
DNA High Binding Solution

(Catalog No. R-1103)

Description

EpiGentek's proprietary DNA High Binding Solution provides a highly efficient method for binding DNA onto a microplate surface. This product allows the DNA to be tightly bound to the wells in a quick manner without the need for overnight incubations or heating microplates.

1. Binding incubation time is only 60 to 90 minutes
2. No blocking is required
3. Wide binding range of 0.1 ng to 200 ng
4. Ability to bind short DNA fragments, as short as 100 base pairs
5. DNA bound onto plastic surfaces is stabilized sufficiently to allow storage for 6 months prior to use

Volume

30 ml

Storage Temp & Conditions

Store at room temperature. DNA bound onto plastic surfaces is sufficiently stabilized to allow 6 months of storage, prior to use.

Application Notes

Used for DNA binding to microplate for various applications

Example Protocol

1. Add 100 µl of DNA high binding solution into the microplate wells and 1-8 µl of DNA solution (0.1-200 ng) for coating.
2. Incubate for 2 hours at 37°C.
3. Remove the solution from the wells and wash with PBS-T 2 times.
4. Dry at room temperature for 10-20 min.

The plate can now be immediately used or stored at 4°C for up to 6 months.

Ordering Information:

Products	Size	Cat. No.
DNA High Binding Solution	30 ml	R-1103 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.