

## PRODUCT INFORMATION

# pUC18 DNA

#SD0051 50 µg

Lot: —

Concentration: 0.5 µg/µL

Store at -20°C



## Description

Double-stranded closed circular high copy plasmid, 2686 base pairs with a molecular weight of  $1.74 \times 10^6$  Da. Isolated from *E.coli* (*dam*<sup>+</sup>, *dcm*<sup>+</sup>). GeneBank/EMBL accession number L09136. For Map and features see [www.thermoscientific.com/reviewer](http://www.thermoscientific.com/reviewer).

## Features

- Purified by chromatography using proprietary patented technology.
- More than 90% in the supercoiled form.

## Applications

- Cloning.
- Sequencing of insert DNA.
- Preparation of DNA molecular weight standards.

## Storage Buffer

10 mM Tris-HCl (pH 7.6) and 1 mM EDTA.

## CERTIFICATE OF ANALYSIS

DNA concentration is confirmed spectrophotometrically. Identity and homogeneity of DNA is confirmed by digestion by HindIII+NdeI and HindIII and analysis of DNA fragmentation patterns.

## Quality authorized by:

 Jurgita Zilinskiene

## PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively for research purposes and in vitro use only. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to [www.thermoscientific.com/onebio](http://www.thermoscientific.com/onebio) for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.