

DNA Coating Solution

CATALOG NUMBER: AKR-5183

QUANTITY: 100 mL

Background

It is known that DNA has a low binding affinity for unmodified plastic surfaces, so positively charged groups such as amino groups are often used to precoat plates. However, these precoating procedures are costly and time-consuming.

Cell Biolabs' DNA Coating Solution is a quick, simple and cost-effective method for immobilizing DNA and Oligonucleotides onto unmodified microplate wells.

Storage

Store at room temperature.

Shelf Life

2 years from date of receipt under proper storage conditions.

Protocol:

1. Add 50-100 μ L of DNA at 1-20 μ g/mL in TE Buffer to each well of a 96-well microplate.
2. Add an equal volume (50-100 μ L) of DNA Coating Solution. Mix well by pipetting and incubate 2 hours or overnight at room temperature on an orbital shaker.
3. Remove unbound DNA with a wash solution compatible with downstream application, such as PBS or TBS.

Example Results

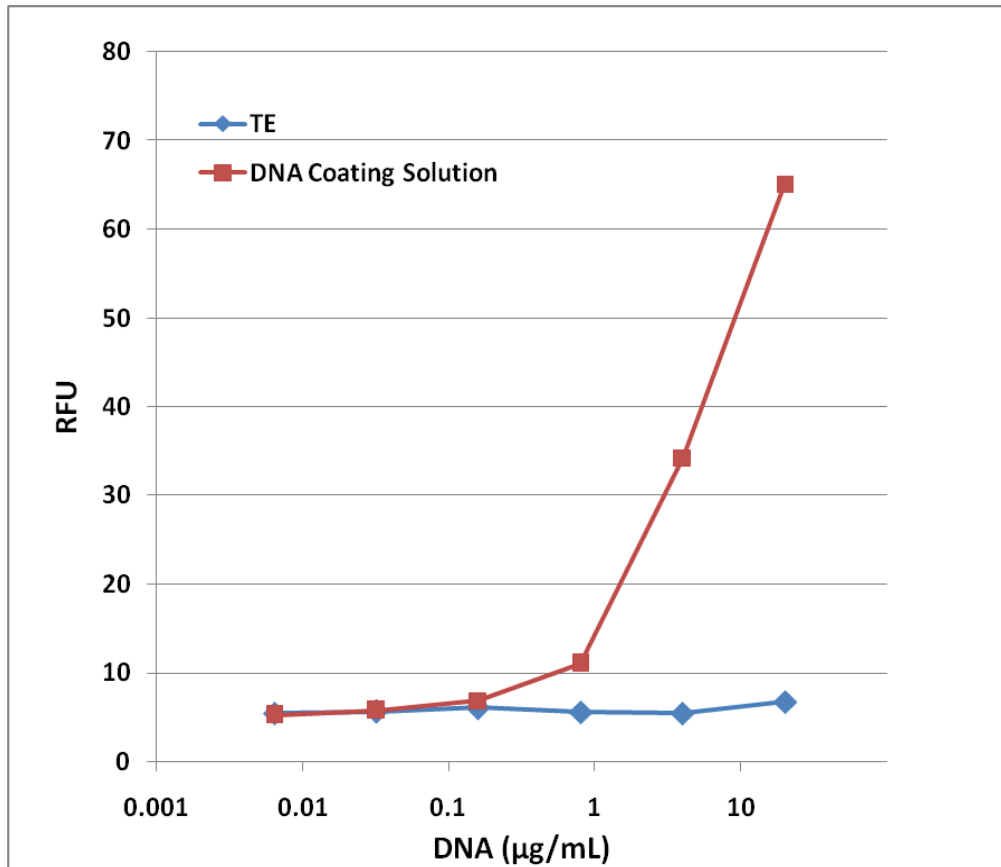


Figure 1: DNA Binding to 96-well Plate. 50 µL of Calf Thymus DNA solution in TE Buffer was mixed with 50 µL of either TE Buffer or DNA Coating Solution in a 96-well microplate. After a 2 hour incubation at room temperature, wells were washed with PBS and the adsorbed DNA in the wells was detected with CyQuant® GR Dye.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.