

resDNASEQ™ Quantitative DNA Kits

For quantification of residual genomic and plasmid DNA from cell lines used in the bioproduction process

Catalog Numbers 4402085, 4458435, A46014, A26366, 4464335, 4458441, 4464336, A46066, A41797, and A50337

Pub. No. MAN0030006 **Rev.** A

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *resDNASEQ™ Quantitative DNA Kits User Guide* (Pub. No. 4469836). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

For information about using resDNASEQ™ Quantitative DNA Kits with AccuSEQ™ Software v2.2 or later, see *resDNASEQ™ Quantitative DNA Kits User Guide* (Pub. No. 4469836).

■ Prepare the control serial dilutions (Genomic DNA)	1
■ Prepare the control serial dilutions (Plasmid DNA - Kanamycin Resistance)	2
■ Prepare the PCR reaction mix	3
■ Prepare the PCR plate	3
■ Setup, run, and analyze samples with AccuSEQ™ Software on the QuantStudio™ 5 Real-Time PCR Instrument	5
■ Limited product warranty	9

Prepare the control serial dilutions (Genomic DNA)

1. Label nonstick 1.5-mL microfuge tubes: **NTC**, **SD1**, **SD2**, **SD3**, **SD4**, **SD5**.
For CHO, Vero, MDCK, and NS0 kits, label an additional tube with **SD6**.
2. Add 50 µL of DNA Dilution Buffer to tube NTC. Put aside.
3. Add 990 µL of DNA Dilution Buffer to tube SD1.
4. Add 450 µL of DNA Dilution Buffer to tubes SD2, SD3, SD4, SD5, and (for CHO, Vero, MDCK, and NS0 only) SD6.
5. Remove the tube of DNA control (30 ng/µL) from the freezer.
6. After the DNA thaws, vortex it gently for 2 seconds, then briefly centrifuge.
7. Perform the serial dilutions:
 - a. Add 10 µL of the DNA control to the tube that is labeled SD1, then vortex thoroughly and briefly centrifuge.
 - b. Transfer 50 µL of the DNA from tube SD1 to tube SD2, then vortex thoroughly and briefly centrifuge.

- c. Continue to transfer 50 μ L of DNA from the previous dilution tube to the next dilution tube until you add DNA to tube SD5 (*E. coli*, HEK293, Human, *Pichia*, and Sf9 and Baculovirus) or SD6 (CHO, Vero, MDCK, and NS0). Final dilutions are shown in the table. After each transfer, vortex thoroughly, then centrifuge briefly.

Serial dilution (SD) tube	Dilution	pg DNA/reaction (10 μ L of the diluted DNA used in final 30 μ L of PCR reaction)
Control	DNA control tube	300,000 pg
SD 1	10 μ L DNA control + 990 μ L DDB	3,000 pg
SD 2	50 μ L SD1 + 450 μ L DDB	300 pg
SD 3	50 μ L SD2 + 450 μ L DDB	30 pg
SD 4	50 μ L SD3 + 450 μ L DDB	3 pg
SD 5	50 μ L SD4 + 450 μ L DDB	0.3 pg
SD 6 (for CHO, Vero, MDCK, and NS0 only)	50 μ L SD5 + 450 μ L DDB	0.03 pg

8. Store the DNA dilution tubes:

Temperature	For use
4°C	Same day
-20°C	≤ 1 week
-20°C	SD1 in single-use aliquots ≤ 6 months

Prepare the control serial dilutions (Plasmid DNA - Kanamycin Resistance)

1. Label nonstick 1.5-mL microfuge tubes: **NTC**, **SD1**, **SD2**, **SD3**, **SD4**, **SD5**, and **SD6**.

The dilution **SD6** will not be used for the standard curve. It will be used to confirm the limit of detection (LOD).

- Add 70 μ L of DNA Dilution Buffer to tube NTC. Put aside.
- Add 990 μ L of DNA Dilution Buffer to tube Dilution 1.
- Add 180 μ L of DNA Dilution Buffer to tubes SD1, SD2, SD3, SD4, and SD5.
- Add 100 μ L of DNA Dilution Buffer to tube SD6.
- Remove the tube of KanR DNA control (3.0×10^7 copies/ μ L) from the freezer.
- After the DNA thaws, vortex it thoroughly for 15-30 seconds, then tap gently to bring contents to the bottom of the tube. Do not centrifuge.
- Perform the serial dilutions:
 - Add 10 μ L of the KanR DNA control (3.0×10^7 copies/ μ L) to the tube that is labeled Dilution 1, vortex thoroughly for 15-30 seconds, then tap gently to bring contents to the bottom of the tube.
 - Transfer 20 μ L of the DNA from tube Dilution 1 to tube SD1, vortex thoroughly for 15-30 seconds, then tap gently to bring contents to the bottom of the tube..
 - Continue to transfer 20 μ L of DNA from the previous dilution tube to the next dilution tube until you add DNA to tube SD5.

- d. Transfer 100 µL of DNA from SD5 to SD6, then vortex thoroughly. Final dilutions are shown in the table. After each transfer, vortex thoroughly for 15-30 seconds, then tap gently to bring contents to the bottom of the tube. Do not centrifuge.

Serial dilution (SD) tube	Dilution	Concentration (copy number/ µL)	Copy number/ PCR reaction
Control	DNA control tube	3.0×10^7	N/A
Dilution 1	10 µL DNA control + 990 µL DDB	300,000	N/A
SD 1	20 µL Dilution 1 + 180 µL DDB	30,000	300,000
SD 2	20 µL SD1 + 180 µL DDB	3,000	30,000
SD 3	20 µL SD2 + 180 µL DDB	300	3,000
SD 4	20 µL SD3 + 180 µL DDB	30	300
SD 5	20 µL SD4 + 180 µL DDB	3	30
SD 6 (LOD)	100 µL SD5 + 100 µL DDB	1.5	15

9. Store the DNA dilution tubes:

Temperature	For use
4°C	≤2 days
-20°C	Dilution 1 ≤1 week
-20°C	Dilution 1 in single-use aliquots ≤6 months

Prepare the PCR reaction mix

- Determine the number of reactions needed for the controls and test samples that you will quantify.
- Thaw all kit reagents completely at room temperature, thoroughly mix reagents, and briefly centrifuge.
- Prepare a PCR reaction mix using the reagents and volumes shown in the table below.
 - Multiply the PCR reaction volume for one reaction (30 µL) by the number of reactions that you need to run.
 - Use 10% excess volume to compensate for pipetting losses.

Note: Use reagents from the same lot for all reactions.

Kit reagents	Volume for 1 30-µL reaction	Volume for 36 30-µL reactions (includes 10% overage)
Negative Control (Water)	2 µL	79.2 µL
10X DNA assay mix appropriate for the cell line being tested	3 µL	118.8 µL
TaqMan™ Environmental Master Mix 2.0	15 µL	594 µL
DNA template	10 µL	Add DNA template to each well separately, not as part of Master Mix
Total	30 µL	792 µL of Master Mix

Prepare the PCR plate

Plate setup differs slightly for each AccuSEQ™ System. See your software user guide for specific instructions. Place samples, NTCs, and standards in different quadrants of the plate.

Note: AccuSEQ™ Software v3.2 or later has a **Factory default/Admin Defined Template** for the Quantitative Sf9 and Baculovirus DNA Assay. Use the **resDNA_5Std** or **resDNA_6Std** template if setting up the Quantitative Sf9 and Baculovirus DNA assay in AccuSEQ™ Software v3.1. For more information about the **Factory default/Admin Defined Template** in AccuSEQ™ Software v3.2 or later, see “Create a Quantitative Sf9 and Baculovirus DNA experiment” on page 6.

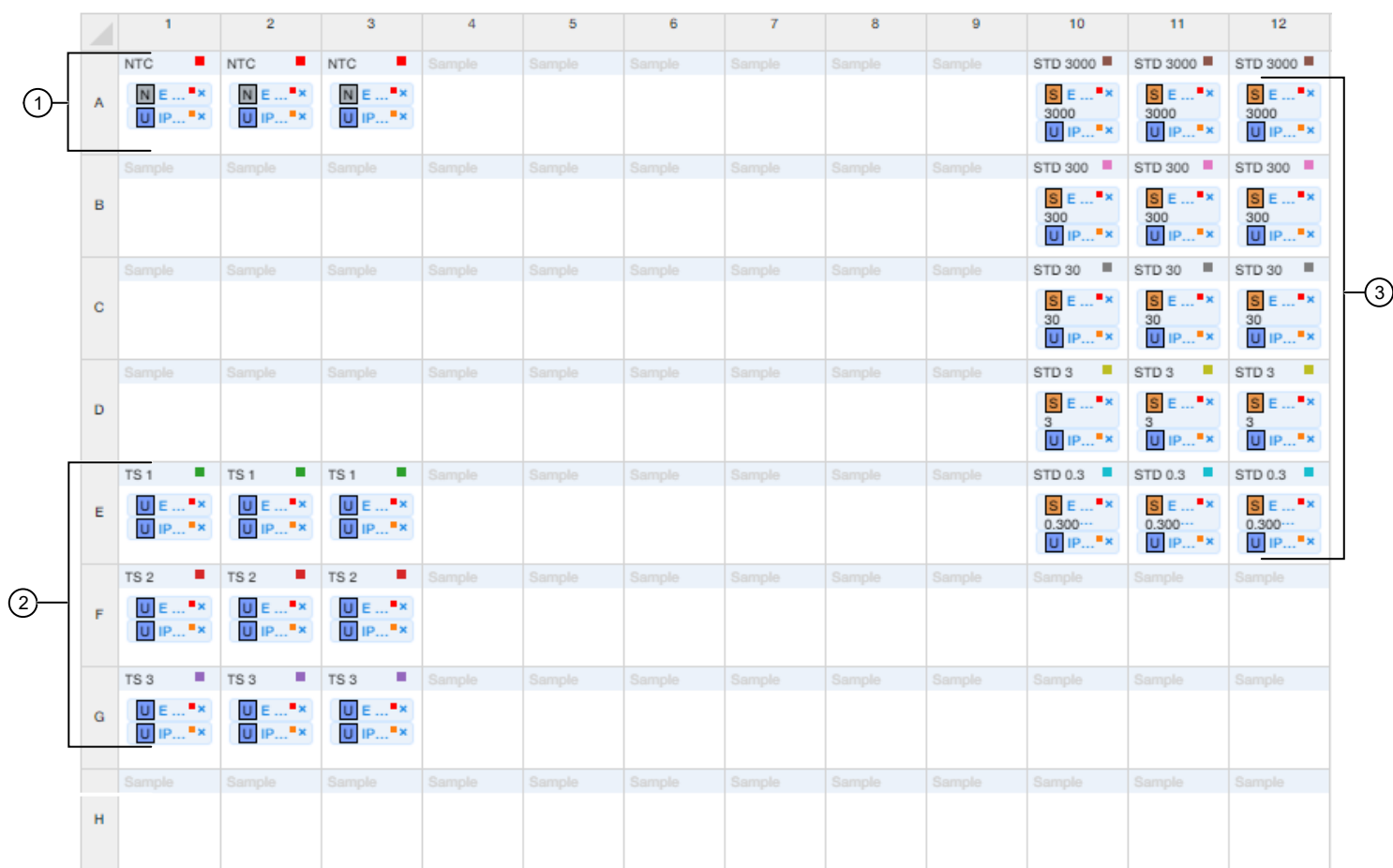


Figure 1 Default plate setup in the resDNA_5Std template in AccuSEQ™ Real-Time PCR Software v3.1 or later

① No template controls

③ Standard curve

② Samples

1. Add 20 μ L PCR reaction mix to each well.

2. Add 10 μ L of PCR NTC to the appropriate wells.

3. Add 10 μ L each of extracted sample DNA to the appropriate wells.

Note: If you prepared samples with the automated protocol, use a multichannel pipette to transfer the extracted sample.

4. Add 10 μ L of standard dilutions to the appropriate wells.

Note: Use different sets of pipettors to dispense test sample and standard curve dilutions to avoid cross-contamination of test samples.

5. Use a film applicator to seal the plate with optical film, then briefly centrifuge with a miniplate centrifuge that is compatible with 96-well plates.

Setup, run, and analyze samples with AccuSEQ™ Software on the QuantStudio™ 5 Real-Time PCR Instrument

Create a resDNASEQ™ experiment

1. In the  **Home** screen, click the **Factory default/Admin Defined Template** tab, then select a **resDNASEQ_5std** or **resDNASEQ_6std** template.

Note: AccuSEQ™ Software v3.2 or later has a **Factory default/Admin Defined Template** for the Quantitative Sf9 and Baculovirus DNA Assay. For more information about the **Factory default/Admin Defined Template** in AccuSEQ™ Software v3.2 or later, see “Create a Quantitative Sf9 and Baculovirus DNA experiment” on page 6.

Serial Dilutions (Standards)	Template	Assays
5	_5Std	<i>E. coli</i> , HEK293, Human, <i>Pichia</i>
6	_6Std	CHO, Vero, MDCK, NS0

2. In the **Experiment Properties** pane of the **Setup** tab:

- a. (Optional) Change the system-generated name of the experiment.

Note: Names must be unique. Deleted experiment names can not be reused.

- b. (Optional) Enter the plate **Barcode**, then add **Comments**.

Note: Comments are not editable post analysis.

Default resDNASEQ™ settings (cannot be changed)

- **Experiment Type** is **Quantitation-Standard Curve**
- **Chemistry** is **TaqMan™ Reagents**
- **Ramp Speed** is **Standard - 2hrs**

- c. Click **Next**.

Note: Experiment names cannot be changed after this step.

3. In the **qPCR Method** pane of the **Setup** tab, view the default volume and cycling conditions (cannot be changed).

4. Click **Next**.

5. In the **Samples** pane of the **Setup** tab, enter the sample **Name** and **Dilution**. **Sample Volume** is not needed. Add more **Samples** if needed.

IMPORTANT! Do not change the **Targets**.

6. In the **Samples** pane of the **Setup** tab, scroll to the right, then enter the spike information.

For more information about plate setup, see the *AccuSEQ™ Real-Time PCR Software v3.2 User Guide* (Pub. No. MAN0029199).

7. Click **Next**.

The **Run** tab is displayed.

8. Experiments are auto-saved in the software. To save, exit the experiment. The software prompts you to save changes. Click **Yes**.


Note: Clicking  **Save As** creates a copy of the experiment.

9. (Optional) Click  **Print** to print the plate layout for use in preparing the reactions.

10. Assemble the PCR reactions using the instructions in this guide.

Create a Quantitative Sf9 and Baculovirus DNA experiment

Note: The **Quantitative Sf9 and Baculovirus DNA** template is available with AccuSEQ™ Software v3.2 or later.



1. In the  **Home** screen, click the **Factory default/Admin Defined Template** tab, then select a **Quantitative Sf9 and Baculovirus DNA** template.
2. In the **Experiment Properties** pane of the **Setup** screen:
 - a. (Optional) Change the system-generated name of the experiment.
Note: Names must be unique. Deleted experiment names can not be reused.
 - b. (Optional) Enter the plate barcode in the **Barcode** field, then add comments in the **Comments** field.
Note: Comments are not editable post analysis.
Default Quantitative Sf9 and Baculovirus DNA settings (cannot be changed)
 - **Experiment Type—Quantitation-Standard Curve**
 - **Chemistry—TaqMan™ Reagents**
 - **Ramp Speed—Standard - 2hrs**
 - c. Click **Next**.
Note: Experiment names cannot be changed after this step.
3. In the **qPCR Method** pane of the **Setup** screen, view the default volume and cycling conditions (cannot be changed).
4. Click **Next**.
5. In the **Samples** table in the **Plate Setup** pane, enter the sample name and the sample dilution. Sample volume is not needed. Click **Add** to add more samples if needed.

IMPORTANT! Do not change the **Targets**.

6. In the **Samples** table of the **Plate Setup** pane, scroll to the right, then enter the spike information.
For more information about plate setup, see *AccuSEQ™ Real-Time PCR Software v3.2 User Guide* (Pub. No. MAN0029199).
Note: The software calculates the pg of DNA in the initial sample and the percentage of recovered DNA

Table 1 Plate well descriptions

Name	Description
SC1 to SC6	Standard curve dilutions in triplicate
Matrix	24 samples in triplicate
NTC	No template control in triplicate

7. (Optional) Double-click wells to add comments. Comments can also be added post-analysis.
8. Click **Next**.
The **Run** screen is displayed.
9. Experiments are auto-saved in the software. To save, exit the experiment. The software prompts you to save changes. Click **Yes**.
Note: Clicking  **Save As** creates a copy of the experiment.
10. (Optional) Click  **Print** to print the plate layout for use in preparing the reactions.
11. Assemble the PCR reactions using the instructions in this guide.

Create a Quantitative Kanamycin experiment

Note: The **Quantitative Kanamycin** template is available with AccuSEQ™ Software v3.2 or later.

For information about creating a Plasmid DNA template using AccuSEQ™ Software v3.1, see *resDNASEQ™ Quantitative DNA Kits User Guide* (Pub. No. 4469836).

1. In the **Home** screen, click the **Factory default/Admin Defined Template** tab, then select a **Quantitative Kanamycin** template.
2. In the **Experiment Properties** pane of the **Setup** screen:
 - a. (Optional) Change the system-generated name of the experiment.
Note: Names must be unique. Deleted experiment names can not be reused.
 - b. (Optional) Enter the plate barcode in the **Barcode** field, then add comments in the **Comments** field.
Note: Comments are not editable post analysis.
Default Quantitative Kanamycin settings (cannot be changed)
 - **Experiment Type—Quantitation-Standard Curve**
 - **Chemistry—TaqMan™ Reagents**
 - **Ramp Speed—Standard - 2hrs**
 - c. Click **Next**.
Note: Experiment names cannot be changed after this step.
3. In the **qPCR Method** pane of the **Setup** screen, view the default volume and cycling conditions (cannot be changed).
4. Click **Next**.
5. In the **Targets** table of the **Plate Setup** pane, enter the molecular weight (g/mol) of the targets.

Targets (2) [Add](#)

Quantity	Molecular Weight (g/mol)
	<input type="text" value="0"/>
	0.0000

6. In the **Samples** table in the **Plate Setup** pane, enter the sample name and the sample dilution. Sample volume is not needed. Click **Add** to add more samples if needed.
7. In the **Samples** table in the **Plate Setup** pane, scroll to the right, then enter the spike information.
 For more information about plate setup, see *AccuSEQ™ Real-Time PCR Software v3.2 User Guide* (Pub. No. MAN0029199).
Note: The software calculates the mass (g) and % recovery.

Table 2 Plate well descriptions

Name	Description
SC1 to SC6	Standard curve dilutions in triplicate
X1 to X3, D1 to D4	24 samples in triplicate
NTC	No template control in triplicate

8. (Optional) Double-click wells to add comments. Comments can also be added post-analysis.
9. Click **Next**.
 The **Run** screen is displayed.
10. Experiments are auto-saved in the software. To save, exit the experiment. The software prompts you to save changes. Click **Yes**.
Note: Clicking **Save As** creates a copy of the experiment.
11. (Optional) Click **Print** to print the plate layout for use in preparing the reactions.
12. Assemble the PCR reactions using the instructions in this guide.

Start the run

Start the run in the AccuSEQ™ Software.

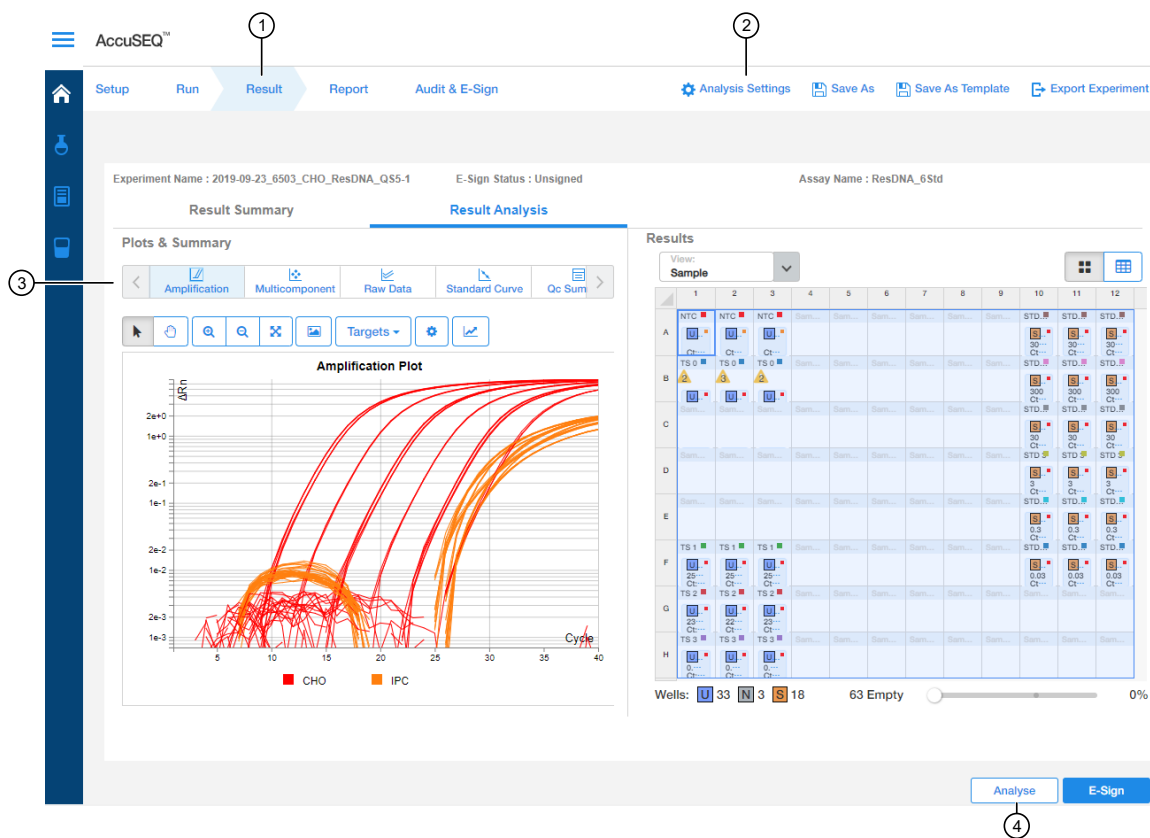
Option	Description
If the experiment is open	Click Start Run .
If the experiment is closed	<ol style="list-style-type: none"> Open the experiment. Click the Run tab. Click Start Run.

A message stating **Run has been started successfully** is displayed when the run has started.

Analyze the results

After the qPCR run is finished, use the following general procedure to analyze the results. For more detailed instructions see the *AccuSEQ™ Real-Time PCR Software v3.2 User Guide* (Pub. No. MAN0029199).

1. In the AccuSEQ™ Real-Time PCR Software, open the experiment, then navigate to the **Result** tab.



① **Result** tab

② **Analysis** Settings

③ Plot horizontal scrollbar

④ **Analyze** button

2. In the **Result Analysis** tab, review the Amplification Curve plots for amplification profiles in the controls, samples, and the standard curve. Ensure that auto threshold is selected.

Note: For AccuSEQ™ Real-Time PCR Software v3.1 Plasmid- KanR assay only, ensure that the threshold was set to 0.04 for Kanamycin target and 0.02 for IPC.

3. In the **Result Analysis** tab, review the **QC Summary** for any flags in wells.

4. In the **Result Analysis** tab, review the **Standard Curve** plot. Verify the values for the Slope, Y-intercept, R^2 , and Efficiency.

Note: The **Standard Curve** efficiency should be between 90-110% and the $R^2 > 0.99$. If these criteria are not met, up to two points, not in the same triplicate, can be removed from the standard curve data, and the analysis repeated.

5. (Optional) Navigate to the **Report** tab to generate a report of the experiment, or to export results.

6. (Optional for AccuSEQ™ Real-Time PCR Software v3.1) To manually convert the copy number output to a mass measurement, multiply the copy number given by the AccuSEQ™ Real-Time PCR Software to the average molecular weight of the plasmid, then divide by the Avogadro constant.

$$\text{Mass (g)} = \frac{\text{Copy number} \times \text{Molecular weight (gmol}^{-1}\text{)}}{6.0221 \times 10^{23} \text{ mol}^{-1}}$$

Note: AccuSEQ™ Real-Time PCR Software v3.2 and later calculates mass (g) and % Recovery for the **Quantitative Kanamycin** template in the **Results** table.

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Ltd | 7 Kingsland Grange | Woolston, Warrington WA1 4SR | United Kingdom

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0030006 A

Revision	Date	Description
A	5 December 2023	MAN0030006 Rev A is updated from 4469837 Rev K Update to include run and analysis information for AccuSEQ™ Real-Time PCR Software v3.2 on the QuantStudio™ 5 Real-Time PCR System and AccuSEQ™ Real-Time PCR Software v2.2 on the 7500 Real-Time PCR System. Remove the resDNASEQ™ Quantitative Synthetic Vero DNA Kit (Cat. No. A53242).
K	8 November 2021	Revision of 4469837, the previous part number for the resDNASEQ™ Quantitative DNA Kits Quick Reference. Update to include the resDNASEQ™ Quantitative Synthetic Vero DNA Kit (Cat. No. A53242).
J	28 March 2021	Revision of 4469837, the previous part number for the resDNASEQ™ Quantitative DNA Kits Quick Reference. Update to the control serial dilutions required for the resDNASEQ™ Quantitative Plasmid DNA - Kanamycin Resistance Gene Kit (Cat. No. A50337).

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021-2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.