

# EpiDyne™ Nucleosome Remodeling Assay Substrate ST601-GATC1, Biotinylated

Catalog No. 16-4111  
Lot No. 17258001  
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 217 base pair DNA sequence that includes the 601 sequence with an added GATC and a 5' biotin-TEG group.

## Formulation:

Purified recombinant mononucleosomes 1.62 mg/ml (DNA + protein weight) in 30.9 µl of 10mM Tris-HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. MW = 243,380 Da. Molarity = 6.67 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 control for nucleosome sliding assays using DpnII to determine accessibility of GATC.

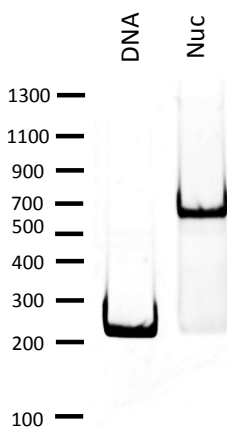
### DNA Sequence:

GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGGTCGTAG  
ACAGCTCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCGCGT  
TTTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCA  
GATATATACATCGATGATGATGGATAGATGGATGATGGATGGAT  
GGATGATGATGGATGAATAGATGGATGGATGAAGCTT

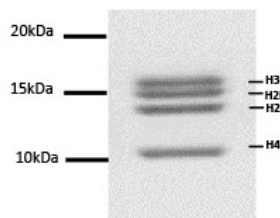
## References:



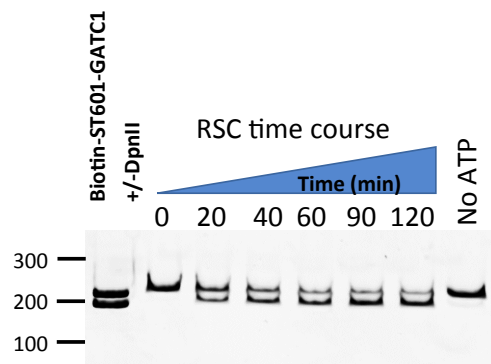
# EpiCypher™



**DNA Gel Data:** Free DNA (DNA, 200 ng) and EpiDyne Nucleosome Remodeling Assay Substrate Biotin ST601-GATC1 (Nuc, 370 ng) run on an acrylamide gel and stained with ethidium bromide to visualize DNA.



**Protein Gel Data:** Coomassie stained PAGE gel of proteins in EpiDyne Nucleosome Remodeling Assay Substrate Biotin ST601-GATC1 (2 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4) are indicated.



**Nucleosome Remodeling Data:** RSC/ATP-dependent nucleosome remodeling reaction in the presence of DpnII restriction enzyme. EpiDyne Nucleosome Remodeling Assay Substrate Biotin ST601-GATC1 nucleosomes were incubated with (10 nM) or without RSC remodeler for up to two hours in the presence of 2 mM ATP and 50U of DpnII. Samples were quenched at specified intervals and run on 8% native PAGE gel.

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