

Nucleosome Assembly 601 Sequence DNA, Biotinylated

Catalog No. 18-0005

Lot No. 21082009-04

Pack Size 50 µg



EpiCypher®

Product Description:

Nucleosome Assembly 601 Sequence DNA, Biotinylated is a 147 base-pair double-stranded DNA fragment that was identified by Lowary and Widom (1998) using the SELEX method. The 601 sequence DNA has high affinity for histone octamers and is useful for *in vitro* nucleosome assembly. There is a biotin group on the 5' end of the fragment, which makes it ideal for use in nucleosome binding assays and pull-down experiments.

NOTE: This product replaces Catalog No. 18-0001.

Formulation:

50 µg lyophilized 601 sequence DNA containing a biotin group on the 5' end.

Storage and Stability:

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

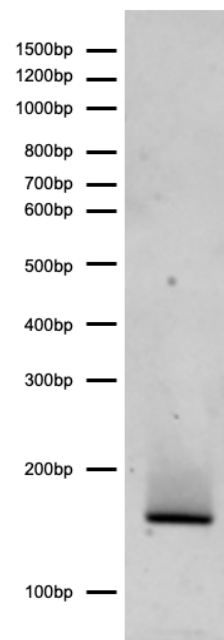
Application Notes:

Nucleosome Assembly 601 Sequence DNA, Biotinylated is useful for assembly of nucleosomes using purified or recombinant histone octamers (Catalog No. 16-0001). See Luger et al. (1999) for recommended nucleosome reconstitution protocol. The biotin group on the DNA allows for immobilization for binding experiments or enzymatic assays. See www.epicypher.com for more information and protocols.

References:

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

Luger et al (1999). *Methods Enzymol* 304:3-19.



DNA Gel Data: Nucleosome Assembly 601 Sequence DNA, Biotinylated (100 ng) resolved via 6% native PAGE in 0.4x TBE. 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.