

PRODUCT INFORMATION

pUC57 DNA

#SD0171 **50 µg**

Lot: —

Concentration: 0.5 µg/µL

Store at -20°C

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 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.

www.thermoscientific.com/onebio

Description

Double-stranded closed circular high copy plasmid, 2710 base pairs with a molecular weight of 1.76×10^6 Da. Isolated from *E.coli* (*dam*⁺, *dcm*⁺).

GeneBank/EMBL accession number Y14837.

For map and features see

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.
- Generation of nested deletions with ExoIII.

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 NdeI+HindIII and HindIII and analysis of DNA fragmentation patterns.

Quality authorized by:

 Jurgita Zilinskiene