

NEBNext Immune Sequencing data analysis user guide

pRESTO tool introduction

The NEBNext Immune Sequencing data analysis workflow is based on the pRESTO tool suite, the Repertoire Sequencing TOolkit. pRESTO performs all stages of raw sequence processing prior to alignment against reference germline sequences. pRESTO is flexible and customizable and is composed of multiple modules. This tutorial is meant to be concise and allow you to understand and run an example workflow that is configured for the standard NEBNext Immune Sequencing workflow. For a more detailed tutorial please consider reading the [pRESTO tutorial](#).

In an example of an Illumina MiSeq paired-end 2x300 cycle run, Figure 1 (adapted from pResto documentation figures) shows the read schematic. Each read was sequenced from one end of the target cDNA so that the two reads together cover the entire variable region of BCR and TCR. The V(D)J reading frame proceeds from the start of read 2 to the start of read 1. Read 1 is in the opposite orientation (reverse complement), contains a partial C-region, and is 300 nucleotides in length. Read 2 contains the 5'RACE template switch site with a 17 nucleotide UMI barcode preceding it and is 300 nucleotides in length.

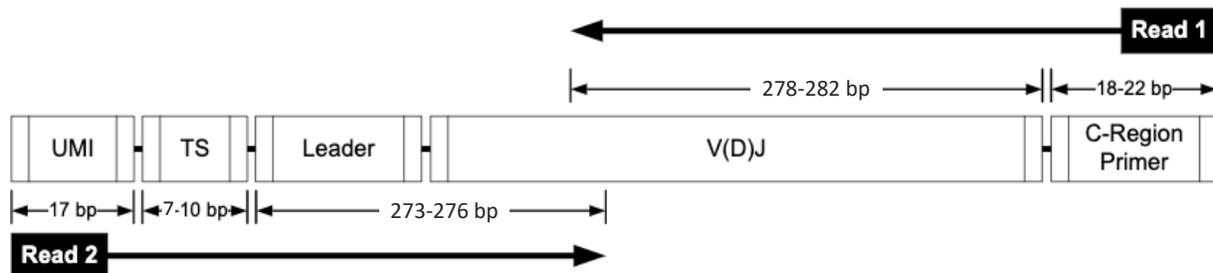


Figure 1. NEBNext Immune Sequencing library read schematic example of a run with Illumina MiSeq paired-end 2x300 reads.

pRESTO workflow overview



1. **Quality filter:** remove reads with mean Phred quality scores (Q) less than a user defined, sensitive value (set to 20%)
2. **Remove primers:** remove PCR primers and annotate reads with UMI (with maximum allowable error rate set to 20%)
3. **UMI consensus*:** generate consensus sequences for each UMI barcode by multiple alignment
4. **Sequence assembly:** assemble paired-reads on the UMI consensus into a long sequence (by default, when the reads do not intercept each other, the full fragment can be inferred using IgBlast)
5. **Filters:** quality-filter sequences
6. **Final repertoire:** obtain the final repertoire output that can be used for VDJ alignment tools like IgBlast. Only unique sequences with more than N (set to 2) representative reads are used in downstream analysis.

*Troubleshooting consensus generation

Workflow parameters have been selected based on our experience with Illumina MiSeq instruments and V3 reagent kits. Since each experiment accumulates experimental errors differently, the parameters may require tuning based on the following factors.

- Filter reads if there is a high rate of mismatches between UMIs in a UMI consensus (**maxerror parameter**).
- Filter reads that do not share common primers within a UMI consensus group (**prcons parameter**).
- Align C-regions to a C-InternalRegions.fasta file, filtering for no more than 30% errors and limited to 100 nucleotides.
- The annotation specifying the number of raw reads used to build each sequence, is updated to the minimum of Fwd and Rev reads.
- Duplicate sequences, sharing the same constant region, are removed.
- Sequences with more than n (20) "N" are also removed.

For more UMI consensus troubleshooting details, please go to the pRESTO tool webpage for [UMI consensus](#) or the following reference.

Dysregulation of B Cell Repertoire Formation in Myasthenia Gravis Patients Revealed through Deep Sequencing.

Vander Heiden JA, et al.

J Immunol. 2017 198(4):1460-1473. doi:10.4049/jimmunol.1601415.

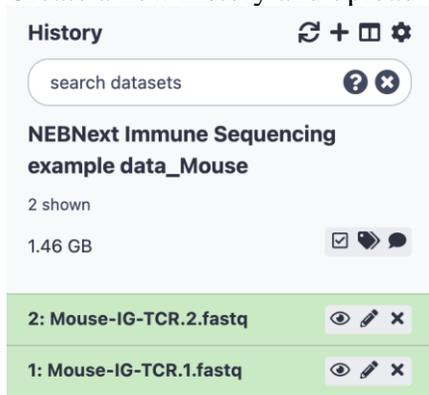
How to use the pRESTO workflow via Galaxy

NEB has implemented a pRESTO workflow on [Galaxy](#) for an easy start to users of immune sequencing data analysis. Galaxy is an open source, web-based platform for data intensive biomedical research. New users to Galaxy can follow the [Galaxy Tours](#) to start learning the platform. The pRESTO workflow published by NEB on Galaxy can be found [here](#).

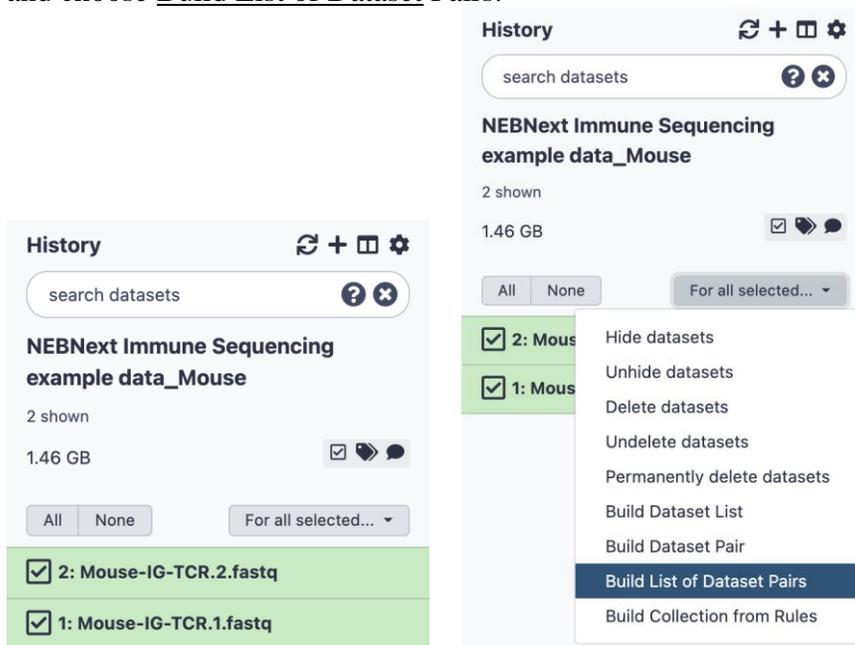
1. Import files before running pRESTO workflow

Input files for running pRESTO and example dataset can be found at an [example history](#) published by NEB. Follow the steps below to import your own files.

1.1 Create a new history and upload fastq files into the history.



1.2 Select all of the fastq files using the check button, click the “For all selected...” button and choose Build List of Dataset Pairs.



1.3 Enter the filters for forward and reverse fastq files, click Auto-pair, type the name of the collection, check that each R1 and R2 are correctly paired, and click Create list.

Create a collection of paired datasets

Could not automatically create any pairs from the given dataset names. You may want to choose or enter different filters and try auto-pairing again, **cancel** and reselect new elements.

Collections of paired datasets are ordered lists of dataset pairs (often forward and reverse reads). These collections can be passed to tools an...

1 unpaired forward - 1 filtered out Clear Filters Auto-pair 1 unpaired reverse - 1 filtered out

1.fastq 2.fastq

Mouse-IG-TCR.1.fastq Pair these datasets Mouse-IG-TCR.2.fastq

0 pairs Unpair all

Collections of paired datasets are ordered lists of dataset pairs (often forward and reverse reads). These collections can be passed to tools an...

0 unpaired forward - 0 filtered out Clear Filters Auto-pair 0 unpaired reverse - 0 filtered out

1.fastq 2.fastq

No datasets were found matching the current filters.

1 pairs Unpair all

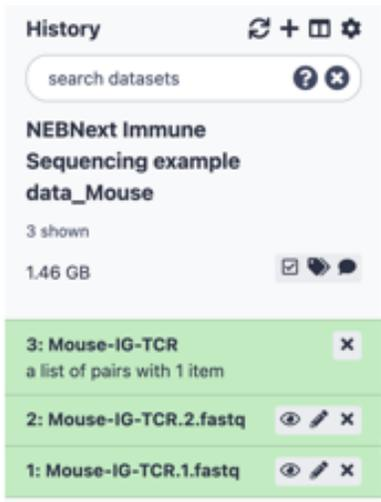
Mouse-IG-TCR.1.fastq → Mouse-IG-TCR. ← Mouse-IG-TCR.2.fastq

Hide original elements? Remove file extensions?

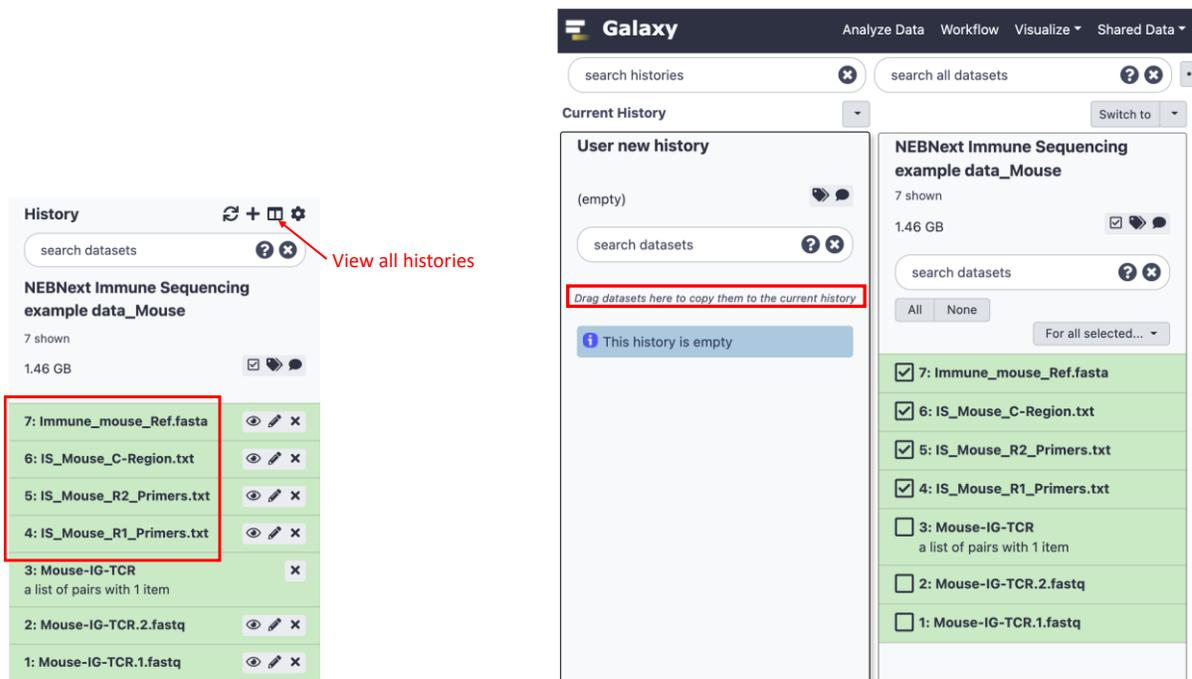
Name: Mouse-IG-TCR

Cancel Create list

the Galaxy Hub Please, help us



1.4 Import input files required to run pRESTO. The input files have been uploaded onto the [example history](#) as file number 4-7. The files can be copied to a new history or downloaded to local computer. To copy the files to a user history, click [view all histories](#) and click [Switch to](#) to make the destination history current, then select the files to be copied and drag them to the current history.



2. Run the pRESTO workflow
 - 2.1. Import the NEB [pRESTO workflow](#) to your Galaxy account.

2.2. Activate the history containing your data and input files. Click Workflow on the top menu and chose the Run Workflow button next to the pRESTO workflow.

The screenshot shows the pRESTO web interface. At the top, the 'Workflow' menu item is highlighted with a red box. Below the navigation bar, there is a search bar for workflows and buttons for '+ Create' and '+ Import'. A table of workflows is displayed with columns for Name, Tags, Updated, Sharing, and Bookmarked. One workflow, 'imported: pRESTO NEBNext Immune Sequencing Kit Workflow v3.2.0', is selected. To its right, a 'Run Workflow' button is highlighted with a red box. On the right side of the interface, a 'History' panel is visible, showing a search bar and a list of datasets including 'Immune_mouse_Ref.fasta', 'IS_Mouse_C-Region.txt', 'IS_Mouse_R2_Primers.txt', 'IS_Mouse_R1_Primers.txt', 'Mouse-IG-TCR', 'Mouse-IG-TCR.2.fastq', and 'Mouse-IG-TCR.1.fastq'.

2.3. Run the pRESTO workflow

- Fill out Workflow Parameters and History Options
 - Num Pairs: number of paired end reads to include in the analysis. Consistent subsampling ensures that each library has the same power to detect transcripts but selecting a very large number will cause all available reads to be used if that is preferred. Example number: 500000.
 - Send results to a new history: select Yes or No
 - 1: Input dataset collection: select the dataset collection in the history
 - 2: R1 Primer FASTA: select input file IS_Mouse_R1_Primers in the history
 - 3: R2 Primer FASTA: select input file IS_Mouse_R2_Primers in the history
 - 4: C-Region FASTA: select input file IS_Mouse_C-Region in the history
 - 5: Immune Ref FASTA: select input file Immune_mouse_Ref in the history
- Click the Run Workflow button.

Workflow: imported: pRESTO NEBNext Immune Sequencing Kit Workflow v3.2.0

Run Workflow

Input Datasets

3: Mouse-IG-TCR

Paired Fastq Dataset Collection

R1 Primer FASTA

4: IS_Mouse_R1_Primer.txt

Read 1 Primer Fasta

R2 Primer FASTA

5: IS_Mouse_R2_Primer.txt

Read 2 Primer Fasta

C-Region FASTA

6: IS_Mouse_C-Region.txt

C-Region Fasta

Immune Ref FASTA

7: Immune_mouse_Ref.fasta

Fasta containing known immune sequences (used to assemble read pairs that do not overlap)

NumReads

500000

decimal 0-0.9999 = fraction of total reads integer 1-N = number of reads

Expand to full workflow form.

History

search datasets

NEBNext Immune Sequencing example history_Mouse_run pRESTO

7 shown, 2 hidden

1.46 GB

- 9: Immune_mouse_Ref.fasta
- 8: IS_Mouse_C-Region.txt
- 7: IS_Mouse_R2_Primer.txt
- 6: IS_Mouse_R1_Primer.txt
- 5: Mouse-IG-TCR
- 2: Mouse-IG-TCR.2.fastq
- 1: Mouse-IG-TCR.1.fastq

3. QC report and output.

The sequencing reads QC metrics are visualized in the pRESTOr AbSeq3 Report. For further reads alignment, the output files included in Unique sequences and Unique sequences (>=2 reads) can be used as input for V(D)J alignment tools, for example IgBlast. An example history after running pRESTO is published [here](#).

History

search datasets

NEBNext Immune Sequencing example history_Mouse_run pRESTO

14 shown, 83 hidden

7.08 GB

- 96: Report on Unique sequences (>=2 reads)
- 93: Unique sequences (>=2 reads)
- 90: Report on Unique Sequences
- 88: Unique sequences
- 86: Report on final sequences
- 84: Final sequences
- 82: pRESTOr AbSeq3 Report on collection 70, collection 54, and others

