japan  
+81.(0)77.543.6116

Clontech Laboratories, Inc.  
A Takara Bio Company  
1290 Terra Bella Ave.  
Mountain View, CA 94043  
Technical Support (US)  
E-mail: [tech@clontech.com](mailto:tech@clontech.com)  
[www.clontech.com](http://www.clontech.com)

(080912; published 09 August 2012)

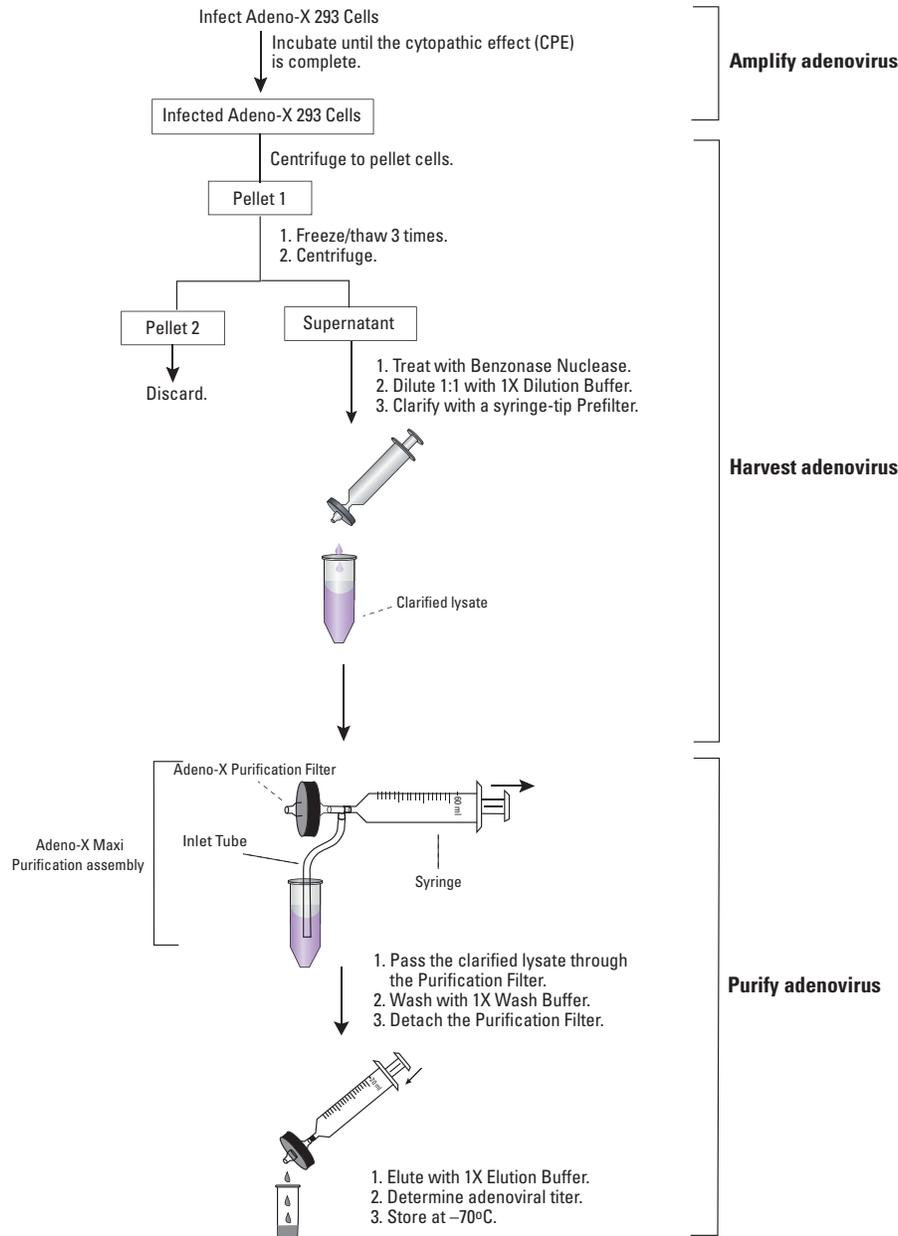


Figure 2. Overview of the Adeno-X Maxi Purification Protocol.

#### Notice to Purchaser

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without written approval of Clontech Laboratories, Inc.

Your use of these products and technologies is subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech, the Clontech logo and Adeno-X are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2012 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.