

Nucleosome Assembly 601 Sequence DNA



EpiCypher®

Catalog No. 18-0006
Lot No. 21032002-54
Pack Size 50 µg

Product Description:

Nucleosome Assembly 601 Sequence DNA is a 147 base-pair double-stranded DNA fragment that was identified by Lowary and Widom using the SELEX method [1]. The 601 sequence DNA has high affinity for histone octamers and is useful for *in vitro* nucleosome assembly.

Formulation:

50 µg lyophilized 601 sequence DNA.

Storage and Stability:

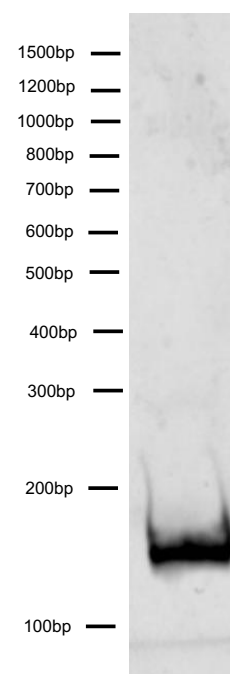
Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C.

Application Notes:

Nucleosome Assembly 601 Sequence DNA is useful for assembly of nucleosomes using purified or recombinant histone octamers (Catalog No. 16-0001). The biotin group on the DNA allows for pull-down of the nucleosomes for nucleosome binding experiments or enzymatic assays.

References:

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



DNA Gel Data: Nucleosome Assembly 601 Sequence DNA (75 ng) resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. Migration positions of DNA molecular weight markers are indicated.

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