

# illustra GenomiPhi V2 DNA Amplification Kit

## Product Specification Sheet

### Introduction

#### Product codes

25660030

25660031

25660032

#### Important

Read these instructions carefully before using the products.

#### Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

#### Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

# Description

illustra™ GenomiPhi™ V2 DNA Amplification Kit is part of the Phi29 DNA polymerase family of products from Cytiva. It contains all of the components necessary for mini-scale whole genome amplification by isothermal multiple strand displacement. Amplification is highly uniform over the entire genome ensuring that locus representation remains extremely close to the original DNA sample. Furthermore, amplification is carried out with very high fidelity due to Phi29 DNA polymerase proofreading activity.

**Note:**      *DO NOT allow reconstituted reagents to warm above 4°C prior to amplification.*

# Short Protocol

**Note:**      *The Short Protocol is appropriate for researchers who have optimized their GenomiPhi reactions. First-time users must review the Full Protocol. The full protocol can be obtained by visiting [cytiva.com](https://www.cytiva.com), by e-mail and fax, or by contacting your local Cytiva Technical Support Group.*

## Heat denaturation of template in sample buffer

| Step | Action                                                                |
|------|-----------------------------------------------------------------------|
| 1    | Mix 1 µL of template DNA (at least 10 ng) with 9 µL of sample buffer. |
| 2    | Heat to 95°C for 3 minutes.                                           |
| 3    | Cool to 4°C on ice.                                                   |
| 4    | Proceed with the next part of the protocol.                           |

## Preparation of amplification reaction

| Step | Action                                                                                           |
|------|--------------------------------------------------------------------------------------------------|
| 1    | For each amplification reaction, combine 9 µL of Reaction Buffer with 1 µL of enzyme mix on ice. |
| 2    | Add this to the cooled sample.                                                                   |
| 3    | Proceed with the next part of the protocol.                                                      |

## Incubation at 30°C

| Step | Action                                                                            |
|------|-----------------------------------------------------------------------------------|
| 1    | Incubate the sample at 30°C for 90 minutes (2 hours suggested for crude lysates). |
| 2    | Proceed with the next part of the protocol.                                       |

## Post-amplification heat inactivation

| Step | Action                                  |
|------|-----------------------------------------|
| 1    | Heat the sample to 65°C for 10 minutes. |
| 2    | Cool to 4°C.                            |

## Optional purification of amplified products

| Step | Action                                  |
|------|-----------------------------------------|
| 1    | Purify amplified material if necessary. |

**Note:** *DO NOT allow reconstituted reagents to warm above 4°C prior to amplification.*



## cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

illustra and GenomiPhi are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

For use only as licensed by Qiagen GmbH. The Phi 29 DNA polymerase may not be re-sold or used except in conjunction with the other components of this kit. See US patent number 6,323,009, and equivalent patents and patent applications in other countries.

All other third-party trademarks are the property of their respective owners.

© 2020 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

29046315 AC V:3 12/2020