

Diversi-Phi Indexed PhiX

Catalog No: 53225

Format: 1 vial

Quality Control: Diversi-Phi Indexed PhiX is a ready-to-use, dual-indexed control library designed for Illumina sequencing platforms. It serves as a positive control to monitor sequencing quality and enhance nucleotide diversity, and is particularly beneficial for low-complexity libraries.

Diversi-Phi Indexed PhiX is a non-denatured double-stranded DNA dual-indexed PhiX library using all combinations of 8 x 8 bp indexes and provided as 10 µL of 10 nMol in 10 mM Tris pH 8.0.

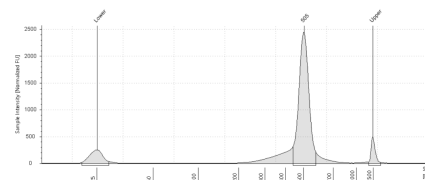
The Diversi-PHI Indexed PhiX Library is 55% Adenine + Thymine, and 45% Guanine + Cytosine.

Please visit www.activemotif.com to download sample sheet information.

Diversi-Phi Indexed PhiX is evaluated on TapeStation to ensure that a 500 bp library is produced. It is quantitated and provided at 10 nMol.

Sequencing recommendations: 2-5% for a typical pool with decent diversity (ATAC-seq, CUT&Tag, CUT&RUN, ChIP-Seq)

Storage and Guarantee: Store at -20°C. This product is guaranteed for 6 months from date of arrival when stored properly.



Diversi-Phi Indexed PhiX run on D1000 TapeStation ScreenTape. The single peak at 500 bp is indicative of the Diversi-Phi Indexed PhiX library.