

EpiDyne™-FRET Remodeling Assay Substrate DNA



EpiCypher™

Catalog No. 18-4201
Lot No. 18213001
Pack Size 50 µg

Product Description:

EpiDyne™-FRET Remodeling Assay Substrate DNA is a 207 base-pair double-stranded Cy3-conjugated DNA fragment. This sequence includes the Lowary 601 nucleosome positioning sequence (see 18-0005), a 5' Cy3 fluorophore, as well as a 3' acceptor sequence to accommodate the histone octamer subsequent to remodeling. The DNA also has a restriction enzyme site embedded in the 601 sequence that is accessible after nucleosome remodeling by the DpnII restriction enzyme.

Formulation:

50 µg EpiDyne™-FRET Remodeling Assay Substrate DNA lyophilized DNA.

Storage and Stability:

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

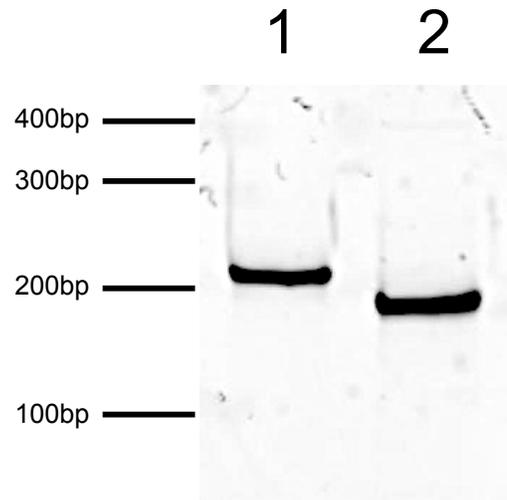
Application Notes:

This product is the DNA component of EpiDyne™-FRET Remodeling Assay Substrate, which can be used for nucleosome remodeling assays using Cy3/Cy5 FRET, or using the restriction enzyme DpnII to determine accessibility of GATC which is masked in its native configuration (prior to remodeling).

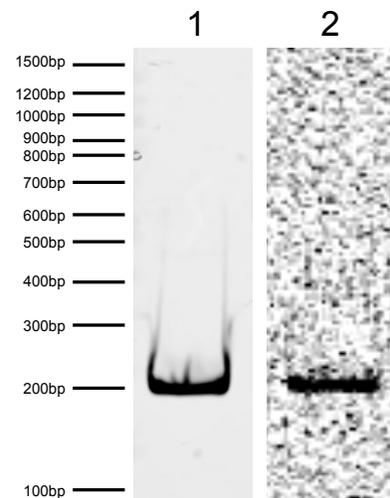
DNA Sequence:

Cy3_CATCAGAATCCCGGTGCCGAGGCCGATCAATTGGTCGT
AGACAGCTCTAGCACCGCTTAAACGCACGTACGCGCTGTCCC
CCGCGTTTTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGG
CACGTGTCAGATATATACATCGATGATGATGGATAGATGGAT
GATGGATGGATGGATGATGATGGATGAATAGATGGATGGAT
GAAGCTT

References:



Restriction Enzyme Data: DNA resolved via native PAGE and stained with ethidium bromide. **Lane 1:** Free DNA (DNA, 200 ng). **Lane 2:** Free DNA incubated with 10U DpnII for 1 hr at 37°C.



Fluorescence Data: DNA resolved via native PAGE and stained with ethidium bromide. **Lane 1:** DNA (200 ng) signal when excited at 312nm. **Lane 2:** DNA (200ng) signal when excited at 520nm.

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