

# EpiDyne<sup>®</sup> Remodeling Assay Substrate DNA ST601-GATC1,2, Biotinylated



## EpiCypher<sup>®</sup>

**Catalog No.** 18-4112  
**Lot No.** 21169004-01  
**Pack Size** 50 µg

### Product Description:

EpiDyne Remodeling Assay Substrate DNA ST601-GATC1,2, Biotinylated is a 217 base-pair double-stranded DNA fragment containing a 5' biotin-TEG group and two GATC restriction enzyme sites. This sequence includes the 601 nucleosome positioning sequence [1], which has high affinity for histone octamers and is useful for nucleosome assembly. The DNA also includes a 3' acceptor sequence to accommodate the histone octamer subsequent to remodeling. When assembled onto a nucleosome, the GATC sites are shielded within the 601 sequence but can be made accessible for restriction enzyme cleavage after nucleosome remodeling. See the EpiDyne Nucleosome Remodeling Assay Tech Note (Restriction Enzyme Assay) for more information:

[epicypher.com/resources/technical-notes/](http://epicypher.com/resources/technical-notes/)

### Formulation:

50 µg lyophilized ST601-GATC1,2, 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.

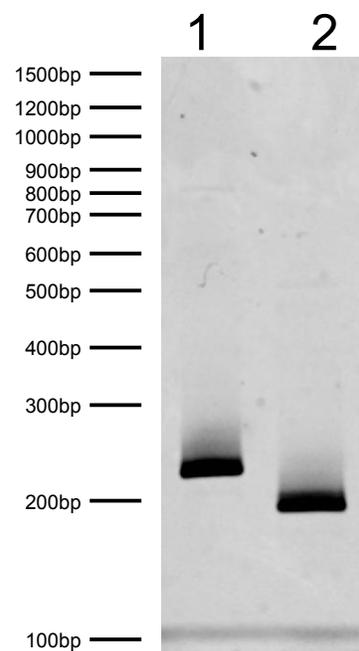
### Application Notes:

ST601-GATC1,2 DNA is useful as naked DNA template control for nucleosome remodeling assays using the EpiDyne Nucleosome Remodeling Assay Substrate ST601-GATC1,2 (Catalog No. 16-4112). See the EpiDyne Restriction Enzyme Assay Technical Note for more information:

[epicypher.com/resources/technical-notes/](http://epicypher.com/resources/technical-notes/).

### References:

[1] Lowary PT and Widom J (1998) *J Mol Biol* 276:19-42.



**DNA Gel Data:** ST601-GATC1,2 DNA resolved via native PAGE gel and stained with ethidium bromide for visualization. **Lane 1:** Free DNA (200 ng) and **Lane 2:** Free DNA digested with 10U DpnII at 37°C for 1 hour (200 ng). Migration patterns of DNA molecular weight markers are indicated.

### DNA Sequence:

```
GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGATCGTAGAC
AGCTCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCCGCGTTTT
AACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCAGATAT
ATACATCGATGATGATGGATAGATGGATGATGGATGGATGGATGATG
ATGGATGAATAGATGGATGGATGAAGCTT
```

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