

resDNASEQ® Quantitative CHO DNA Kit

Catalog Numbers 4402085 (kit), 4442731 (kit with protocol and quick reference card), 4413713 (combo kit†), and 4415413 (combo kit with protocol and quick reference card)

Publication Number 4415257 **Revision** B

Note: For safety and biohazard guidelines, refer to the “Safety” section in the *resDNASEQ® Quantitative CHO DNA Kit User Guide* (Pub. no. 4415260). For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product overview

The resDNASEQ® Quantitative CHO DNA Kit is used to quantitate host-cell residual DNA from Chinese hamster ovary (CHO) cells, a widely used cell line for production of biopharmaceutical products. Use the kit after you extract host-cell DNA from test samples. The resDNASEQ® Quantitative CHO DNA Kit uses TaqMan® quantitative PCR to perform rapid, specific quantitation of sub-picogram levels of residual CHO host-cell DNA. The assay is accurate and reliable across a broad range of sample types, from in-process samples to final product.

Kit contents and storage

Reagent	Description	Storage
resDNASEQ™ CHO DNA Control		
CHO DNA Control	1 tube, 40 µL	Store at –15 to –25°C.
DNA Dilution Buffer (DDB)	1 bottle, 7 mL	Store at –15 to –25°C before first use. Store at 2–8°C after first use.
resDNASEQ™ CHO Real-Time PCR Reagents		
2X Environmental Master Mix	2 tubes, 0.75 mL/tube	Store at –15 to –25°C before first use, protected from light. Store at 2–8°C after first use.
10X CHO DNA Real-Time PCR Assay Mix	1 tube, 300 µL	Store at –15 to –25°C, protected from light.
Negative Control	1 tube, 1.0 mL	Store at –15 to –25°C before first use. Store at 2–8°C after first use.

Prepare for quantitative CHO DNA analysis

Guidelines

When you prepare the serial dilutions of CHO DNA, follow these rules to *avoid carryover contamination* and to ensure proper sample preparation and quantitative PCR (qPCR) of samples:

- Use Ambion® nonstick 1.5-mL tubes.
- Label the top of each tube for identification.
- Use pipettes for the serial dilutions and standard curve that are separate from those used for sample preparation or PCR set up.
- Prepare the serial dilutions in an area physically separate from the test-sample preparation area.

IMPORTANT! To assure accurate quantitative results, Life Technologies protocols call for true triplicate sample preparation and analysis. Extract each test sample in triplicate and perform a single PCR for each extraction. The instrument software then calculates a mean quantity and standard deviation for the triplicate samples, followed by a percent coefficient of variation from this data ($SD/ \text{Mean Quantity} \times 100 = \%CV$). Based on the method qualification results, you can then assign a %CV to ensure accurate results from each sample tested.

† combo = PrepSEQ® Residual DNA Sample Preparation Kit + resDNASEQ® Quantitative CHO DNA Kit

Prepare CHO DNA serial dilutions

Prepare serial dilutions of CHO DNA control from the same experiment to create a standard curve and determine sample recovery rate.

To prepare the tubes for the CHO DNA standard curve, use the following table:

Standard curve (SC) tube	Volume of indicated serial dilution (SD) tube (µL)	Volume of PCR mix (µL)	Amount of DNA (pg)
SC1	SD1: 33	66	3000
SC2	SD2: 33	66	300
SC3	SD3: 33	66	30
SC4	SD4: 33	66	3
SC5	SD5: 33	66	0.3
SC6	SD6: 33	66	0.03

To prepare the reaction master mix, use the following table:

Kit Reagents	Volume for 1 30-µL reaction (µL)	Volume for 36 30-µL reactions (µL)
Negative Control	2	72
10X Primer/probe Mix	3	108
2X Environmental Master Mix	15	540
DNA Template	10	NA
Total	30	720

To set up a 96-well PCR reaction plate, use the following table:

	1	2	3	4	5	6	7	8	9	10	11	12
										Standard Curve (pg)		
A	NTC [†]	NTC	NTC							3000	3000	3000
B	NEG [‡]	NEG	NEG							300	300	300
C										30	30	30
D										3	3	3
E	TS [§] -1	TS-1	TS-1							0.3	0.3	0.3
F	TS-2	TS-2	TS-2							0.03	0.03	0.03
G	TS-3	TS-3	TS-3									
H												

[†] NTC = no template control

[‡] NEG = negative control

[§] TS = test sample

Workflow to prepare the serial dilutions and the standard curve

Prepare CHO DNA serial dilutions for the standard curve

Step 1: Label 7 Ambion® nonstick 1.5-mL tubes: **SD1, SD2, SD3, SD4, SD5, SD6,** and **NTC.**



Step 2: Add 990 µL of DNA dilution Buffer (DDB) to tube SD1.



Step 3: Add 450 µL of DDB to tubes SD2, SD3, SD4, SD5, SD6, and NTC.



Step 4: Remove the tube of CHO DNA Control (30 ng/µL) from the freezer.



Step 5: Thaw on ice, vortex gently for 2 sec, and quick-spin.

Step 6: Add 10 µL of CHO DNA Control to the tube that is labeled SD1, then vortex thoroughly.



Step 7: Transfer 50 µL of the DNA from tube SD1 to tube SD2.



Step 8: Transfer 50 µL of DNA from the previous dilution tube to the next dilution tube, until you add DNA to tube SD6. After each transfer, vortex thoroughly.



Step 9: Store the CHO DNA dilution tube at 4°C for use on the day of preparation. Otherwise, store the tubes at -20°C and use within 1 week.



Prepare the reaction master mix

Step 1: Determine the number of controls and test samples to quantify.



Step 2: Thaw all kit reagents completely at room temperature.



Step 3: Prepare a PCR mix.



Add the master mix to the test and standard curve samples

Step 1: Prepare tubes for the CHO DNA standard curve.



Step 2: Prepare the negative control samples (1 tube labeled NEG).



Step 3: Prepare the test samples.

Step 4: Prepare the No Template Control samples (1 tube labeled NTC).



Step 5: Set up a 96-well PCR plate.



Step 6: Seal the plate with optical film, then quick-spin with a centrifuge rotor that is compatible with 96-well plates.

To run the preparation on the 7500 Fast instrument, go to the next page.

Create, run, and analyze a plate on the 7500 Fast instrument

To set up a plate layout, use the following plate example:

	1	2	3	4	5	6	7	8	9	10	11	12
										Standard Curve (pg)		
A	NTC	NTC	NTC							3000	3000	3000
B	NEG	NEG	NEG							300	300	300
C										30	30	30
D										3	3	3
E	TS-1	TS-1	TS-1							0.3	0.3	0.3
F	TS-2	TS-2	TS-2							0.03	0.03	0.03
G	TS-3	TS-3	TS-3									
H												

To set up the standard curve, use the following table:

Tube label	Row-wells	Task	Quantity	Label (pg)
SC1	A-10, 11, 12	Standard	3000	3000
SC2	B-10, 11, 12	Standard	300	300
SC3	C-10, 11, 12	Standard	30	30
SC4	D-10, 11, 12	Standard	3	3
SC5	E-10, 11, 12	Standard	0.3	0.3
SC6	F-10, 11, 12	Standard	0.03	0.03

To set up tests and controls, use the following table:

Tube label	Row-wells	Task	Quantity (pg)	Label
NTC	A-1, 2, 3	NTC	NA	NTC
NEG	B-1, 2, 3	NTC	NA	NEG
TS-1	E-1, 2, 3	Unknown	NA	Sample-1
TS-2	F-1, 2, 3	Unknown	NA	Sample-2
TS-3	G-1, 2, 3	Unknown	NA	Sample-3

To set the thermal cycling temperature and time, use the following table:

	AmpliQaq Gold® enzyme activation	PCR	
	Hold	40 Cycles	
		Denature	Anneal/extend
Temp (°C)	95	95	60
Time (min:sec)	10:00	0:15	1:00

Workflow to create the plate document, run the plate, and analyze the results

On the 7500 Fast instrument:

Create a plate document

Step 1: Select **Absolute Quantification** from the Template Assay drop-down list.



Step 2: Select **Standard 7500** from the Run Mode drop-down list.



Step 3: Enter **resDNAassay** in the Plate name field, then click **Next**.



Step 4: Click **New Detector** and enter/select variables (name = **CHO**, report dye = **FAM**, quencher dye = **none**, select detector color), then click **Create Another**.



Step 5: Click **New Detector** and enter/select variables (name = **IPC**, report dye = **VIC**, quencher dye = **none**, select detector color).



Step 6: Select the applicable set of wells for the samples, then select FAM/IPC detectors for each well.

Step 7: Set up the standard curve (select the wells, assign the tasks, and label the appropriate wells).



Step 8: Set up the test and controls (select the wells, assign the tasks, and label the appropriate wells), then click **Finish**.



Step 9: Select the **Setup** tab ► **Plate** tab, select and label the appropriate wells.



Step 10: In the **Instrument** tab, set thermal-cycling conditions (reaction volume = **30 µL**, reaction = **standard**, set temperature and time).



Step 11: Select **File** ► **Save as** and confirm that the file is named "resDNAassay", then select **SDS Templates (*.sdt)** in the "Save as type" drop-down list and close the plate document (template).



Run the plate

Step 1: In the SDS software, select **File** ► **New**, then navigate to the SDS Documents folder.



Step 2: Select the **resDNAassay** template file, then click **Open**.



Step 3: In the Plate Name field, enter **ResDNA_date of assay**, then click **Finish** in the New Document Wizard.

Step 4: Make necessary changes to the test sample labels.



Step 5: Load the plate on the instrument.



Step 6: Select the **Instrument** tab, save the document, then click **Start** to start the real-time qPCR run.



Analyze the results

Step 1: Select **Analysis** ► **Analysis Settings** from the **Results** tab.



Step 2: In the Analysis Settings window:

- Select **Manual Ct**,
- Enter **0.2** in the Threshold field,
- Select **Manual Baseline**,
- Enter **3** for Start (cycle),
- Enter **15** for End (cycle),
- Click **OK**.



Step 3: Click  in the toolbar to analyze the plate.

Step 4: Select the **Results** tab ► **Standard Curve** tab, then verify the Slope, Intercept, and R2 values.



Step 5: Right-click the **Standard Curve**, select **Export as JPEG**, then click **OK**. Alternatively, press **PrintScreen**, then paste the image in a WordPad file.



Step 6: Select **Report** tab ► **Report**, then review the mean quantity and standard deviation for each of the samples.



Step 7: Select **File** ► **Export** ► **Results**. In the "Save as type" drop-down list, select **Results Export Files (*.csv)**, then click **Save**.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.



For Research Use Only. Not for use in diagnostic procedures.

The information in this guide is subject to change without notice.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

NOTICE TO PURCHASER: PLEASE REFER TO THE resDNASEQ® QUANTITATIVE CHO DNA KIT PRODUCT INSERT AND USER GUIDE FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners. TaqMan and Ampliqa Gold are registered trademarks of Roche Molecular Systems, Inc., used under permission and license.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit lifetechnologies.com/support

lifetechnologies.com

17 December 2011

