

NEXTFLEX®

Small RNA-Seq Kit v4

Product number: NOVA-5132-31; -32; -41-44

NEXTFLEX SMALL RNA TRIMMING INSTRUCTIONS

Sequencing reads generated with the NEXTFLEX Small RNA Seq Kit v4 require trimming of adapter sequences prior to alignment when using and end-to-end alignment mode. These trimming steps can be accomplished using cutadapt (<https://cutadapt.readthedocs.io/en/stable/>), which is free to download.

The following commands require the latest version of cutadapt, which can be installed using the directions found at <http://cutadapt.readthedocs.io/en/stable/installation.html>.

1. Trim 3' adapter:

```
cutadapt \
--cores 0 \
--quality-cutoff 20 \
--adapter TGGAATTCTCGGGTGCCAAGG \
--output YOUR_FILE.cut.fastq \
--minimum-length 16 \
YOUR_FILE.fastq.gz
```

This command excludes any inserts less than 16 bases and removes bases from the 3' end of the insert with a quality score less than 20. Adjust values for minimum length and quality score as desired. It takes as input "YOUR_FILE.fastq.gz" and writes out to "YOUR_FILE.cut.fastq".

2. Trim 3' and 5' adapter from paired-end reads:

```
cutadapt \
--cores 0 \
--pair-adapters \
--quality-cutoff 20 \
--adapter TGGAATTCTCGGGTGCCAAGG \
-A AGATCGGAAGAGCGTCGTGTAGGGAAAGA \
--output YOUR_FILE.cut.R1.fastq \
--paired-output YOUR_FILE.cut.R2.fastq \
--minimum-length 16 \
YOUR_FILE.R1.fastq.gz \
YOUR_FILE.R2.fastq.gz
```

This command is similar to the single end version but requires paired reads to contain both adapters. If that is not the case, remove the "--pair-adapters" flag. If you would like to set a different quality score cutoff for read 2, use "--Q [value]" with "value" equal to the desired read 2 quality cutoff.

From here, you can align your trimmed fastqs to an appropriate reference using an aligner such as bowtie2.

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

