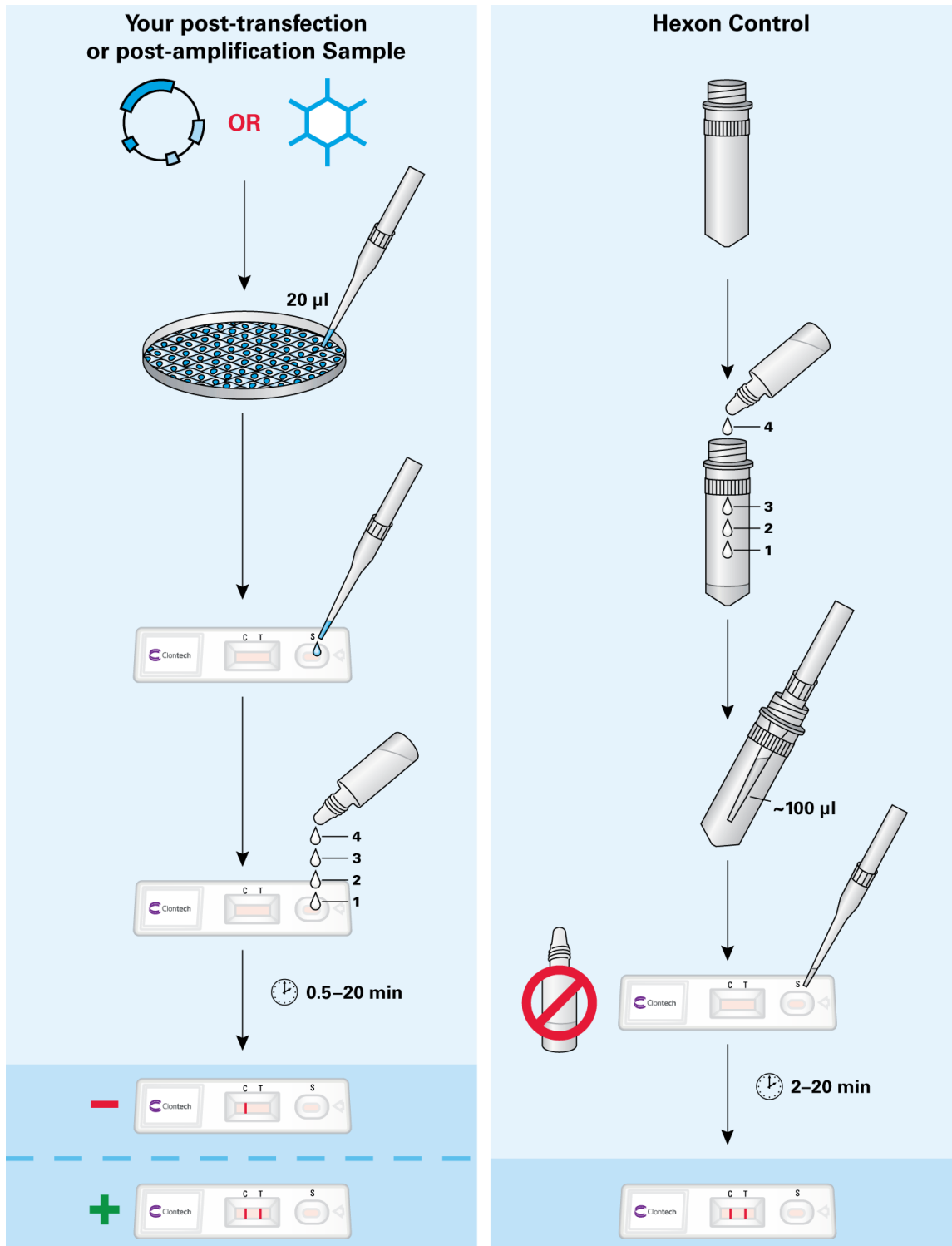


I. Introduction

Adeno-X GoStix (Cat. No. 632270) are designed to instantly confirm the presence of adenovirus in cell culture supernatants or lysates. The test, which detects adenoviral hexon proteins, requires just 20 µl of adenoviral supernatant to determine if infectious virus was produced posttransfection, or to select the best time to harvest your virus following amplification. The included Hexon Control provides confirmation of GoStix test function.



II. Your Sample Test—After Transfection of Adenoviral DNA

1. Take 20 µl of your adenoviral supernatant and apply it to the sample well (S) of the GoStix cassette.
2. Add 4 drops of Chase Buffer 2 to the sample well and wait for the bands to develop.
3. A control band will always appear (C), and a test band (T) will start to appear within 2–20 min. The test band will reach maximum intensity at 20 min if your sample contains sufficient adenovirus. **DO NOT** exceed 20 min. A positive test band indicates your sample contains $> 2.0 \times 10^5$ IFU/ml.

III. Your Sample Test—After Amplification of Adenovirus

1. Take 20 µl of your adenoviral supernatant (or cell lysate) and apply it to the sample well (S) of the GoStix cassette.
2. Add 4 drops of Chase Buffer 2 to the sample well and wait for the bands to develop.
3. A control band will always appear, and a test band will start to appear within 30 sec–20 min. The test band will reach maximum intensity at 5 min if your sample contains sufficient adenovirus. **DO NOT** exceed 20 min, since the likelihood of a false positive will increase.

NOTE: If the viral titer is expected to exceed 10^8 IFU/ml, test a series of sample dilutions, e.g., 1/10, 1/100, 1/1,000, on GoStix cassette.

IV. Hexon Control Test

1. Resuspend the lyophilized Hexon Control in 4 drops of Chase Buffer 2 and pipette up and down to mix.
2. Apply 100 µl of the Hexon Control to the sample well (S) of the GoStix cassette.
3. 2 bands (control and test) will appear in 2–20 min. **DO NOT** exceed 20 min.

Contact Us For Assistance	
Customer Service/Ordering	Technical Support
Telephone: 800.662.2566 (toll-free)	Telephone: 800.662.2566 (toll-free)
Fax: 800.424.1350 (toll-free)	Fax: 800.424.1350 (toll-free)
Web: www.clontech.com	Web: www.clontech.com
E-mail: orders@clontech.com	E-mail: tech@clontech.com

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 prior written approval of Clontech Laboratories, Inc.

Your use of this product is also 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, Adeno-X, and GoStix 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.