Cogent™ NGS Analysis Pipeline Quick Start Guide

The following information is provided as a high-level introduction to the software, also referred to as CogentAP. For more detailed information, please see the Cogent NGS Analysis Pipeline User Manual.

Before You Begin

- A. Supported operating systems
 - Linux: CentOS 8 or higher, RedHat 8 or higher, Ubuntu 18.04 or higher
- B. Hardware minimum requirements
 - CPU: 24 cores
 - Memory: 64 GB RAM
 - Free disk space: 1 TB

NOTE: If analyzing data generated with the ShastaTM Total RNA-Seq Kit or Shasta Whole-Genome Amplification Kit, at least 6 times the size of the input FASTQ files of free disk space is needed

- C. Additional dependencies
 - Internet connectivity on the server
 - Conda 24.4.0 or higher
 - Bash UNIX shell
 - bcl2fastq/BCL Convert
- D. Required input files
 - FASTQ files generated by an Illumina® sequencing platform.

NOTE: For a list of supported Takara Bio chemistries, please refer to our <u>bioinformatics</u> portal web page.

• A well-list text file, Illumina sample sheet, or similar TDT/CSV format file

Confirm Conda version

 Verify Conda is installed and meets or exceeds the required version by typing the following into a terminal window:

conda -V

If Conda is successfully installed, it should return text with the version number.

Example:

conda 24.4.0

 Verify the base Conda environment can be activated by typing:

conda activate

Type the following command to return to the default Linux prompt.

conda deactivate

- Verify the install location of miniforge3 is configured in the file .bash profile
 - a. For an individual user account, type:

```
more ~/.bash profile
```

b. Confirm something similar to the following is showing in the file (all on one line):

```
export
PATH="/home/<USERNAME>/miniforge3
/
bin:$PATH"
```

where <username > is replaced by the username of the account that installed Conda.

If no .bash_profile file exists or the line isn't displaying, it will need to be manually created and populated.

Installation

- 1. <u>Sign up</u> to download the installation package from our website.
- 2. Move or copy the ZIP file downloaded from Step 1 onto the Linux server into the directory location where you want to install.
- 3. Unzip the installation package by running the following two commands in the order listed:

```
unzip Cogent_NGS_Analysis_Pipeline_v3.1.zip \
&& mv Cogent_NGS_Analysis_Pipeline_v3.1 \
CogentAP
```

cd CogentAP

4. Run the following command to install CogentAP and its dependencies:

```
bash CogentAP setup.sh install
```

5. Install the human genome build (for the mouse genome build, use mm39 rather than hg38):

```
bash CogentAP_setup.sh genome_install
hg38
```

NOTE: See Section IV.D of the Cogent NGS Analysis Pipeline v3.1 User Manual for how to set up the \$COGENT AP HOME variable.



Generation of raw-fastq files

- 1. Log in to a server that stores the run folder from Illumina sequencing and has the bcl2fastq program installed.
- 2. Change to a working folder where you want the raw-fastq files to be located after being generated.
- 3. To convert BCL files to raw-fastq files using bcl2fastq, go to Step 3a. If using BCL Convert, go to Step 3b.
 - a. Run bcl2fastq with the following syntax template:

```
bcl2fastq -R <RUN_FOLDER> \
  -o <RUN_ID> \
  --no-lane-splitting \
  --sample-sheet \
  $COGENT_AP_HOME/config/SampleSheet_du
  mmy.csv > <RUN_ID>.stdout \
  2 > <RUN_ID>.stderr
```

where:

- <RUN_FOLDER> is the path to the sequencing run folder and
- <RUN_ID> is the ID number automatically generated by the Illumina sequencer

The file SampleSheet_dummy.csv is stored in the CogentAP config folder

Continue to Step 4.

b. Run BCL Convert with the following syntax template:

Templates for the DummySampleSheet for BCL Convert are stored in the CogentAP config/ folder. Continue to Step 4.

4. Move the raw-fastq files to your preferred storage location. They are typically generated in the <RUN ID> folder and named similar to:

```
Undetermined_ S0_R1_001.fastq.gz
Undetermined_ S0_R2_001.fastq.gz
```

RNA-Seq Analysis

• To demultiplex (demux) RNA-seq data:

```
$COGENT_AP_HOME/cogent rna demux \
-f <FASTQ_R1> \
-p <FASTQ_R2> \
-b <WELL-LIST> \
-t <EXP_TYPE> \
-0 <DEMUX_OUTPUT_DIR>
```

• To analyze RNA-seq data:

```
$COGENT_AP_HOME/cogent rna analyze \
-i <DEMUX_OUTPUT_DIR> \
-g <GENOME> \
-t <EXP_TYPE> \
-o <ANALYSIS_OUTPUT_DIR>
```

where:

- <FASTQ_R1> and <FASTQ_R2> are the full paths to the FASTQ files generated by an Illumina sequencing platform.
- <WELL-LIST> is the full path to the Shasta or ICELL8® system WellList, Illumina's sample sheet, or TDT/CSV format file
- <EXP_TYPE> is the experiment type used (e.g., icel18_fla, refer to the user manual for options)
- <DEMUX_OUTPUT_DIR> is the full path of the demultiplex results directory
- <GENOME> is a name of genome build (e.g., hg38)
- <ANALYSIS_OUTPUT_DIR> is the full path to the output directory created for the analysis results

NOTE: Additional commands are needed if analyzing sequencing data generated with the Shasta Total RNA-Seq Kit. See Section V.B of the Cogent NGS Analysis Pipeline v3.1 User Manual for details.

DNA-Seq or WGA Analysis

• To demultiplex (demux) DNA-seq data:

```
$COGENT_AP_HOME/cogent dna demux \
-f <FASTQ_R1> \
-p <FASTQ_R2> \
-b <WELL-LIST> \
-t <EXP_TYPE> \
-0 <DEMUX OUTPUT DIR>
```



To analyze DNA-seq data:

where:

- <COGENT_AP_HOME> is the path to the directory where CogentAP is installed
- <FASTQ_R1> and <FASTQ_R2> are the full paths to the FASTQ files generated by an Illumina sequencing platform.
- <WELL-LIST> is the full path to the Shasta, ICELL8 cx, or ICELL8 system WellList, Illumina's sample sheet, or TDT/CSV format file

- <EXP_TYPE> is the experiment type used (e.g., shasta_wga, refer to the user manual for more options)
- <DEMUX_OUTPUT_DIR> is the full path of the demultiplex results directory
- <GENOME> is a name of genome build (e.g., hg38)
- <BIN_SIZE> is the bin size used for CNV analysis using Ginkgo. Must be either 500kb or 1mb
- <READ_LENGTH> is the read length of the input data. Must be either 76bp or 151bp
- <READ_FILTER> is the minimum number of PE reads required per barcode to be kept in downstream analysis.
- <analysis_output_directory created for the analysis results
- <BARCODES_FILE> is the full path to the Shasta or ICELL8 system WellList, Illumina's sample sheet, or TDT/CSV format file.

Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: takarabio.com/service	web: takarabio.com/support
e-mail: ordersUS@takarabio.com	e-mail: technical_support@takarabio.com

Notice to Purchaser

Our products are to be used for Research Use Only. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2025 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara B io Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

Takara Bio USA, Inc.

2560 Orchard Parkway, San Jose, CA 95131, USA

United States/Canada: +1.800.662.2566 • Asia Pacific: +1.650.919.7300 • Europe: +33.(0)1.3904.6880 • Japan: +81.(0)77.565.6999

01.25 US

