

EpiDyne® Remodeling Assay Substrate DNA ST601-GATC1,2,3, 50-N-66, Biotinylated

Catalog No. 18-4116
Lot No. 19246001
Pack Size 50 µg



EpiCypher®

Product Description:

EpiDyne® Remodeling Assay Substrate DNA ST601-GATC1,2,3, 50-N-66, Biotinylated is a 263 base-pair double-stranded DNA fragment containing a 5' biotin-TEG group. This sequence includes the Lowary 601 nucleosome positioning sequence (see EpiCypher Catalog No. 18-0005), as well as both a 5' and 3' acceptor sequence to accommodate the histone octamer subsequent remodeling. ST601-GATC1,2,3, 50-N-66 DNA has restriction enzyme sites embedded in the 601 sequence that are accessible after nucleosome remodeling.

Formulation:

50 µg lyophilized ST601-GATC1,2,3, 50-N-66 Biotinylated DNA.

Storage and Stability:

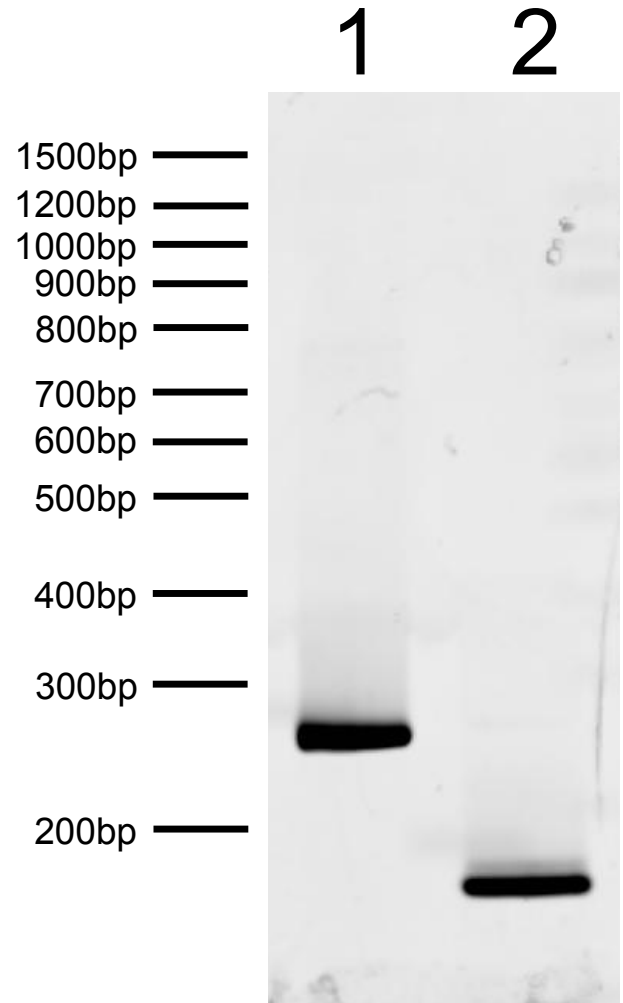
Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C.

Application Notes:

EpiDyne Remodeling Assay Substrate DNA ST601-GATC1,2,3, 50-N-66 Biotinylated is useful as a control for restriction enzyme accessibility nucleosome remodeling assays using the Biotinylated EpiDyne Remodeling Assay Substrate.

DNA Sequence:

```
GAACCAATGGGACCATGCTTCACACCGATATCATCGCTTATGTGTT  
GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGATCGTAG  
ACAGATCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCGCG  
TTTTAACCGCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCA  
GATATATACATCGATGATGATGGATAGATGGATGATGGATGGAT  
GGATGATGATGGATGAATAGATGGATGGATGAAGCTT
```



DNA Gel Data:

Lane 1: Free DNA (DNA, 100 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.