

Nucleosome Assembly 601 Sequence DNA, 199 bp, Biotinylated

Catalog No	18-2044	Tag	Biotinylated
Lot No	24078002-01	MW	123,336.7 Da
Pack Size	50 µg	Source	Synthetic DNA

DESCRIPTION

Nucleosome Assembly 601 Sequence DNA, 199 bp, Biotinylated is a double-stranded DNA fragment with 26 bp linkers on either end of a 147 bp 601 Widom sequence, 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. The DNA contains a 5' biotin-TEG group.

TECHNICAL INFORMATION

Storage	Stable for 2 years at -20°C from date of receipt. After resuspending, aliquots should be stored at -80°C
Formulation	50 µg lyophilized 601 sequence DNA

APPLICATION NOTES

Nucleosome Assembly 601 Sequence DNA, 199 bp, Biotinylated is useful for assembly of nucleosomes using purified or recombinant histone octamers (EpiCypher 16-0001). See Luger et al. [2] for recommended nucleosome reconstitution protocol. This product also serves as a DNA only control for experiments utilizing the 199x601 nucleosome (EpiCypher 16-2044). These products are part of the Methyl DNA designer nucleosome set (epicypher.com/products/nucleosomes/methyl-dna-designer-nucleosomes).

VALIDATION DATA

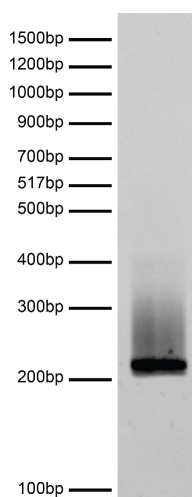


FIGURE 1 DNA gel data. Nucleosome Assembly 601 Sequence DNA, 199 bp, Biotinylated (100 ng) resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. Migration positions of DNA molecular weight markers are indicated.

REFERENCES

- [1] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715
- [2] Luger et al. *Methods Enzymol.* (1999). PMID: 10372352