

EpiDyne[®] Remodeling Assay Substrate DNA ST601-GATC0, Biotinylated



EpiCypher[®]

Catalog No 18-4110
Lot No 17258001
Pack Size 50 µg

Product Description:

EpiDyne[®] Remodeling Assay Substrate DNA ST601-GATC0 is a 217 base-pair double-stranded DNA fragment containing 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. This negative control does not contain DpnII restriction sites. When paired with the positive control DNA (Catalog No. 18-4111) these controls illustrate the migration range for the Restriction Enzyme Assay. See the EpiDyne Nucleosome Remodeling Assay Tech Note (Restriction Enzyme Assay) for more information:

epicypher.com/resources/technical-notes/

Formulation:

50 µg lyophilized ST601-GATC0 DNA.

Storage and Stability:

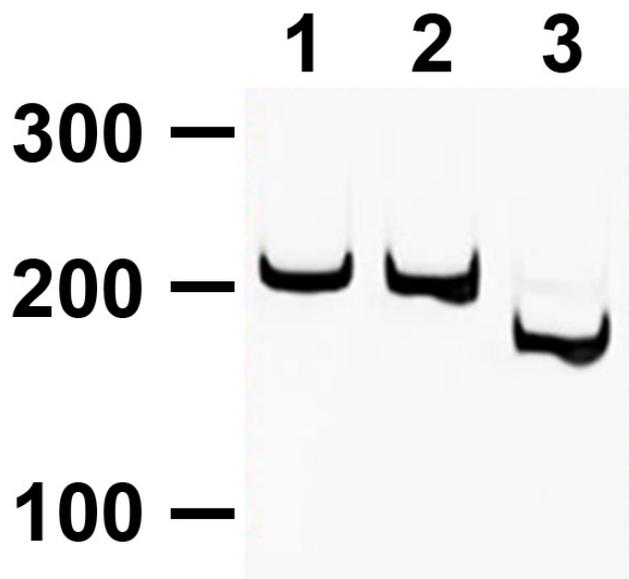
Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C.

Application Notes:

ST601-GATC0 DNA is useful as a negative control for restriction enzyme accessibility nucleosome remodeling assays using the EpiDyne Remodeling Assay Substrate. No DpnII restriction enzyme site is present, while the naturally occurring MfeI (AATTGG in bold) remains present within the 601 sequence.

References:

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



DNA Gel Data: ST601-GATC0 DNA resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (200 ng). **Lane 2:** Free DNA incubated with 10U DpnII for 1 hr at 37°C (200 ng). **Lane 3:** Free DNA incubated with 10U MfeI for 1 hr at 37°C (200 ng). Migration patterns of DNA molecular weight markers are indicated.

DNA Sequence:

```
GAATTCATCAGAATCCCGGTGCCGAGGCCGCTCAAATTGGTCGTAGAC  
AGCTCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCCGCTTTT  
AACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCAGATAT  
ATACATCGATGATGATGGATAGATGGATGATGGATGGATGGATGATG  
ATGGATGAATAGATGGATGGATGAAGCTT
```

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