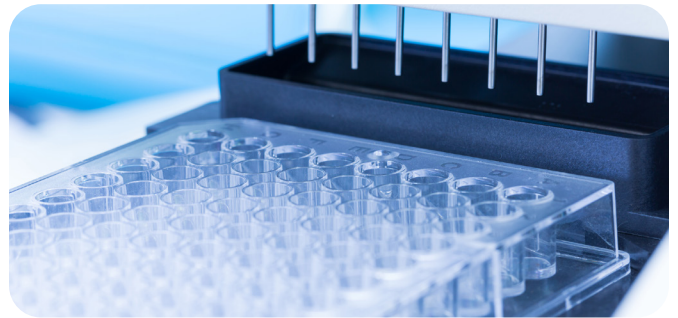




# NEXTFLEX Unique Dual Index Barcodes (set of 1,536)

## You Asked, We Listened!!

The new set of 1,536 NEXTFLEX® Unique Dual Index barcodes were designed to maximize flow cell capacity and minimize risk. The cost per sample is greatly reduced while mitigating chances of sample mix-up by means of unprecedented multiplexing options and enhanced barcode rotation capability between runs. The 1,536 adapters have unique 10 nucleotide indexes on both ends and each lot is functionally validated to increase data security by mitigating sample mis-assignment caused through index hopping and signal spreading, especially on sequencers with patterned flow cells such as the Illumina NovaSeq® sequencer or other Illumina® NGS platforms.



### Key features

- Save time and sequencing costs by multiplexing up to 1,536 samples in a single run on the high-throughput Illumina® NovaSeq® sequencer or any other Illumina® NGS platform
- Avoid costly mistakes due to sample swapping by having ready access to 1,536 unique dual index barcodes off-the-shelf
- Boost your confidence in sample identification using functionally validated unique dual indices

For research use only. Not for use in diagnostic procedures.

ADAPTERS

96 1536

MULTIPLY  
**16x**  
MORE LIBRARIES

### Get the Answer Faster

Samples: 1,536 Libraries | Sequencer: Illumina® NovaSeq® Sequencer | Run Type: S4 Reagent Kit

NOVASEQ®



384 BARCODES



TOTAL RUN TIME



1,536 BARCODES



### Reduce Sequencing Reagent Cost

Samples: 1,536 Libraries | Sequencer: Illumina® NovaSeq® Sequencer | Run Type: S4 Reagent Kit

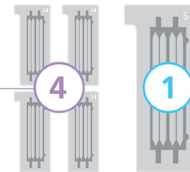
NOVASEQ®



384 SAMPLES



S4 FLOW CELLS



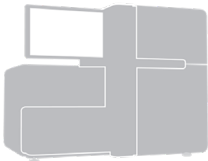
1,536 SAMPLES



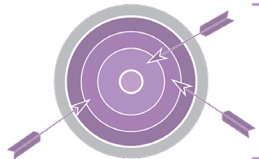
### Improve the Accuracy of Data Generated with Patterned Flow Cells

Sequencer: Illumina® HiSeq® Sequencer | Reagents: NEXTFLEX® Rapid XP DNA-seq Kit & 384 UDI Barcodes or Single Index Barcodes

HISEQ®

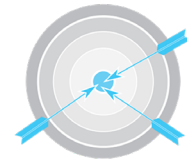


SINGLE INDEX



**0.2966%**  
INDICES MISASSIGNED

UNIQUE DUAL-INDEXED



**0.0013%**

## Ordering information

Catalog #	Kit name	Quantity
Nova-534100	NEXTFLEX® Unique Dual Index Barcodes (1,536 barcodes, 3,072 rxn)	3,072 rxn
Custom	NEXTFLEX® Unique Dual Index Pre-Plated Automation Barcodes (1,536 barcodes, 1,536 rxn)	1,536 rxn

Think of the possibilities! Learn more at [www.revvy.com](http://www.revvy.com)