

Human Genomic DNA

Cat. No. 11 691 112 001

100 µg (500 µl)

 **Version 09**

Content version: April 2016

Store at +2 to +8°C

1. Product Overview

Concentration

0.2 mg/ml in 10 mM Tris-HCl, 1 mM EDTA, pH 8.0.

Source

High molecular weight genomic DNA is isolated from human blood (buffy coat).

Application

This DNA is suitable for Southern hybridization analysis, genomic library construction and amplification of large DNA targets.

Storage and Stability

The unopened reagent is stable at +2 to +8°C until the expiration date printed on the label.

- Ⓢ Do not freeze. Due to logistic reasons this product has to be delivered on dry ice. This one time freezing does not affect the quality of the DNA as checked by our Quality Control.
- Ⓢ After delivery store at +2 to +8°C

2. Quality Control

Purified genomic DNA was found to have a molecular weight greater than 50 kb as determined by visualisation on a 0.5% ethidium bromide stained agarose gel.

PCR amplification products up to 23 kb could be obtained using the Roche Diagnostics Expand 20kb^{PLUS} PCR System*.

- Ⓢ The blood of the donors for this preparation was tested for Hbs antigen and for the presence of anti-bodies to HIV-1, HIV-2, HCV and found to be negative

Changes to Previous Version

- Editorial Changes

Trademarks

EXPAND is a trademark of Roche.

All other brands or product names are the trademarks of their respective holders.

Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

* available from Roche Diagnostics

Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit sigma-aldrich.com, and select your home country. Country-specific contact information will be displayed.