

# EpiDyne® Nucleosome Remodeling Assay Substrate ST601-GATC1, 50-N-66, Biotinylated



## EpiCypher®

**Catalog No.** 16-4114  
**Lot No.** 19241001  
**Pack Size** 50 µg

### Product Description:

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4. Accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805) wrapped by provided 263 base pair DNA sequence that includes the 601 sequence with an added GATC sequence and a 5' biotin-TEG group.

### Formulation:

Purified recombinant mononucleosomes 1.62 mg/ml (DNA + protein weight) in 46.3 µl of 10mM Tris-HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. MW = 271,084.4 Da. Molarity = 3.98 µM.

### Storage and Stability:

Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

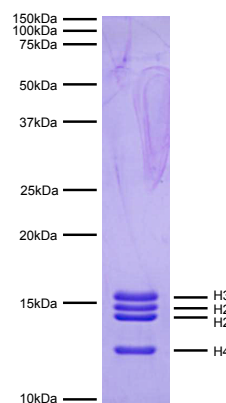
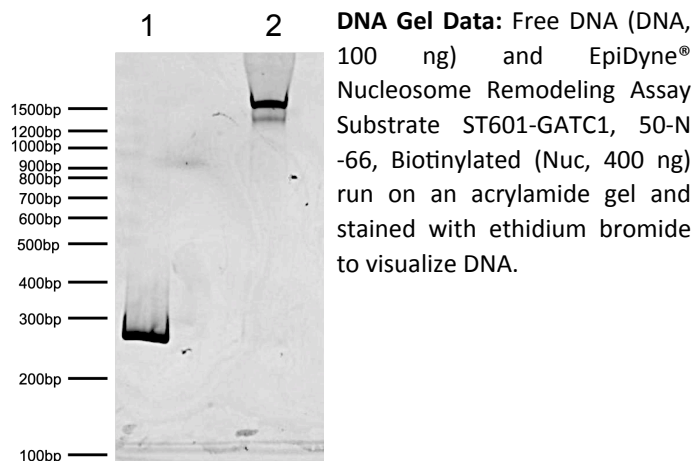
### Application Notes:

This product is a template for nucleosome remodeling assays using the restriction enzyme DpnII to determine accessibility of GATCs which is masked in its native configuration (prior to remodeling).

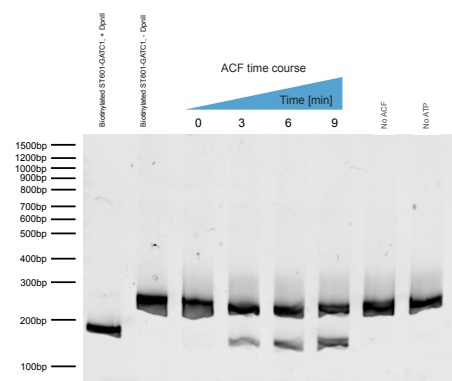
### DNA Sequence:

GAACCAATGGGACCATGCTTCACACCGATATCATCGCTTATGTGTT  
GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGGTCGTAG  
ACAGCTCTAGCACCGCTTAAACGCACGTACGCGCTGCCCCGCGT  
TTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCA  
GATATATACATCGATGATGATGGATAGATGGATGATGGATGGAT  
GGATGATGATGGATGAATAGATGGATGGATGAAGCTT

### References:



**Protein Gel Data:** Histone proteins in EpiDyne® Nucleosome Remodeling Assay Substrate ST601-GATC1, 50-N-66 Biotinylated resolved via SDS-PAGE and coomassie staining (2 µg). Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4) are indicated.



**Nucleosome Remodeling Data:** ACF/ATP-dependent nucleosome remodeling reaction in the presence of DpnII restriction enzyme. EpiDyne® Nucleosome Remodeling Assay Substrate ST601-GATC1, 50-N-66, Biotinylated nucleosomes incubated with (10 nM) or without ACF for up to 9 minutes in the presence of 2 mM ATP and 50U of DpnII. Samples were quenched at specified intervals and resolved via

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