

Nucleosome Assembly Control DNA

Catalog No: 53502

Format: 50 µg

Quality Control: A 187 base pair DNA fragment was PCR amplified from plasmid p601 (Widom, 1998). Gel analysis shows a band running at 187 base pairs with little or no contaminating primer dimer at approx. 50 base pairs.

Contents

Nucleosome Assembly Control DNA is provided lyophilized. Please resuspend DNA in sterile water or appropriate buffer as needed for nucleosome assembly.

Storage and Guarantee: Lyophilized DNA can be stored at -20°C . This product is guaranteed stable for 6 months from date of receipt when stored properly.

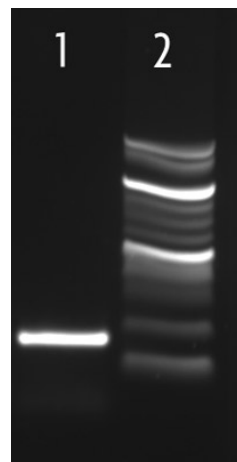


Figure 1:
5 ul of purified PCR product was loaded on a 2.5% agarose gel.
Lane 1: Nucleosome Assembly Control DNA
Lane 2: 100 bp DNA ladder