

# EpiDyne® Remodeling Assay Substrate DNA ST601-GATC1,2,3, Biotinylated



EpiCypher®

**Catalog No.** 18-4113  
**Lot No.** 19004001  
**Pack Size** 50 µg

## Product Description:

EpiDyne® Remodeling Assay Substrate DNA ST601-GATC1,2,3, Biotinylated is a 217 base-pair double-stranded DNA fragment. This sequence includes the Lowary 601 nucleosome positioning sequence (see 18-0005) as well as a 3' acceptor sequence to accommodate the histone octamer subsequent to remodeling. ST601-GATC1,2,3 DNA has restriction enzyme sites embedded in the 601 sequence that is accessible after nucleosome remodeling.

## Formulation:

50 µg lyophilized ST601-GATC1,2,3 sequence DNA.

## Storage and Stability:

Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C.

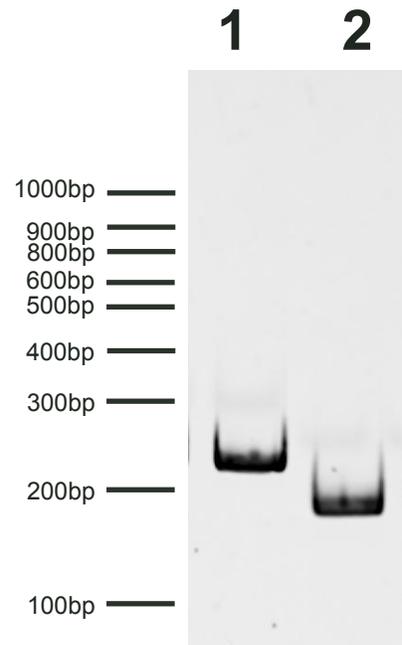
## Application Notes:

EpiDyne Remodeling Assay Substrate DNA ST601-GATC1,2,3, Biotinylated is useful as a negative control for restriction enzyme accessibility nucleosome remodeling assays using the Biotinylated- EpiDyne Remodeling Assay Substrate.

## DNA Sequence:

```
GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGATCGTAGA  
CAGATCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCGCGTT  
TTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCAGA  
TATATACATCGATGATGATGGATAGATGGATGATGGATGGATGGA  
TGATGATGGATGAATAGATGGATGGATGAAGCTT
```

## References:



## DNA Gel Data:

**Lane 1:** Free DNA (DNA, 200 ng). **Lane 2:** Free DNA incubated with 10U DpnII for 1 hr at 37°C.

This product is for *in vitro* research use only and is not intended for use in humans or animals.