

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



EpiCypher®

Catalog No. 16-4116
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 three added GATC sequences and a 5' biotin-TEG group.

Formulation:

Purified recombinant mononucleosomes 0.97 mg/ml (DNA + protein weight) in 34.5 µl of 10mM Tris-HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. MW = 271,082.4 Da. Molarity = 5.35 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.

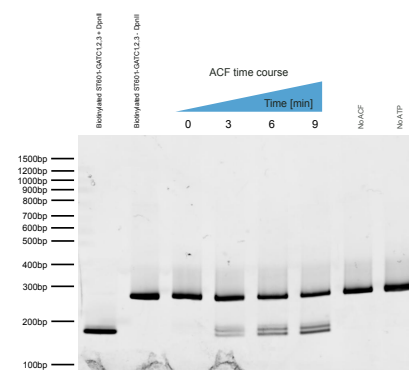
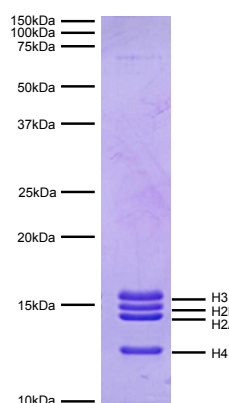
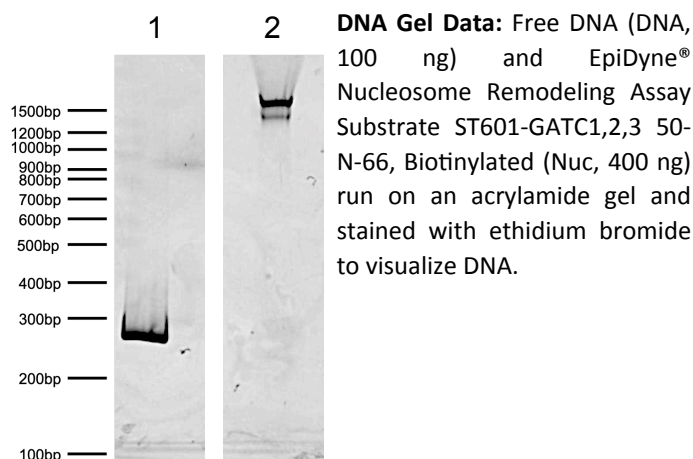
Application Notes:

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
GAATTCATCAGAATCCCGGTGCCGAGGCCGATCAATTGATCGTAG
ACA**GATC**TAGCACCGCTTAAACGCACGTACGCGCTGTCCCCGCG
TTTTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCA
GATATATACATCGATGATGATGGATAGATGGATGATGGATGGAT
GGATGATGATGGATGAATAGATGGATGGATGAAGCTT

References:



Nucleosome Remodeling Data: ACF/ATP-dependent nucleosome remodeling reaction in the presence of DpnII restriction enzyme. EpiDyne® Nucleosome Remodeling Assay Substrate ST601-GATC1,2,3, 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 SDS-PAGE.

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

8% native PAGE.



EpiCypher®

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

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com