

Nucleosome Assembly 601 Sequence DNA, Biotinylated

Catalog No	18-0005	Tag	Biotinylated
Lot No	23129011-01	MW	91197.5 Da
Pack Size	50 µg	Source	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. 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 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. The biotin group on the DNA allows for pull-down of the nucleosomes for nucleosome binding experiments or enzymatic assays.

VALIDATION DATA

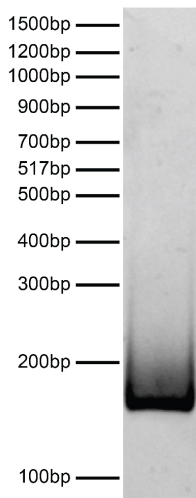


FIGURE 1 DNA gel data. Nucleosome Assembly 601 Sequence DNA, Biotinylated (75 ng) resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. Migration positions of DNA molecular weight markers are indicated.

REFERENCES

- [1] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715
- [2] Luger et al. *Methods Mol. Biol.* (1999). PMID: 10804500