

## Nucleosome Assembly 601 Sequence DNA

<b>Catalog No</b>	18-0006	<b>Tag</b>	None
<b>Lot No</b>	25351006-02	<b>MW</b>	90,716 Da
<b>Pack Size</b>	100 µg	<b>Source</b>	Synthetic DNA

### 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.

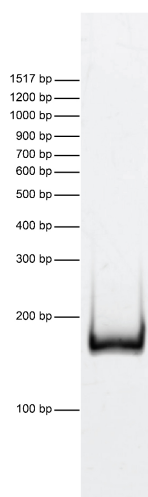
### TECHNICAL INFORMATION

<b>Storage</b>	Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C.
<b>Formulation</b>	100 µg lyophilized 601 sequence DNA.

### APPLICATION NOTES

Nucleosome Assembly 601 Sequence DNA is useful for assembly of nucleosomes using purified or recombinant histone octamers (EpiCypher 16-0001).

### VALIDATION DATA



**FIGURE 1 DNA gel data.** Nucleosome Assembly 601 Sequence DNA (100 ng) resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. Migration positions of DNA molecular weight markers are indicated.

### REFERENCES

[1] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715