

Version 01

Update: 07/18/2023

Product Manual

Pertuzumab Pharmacokinetic ELISA Kit

Cat. No. L00978

For Research Use Only. Not for Use in Diagnostic Procedures.

The operator should read the technical manual carefully before using this product.

Contents

| | |
|--|----|
| I. DESCRIPTION | 2 |
| II. ASSAY PRINCIPLE | 2 |
| III. ANALYTICAL CHARACTERISTICS | 2 |
| IV. KIT CONTENTS | 3 |
| V. STORAGE | 3 |
| VI. REAGENTS/EQUIPMENT NEEDED BUT NOT SUPPLIED | 3 |
| VII. PRECAUTIONS | 4 |
| VIII. SPECIMEN COLLECTION AND STORAGE | 4 |
| IX. PROTOCOL | 5 |
| ● Reagent Preparation | 5 |
| ● Capture Plate Preparation | 6 |
| ● Test Procedure | 7 |
| X. ASSAY PROCEDURE SUMMARY | 8 |
| XI. ANALYTICAL PERFORMANCE | 9 |
| ● Linearity and Limit of Detection | 9 |
| ● Precision and Accuracy | 10 |
| ● Selectivity | 10 |
| ● Dilutional Linearity and Hook Effect | 11 |
| ● Specificity | 11 |
| XII. TROUBLESHOOTING | 12 |
| XIII. REFERENCES | 13 |

I. DESCRIPTION

Pertuzumab, marketed under the brand name Perjeta, is a HER2/neu receptor antagonist. It works by preventing the formation of the HER2/HER3 dimer, which in turn blocks signaling by the dimer. Pertuzumab is used in combination with trastuzumab for the treatment of patients with HER2-positive tumors. It is a humanised IgG1 monoclonal antibody produced in mammalian cells.

GenScript's Pertuzumab Pharmacokinetic ELISA Kit had been comprehensively developed and validated for quantitative measurement of Pertuzumab in cynomolgus monkey serum and plasma, based on the ICH M10 and the FDA bioanalytical method validation guidance for industry. Its precision, accuracy, dilutional linearity, specificity, selectivity, stability, and hook effect were acceptable according to the guidances^[1-4]. The Pertuzumab ELISA kit is a validated tool for whole Pertuzumab and its biosimilar quantification in biological matrices for drug research and development.

II. ASSAY PRINCIPLE

Pertuzumab Pharmacokinetic ELISA Kit is a sandwich ELISA assay with a pair of anti-idiotypic monoclonal capture and detection antibodies. When standards or samples are added to the capture plate, the Pertuzumab in the sample can be captured on the plate coated with the Pertuzumab capture antibody. Then the Biotin Anti-Pertuzumab Antibody is added to interact with the Pertuzumab bound on the plate. Streptavidin-HRP (Streptavidin-Horseradish Peroxidase conjugate) is added to interact with the Biotin Anti-Pertuzumab Antibody. After the washing steps, TMB Solution is added, resulting in the formation of blue color. The reaction is stopped by adding Stop Solution. Adding the Stop Solution changes the color from blue to yellow. The intensity of the color can be read at 450 nm and 630 nm by a microplate reader.

The quantity of Pertuzumab in the sample is precisely quantified against a Pertuzumab standard curve.

III. ANALYTICAL CHARACTERISTICS

| Features | Specifications |
|---------------------------------|--|
| LLOQ | 6.25 ng/mL |
| ULOQ | 400 ng/mL |
| Intra-assay | CV ≤ 10% |
| Inter-assay | CV ≤ 15% |
| Minimum required dilution (MRD) | 1:50, validated non-human primate plasma |

IV. KIT CONTENTS

The kit provides the following reagents and solutions for the quantitative measurement of Pertuzumab and its biosimilar in biological matrices.

Table 1. Components of the kit

| Component | Quantity/Size | Part No. |
|--|----------------------|----------|
| Capture Plate | 1 plate | L1-80 |
| Standard Stock | 1 vial (50 μ L) | L1-10 |
| 100 \times Biotin Anti-Pertuzumab Antibody | 1 vial (160 μ L) | L1-20 |
| Streptavidin-HRP | 1 bottle (12 mL) | L1-30 |
| Sample Dilution Buffer | 1 bottle (60 mL) | L1-60 |
| Assay Dilution Buffer | 1 bottle (15 mL) | L1-90 |
| 20 \times Wash Solution | 1 bottle (60 mL) | L1-70 |
| Stop Solution | 1 bottle (6 mL) | A1-50 |
| TMB Solution | 1 bottle (12 mL) | A1-40 |
| Plate Sealer | 2 pieces | N/A |
| User Manual | 1 piece | N/A |

- Capture Plate: 96 well microplates (8 wells x 12 strips); 12 strips are configured in plate; plate is sealed in a foil pouch with a desiccant.
- Standard Stock contains 40 μ g/mL of Pertuzumab.

V. STORAGE

The unopened kit is stable for at least 12 months from the date of manufacture at 2°C to 8°C, and the opened kit is stable for up to 21 days from the date of opening at 2°C to 8°C.

VI. REAGENTS/EQUIPMENT NEEDED BUT NOT SUPPLIED

- Fresh matrix (normal serum or plasma from human or non-human primate)
- Microplate reader capable of measurement at 450 nm with the correction wavelength set at 630 nm
- Data analysis and graphing software. It is recommended to use software which is capable of generating a four-parameter logistic (4-PL) curve-fit
- Automated microplate washer
- Deionized or distilled water
- Graduated cylinder
- Plastic container
- Tubes to aliquot and dilute samples
- 10 μ L, 200 μ L, and 1000 μ L precision pipettes and pipette tips

- Multichannel pipettes
- Disposable reagent reservoir
- Absorbent paper
- Laboratory timer
- Refrigerator
- Centrifuge
- 25 ± 2 °C and 37 ± 2 °C incubator
- Rotary shaker
- Vortex Mixer

VII. PRECAUTIONS

1. All reagents containing human material should be handled as potentially infectious. Operators should wear gloves and protective clothing when handling any patient sera or serum based products.
2. Reagents that contain preservatives may be toxic if ingested, inhaled, or spilled on the skin.
3. Avoid contact with skin, eyes, or clothing with Stop Solution or TMB Substrate. Keep the container tightly closed. In case of an accident, please seek medical advice immediately.
4. Do not use the kit if there is any visible damage to the packaging or kit contents.
5. Do not mix components from different batches. Do not mix with components from other manufacturers.
6. Do not use reagents beyond the stated expiry date.
7. All reagents must be equilibrated to room temperature (20°-25°C) before running the assay. Only take an appropriate amount of reagents at once. Do not put unused reagents back into the vials as reagent contaminations may occur.
8. Before opening the Standard Stock, quickly span the vial to ensure that all the liquid has collected at the bottom, and prevent the liquid from splashing when opening the lid.
9. Use only distilled or deionized water and clean glassware.
10. Do not let wells dry during the test, add reagents immediately after completing washing steps.

VIII. SPECIMEN COLLECTION AND STORAGE

1. The handling and storage information provided here is intended to be used as a general guideline. Sample stability has not been evaluated. When samples need to be stored for a long time, users need to evaluate the stability of the samples. It is the responsibility of the individual laboratory to use all available references and/or its own studies when establishing alternate

stability criteria that meet their needs.

2. Store specimens at -20°C or lower if not tested immediately. Avoid repeated freeze-thaw cycles.

IX. PROTOCOL

- **Reagent Preparation**

All reagents must be equilibrated to room temperature before use (20°C-25°C). All samples and reagents should be vortexed before use. Store all reagents back in the refrigerator promptly after use.

1× Wash Solution: Dilute the 20× Wash Solution with deionized or distilled water with a volume ratio of 1:19. For example, dilute 40 mL of 20× Wash Solution with 760 mL of deionized or distilled water to make 800 mL of 1× Wash Solution. Store the solution at 2°C to 8°C when not in use.

Note: If any precipitate is found in the 20× Wash Solution, incubate the bottle in a water bath (up to 50°C) with occasional mixing until all the precipitate is dissolved.

1×Biotin Anti-Pertuzumab Antibody: Dilute the 100×Biotin Anti-Pertuzumab Antibody with Assay Dilution Buffer with a volume ratio of 1:99. For example, dilute 120 µL of 100×Biotin Anti-Pertuzumab Antibody with 11.88 mL of Assay Dilution Buffer to make 12 mL of 1×Biotin Anti-Pertuzumab Antibody. Store the solution at 2°C to 8°C when not in use.

Calibration Standard Preparation: Calibration standards should be prepared with a fresh matrix (The reagents are not provided in the kit) to generate eight Pertuzumab concentrations: fresh matrix (NC), 6.25, 12.5, 25, 50, 100, 200, and 400 ng/mL. Preparation of a whole set of standards is recommended in table 2. S-Int1 preparation is described below as an example.

Table 2. Recommended standard preparation

| Standard ID | Dilution Factor | Source | Source Volume (µL) | Matrix Volume (µL) | Final Volume (µL) | Final Conc. (ng/mL) |
|-------------|-----------------|------------------------|--------------------|--------------------|-------------------|---------------------|
| S-Int1 | 10 | Standard Stock (L1-10) | 10 | 90 | 100 | 4000 |
| Std1 | 10 | S-Int1 | 10 | 90 | 100 | 400 |
| Std2 | 2 | Std1 | 30 | 30 | 60 | 200 |
| Std3 | 2 | Std2 | 30 | 30 | 60 | 100 |
| Std4 | 2 | Std3 | 30 | 30 | 60 | 50 |
| Std5 | 2 | Std4 | 30 | 30 | 60 | 25 |
| Std6 | 2 | Std5 | 30 | 30 | 60 | 12.5 |
| Std7 | 2 | Std6 | 30 | 30 | 60 | 6.25 |
| NC | 0 | / | / | 60 | 60 | / |

Quality Control Preparation: QCs should be prepared with fresh matrix to generate five Pertuzumab concentrations: 6.25 (LLOQ), 18.75 (LQC), 75 (MQC), 300 (HQC) and 400 (ULOQ) ng/mL. Preparation of a whole set of standards is recommended in table 3. Q-Int1 preparation is described below as an example.

Note: QC is quality control. LQC is low quality control. MQC is medium quality control. HQC is high quality control. LLOQ is lower limit of quantification. ULOQ is upper limit of quantification.

Q-Int1 preparation: Vortex and centrifuge Standard Stock for several seconds. Dilute Standard Stock with a fresh matrix with a volume ratio of 1:10. For example, add 10 µL of Standard Stock to 90 µL of fresh matrix and mix it well to make 100 µL of Q-Int1.

Table 3. Recommended quality control preparation

| QC ID | Dilution Factor | Source | Source Volume (µL) | Matrix Volume (µL) | Final Volume (µL) | Final Conc. (ng/mL) |
|--------|-----------------|---------------|--------------------|--------------------|-------------------|---------------------|
| Q-Int1 | 10 | Stock (L1-10) | 10 | 90 | 100 | 4000 |
| ULOQ | 10 | S-Int1 | 10 | 90 | 100 | 400 |
| HQC | 1.3 | ULOQ | 60 | 20 | 80 | 300 |
| MQC | 4 | HQC | 15 | 45 | 60 | 75 |
| LQC | 4 | MQC | 15 | 45 | 60 | 18.75 |
| LLOQ | 3 | LQC | 20 | 40 | 60 | 6.25 |

- **Capture Plate Preparation**

It is recommended that all standards, quality controls, and samples be prepared in duplicate at least. Table 4 is an example for the setup of Pertuzumab standards and samples.

Count the strips according to the number of test samples and install the strips. Make sure the strips are tightly snapped into the plate frame.

Leave the unused strips in the foil pouch and store at 2°C to 8°C. The strips must be stored in the closed foil pouch to prevent moisture from damaging the Capture Plate.

Table 4. Setup of standards, quality controls and samples on Capture Plate

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|----------|------|------|------|------|-----|-----|-----|-----|-----|-----|-----|-----|
| A | Std1 | Std1 | ULOQ | ULOQ | S4 | S4 | S12 | S12 | S20 | S20 | S28 | S28 |
| B | Std2 | Std2 | HQC | HQC | S5 | S5 | S13 | S13 | S21 | S21 | S29 | S29 |
| C | Std3 | Std3 | MQC | MQC | S6 | S6 | S14 | S14 | S22 | S22 | S30 | S30 |
| D | Std4 | Std4 | LQC | LQC | S7 | S7 | S15 | S15 | S23 | S23 | S31 | S31 |
| E | Std5 | Std5 | LLOQ | LLOQ | S8 | S8 | S16 | S16 | S24 | S24 | S32 | S32 |
| F | Std6 | Std6 | S1 | S1 | S9 | S9 | S17 | S17 | S25 | S25 | S33 | S33 |
| G | Std7 | Std7 | S2 | S2 | S10 | S10 | S18 | S18 | S26 | S26 | S34 | S34 |
| H | NC | NC | S3 | S3 | S11 | S11 | S19 | S19 | S27 | S27 | S35 | S35 |

S: Sample number

- **Test Procedure**

Standards and Samples Incubation

1. Dilute standards, QCs and samples with Sample Dilution Buffer with a volume ratio of 1:50.

Note: Both standards and QCs are working solutions that have been diluted in the matrix, see PROTOCOL.Reagent Preparation for step details.

2. Add 100 μ L of the diluted standard solutions, controls and samples to the corresponding wells in the Capture Plate.
3. Cover the plate with Plate Sealer and incubate at 37°C for 60 minutes.
4. Remove the Plate Sealer and wash the plate with 260 μ L of 1 \times Wash Solution five times.
5. Tap the inverted plate onto absorbent paper to remove residual liquid in the wells after the wash steps.

Detection Antibody Incubation

6. Add 100 μ L of 1 \times Biotin Anti-Pertuzumab Antibody to all the testing wells.
7. Cover the plate with Plate Sealer and incubate at 37°C for 30 minutes.
8. Remove the Plate Sealer and wash the plate with 260 μ L of 1 \times Wash Solution five times.
9. Tap the inverted plate onto absorbent paper to remove residual liquid in the wells after the wash steps.

Enzyme Conjugate Incubation

10. Add 100 μ L of Streptavidin-HRP to all the testing wells.
11. Cover the Plate with Plate Sealer and incubate at 37°C for 10 minutes.
12. Remove the Plate Sealer and wash the plate with 260 μ L of 1 \times Wash Solution five times.
13. Tap the inverted plate onto absorbent paper to remove residual liquid in the wells after the wash steps.

Absorbance Measurement and Calculation

14. Add 100 μ L of TMB Solution to each well and incubate the plate in the dark at 25°C for 15 minutes (start timing after the addition of TMB Solution to the first well).

Note: TMB incubation time could extend to 20 minutes based on test signals.

15. Add 50 μ L of Stop Solution to each well to stop the reaction.
16. Read the absorbance in the microplate reader at 450 nm against 630 nm as a reference filter.
17. Plot the standard curve with the Pertuzumab concentration (ng/mL) on the x-axis and the corresponding mean absorbance value on the y-axis.
18. Using a 4- or 5-parameter logistic curve fitting program, calculate the best-fitting linear line through the points of the standard curve.

X. ASSAY PROCEDURE SUMMARY

- 1**
 - Prepare 1× Wash Solution, 1× Biotin Anti-Pertuzumab Antibody, and Capture Plate.
 - Dilute Standard Stock with **fresh matrix** to generate calibration standards and QCs
- 2**
 - Dilute the test samples and a set of standards and QCs with **Sample Dilution Buffer** with a volume ratio of 1:50.
- 2**
 - Add 100 µL of the diluted standard solutions, controls and samples to the corresponding wells. Incubate the plate at 37°C for 60 minutes.
- 3**
 - Wash the plate with 260 µL of 1× Wash Solution per well five times.
- 4**
 - Add 100 µL of 1× Biotin Anti- Pertuzumab Antibody to the well and incubate at 37°C for 30 minutes.
- 5**
 - Wash the plate with 260 µL of 1× Wash Solution per well five times.
- 6**
 - Add 100 µL of the Streptavidin-HRP and incubate at 37°C for 10 minutes.
- 7**
 - Wash the plate with 260 µL of 1× Wash Solution per well five times.
- 8**
 - Add 100 µL of TMB Solution and incubate the plate in dark at 25°C for 15 minutes.
- 9**
 - Add 50 µL of Stop Solution to each well to stop the reaction.
- 10**
 - Read the plate immediately.

XI. ANALYTICAL PERFORMANCE

- **Linearity and Limit of Detection**

A set of Pertuzumab calibration standards were freshly prepared and analyzed. Standard curves were constructed using a four- or five-parameter logistic curve. The typical dynamic range of the kit is 6.25-400 ng/mL (0.125-8 ng/mL diluted) and its detection limit is 6.25 ng/mL (Table 5 & Figure 1).

Table 5. Sample data for standard curve

| Pertuzumab (ng/mL) | Absorbance (OD 450/630nm) | | | Measured Pertuzumab (ng/mL) | CV % | Accuracy % |
|-----------------------|---------------------------|-------------|---------|-----------------------------------|---------|---------------|
| | Duplicate 1 | Duplicate 2 | Average | | | |
| 400 | 2.499 | 2.421 | 2.460 | 399.96 | 2.2 | 100.0 |
| 200 | 1.264 | 1.362 | 1.313 | 200.44 | 5.3 | 100.2 |
| 100 | 0.674 | 0.658 | 0.666 | 99.06 | 1.7 | 99.1 |
| 50 | 0.342 | 0.348 | 0.345 | 50.80 | 1.1 | 101.6 |
| 25 | 0.171 | 0.175 | 0.173 | 25.16 | 1.7 | 100.6 |
| 12.5 | 0.086 | 0.088 | 0.087 | 12.23 | 2.1 | 97.8 |
| 6.25 | 0.044 | 0.050 | 0.047 | 6.11 | 8.5 | 97.8 |
| N/A | 0.009 | 0.008 | 0.008 | / | 3.4 | / |

Pertuzumab Standard Curve

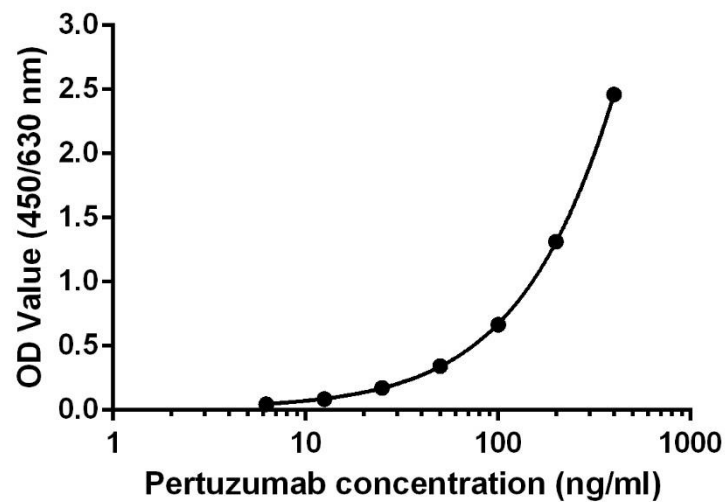


Figure 1: Pertuzumab ELISA kit standard curve.

A set of Pertuzumab calibration standards from 400 ng/mL to 6.25 ng/mL was then diluted with Sample Dilution Buffer with a volume ratio of 1:50.

- **Precision and Accuracy**

Pertuzumab Quality Controls at three concentrations (HQC of 300 ng/mL, MQC of 75 ng/mL, and LQC of 18.75 ng/mL) were tested for precision and accuracy in three batches of the kits, with each sample tested three times.

Table 6. Intra-assay and inter-assay precision and accuracy of the kit

| Quality Control | Pertuzumab (ng/mL) | Intra-assay (n=10) | | | Inter-assay (n=30) | | |
|-----------------|--------------------|-----------------------------|------|------------|-----------------------------|------|------------|
| | | Measured Pertuzumab (ng/mL) | CV % | Accuracy % | Measured Pertuzumab (ng/mL) | CV % | Accuracy % |
| HQC | 300 | 279.02 | 5.7 | 93.0 | 279.12 | 4.5 | 93.0 |
| MQC | 75 | 74.01 | 3.6 | 98.7 | 74.50 | 4.3 | 99.3 |
| LQC | 18.75 | 19.73 | 4.3 | 105.2 | 21.09 | 12.9 | 112.5 |

- **Selectivity**

Selectivity was tested by spiking plasma of ten different samples from non-human primate with Pertuzumab Quality Controls at two concentrations (HQC of 300 ng/mL and LLOQ of 6.25 ng/mL). Based on the ICH M10 guidance, the mean accuracy for LLOQ was required to be within 75%-125% of the low spiked concentration in at least 90% of the evaluated matrices. The mean accuracy for HQC was required to be within 80%-120% of the high spiked concentration in at least 100% of the evaluated matrices (Table 7).

Table 7. Selectivity analysis of the Kit

| HQC-Selectivity | | | LLOQ-Selectivity | | |
|-----------------------------|-----|-----------|-----------------------------|------|-----------|
| Measured Pertuzumab (ng/mL) | CV% | Accuracy% | Measured Pertuzumab (ng/mL) | CV% | Accuracy% |
| 304.62 | 1.5 | 101.5 | 5.14 | 1.2 | 82.2 |
| 296.90 | 5.3 | 99.0 | 5.07 | 6.3 | 81.1 |
| 285.26 | 2.1 | 95.1 | 4.69 | 1.9 | 75.0 |
| 293.24 | 0.2 | 97.7 | 4.70 | 5.2 | 75.2 |
| 294.98 | 1.1 | 98.3 | 5.70 | 4.4 | 91.2 |
| 309.41 | 0.7 | 103.1 | 5.66 | 0.8 | 90.6 |
| 305.39 | 5.6 | 101.8 | 5.76 | 1.7 | 92.2 |
| 315.36 | 3.0 | 105.1 | 5.37 | 5.6 | 85.9 |
| 284.85 | 4.8 | 95.0 | 5.15 | 19.5 | 82.4 |
| 277.20 | 8.5 | 92.4 | 4.63 | 5.1 | 74.1 |

- **Dilutional Linearity and Hook Effect**

Samples with high concentrations of Pertuzumab were used for the assessment of dilutional linearity (Table 9). The presence of a hook effect was investigated simultaneously. The hook effect was not observed in the assayed range (Table 8).

Table 8. Hook effect analysis of the kit

| Pertuzumab (ng/mL) | Absorbance (OD 450/630nm) | | | CV% |
|--------------------|---------------------------|-------------|---------|-----|
| | Duplicate 1 | Duplicate 2 | Average | |
| 20000 | 5.90 | 5.90 | 5.895 | 0.0 |
| 2000 | 5.91 | 5.30 | 5.604 | 7.7 |
| 1000 | 4.10 | 4.19 | 4.146 | 1.6 |

Table 9. Dilutional linearity analysis of the kit

| Dilution Factor | Expected Pertuzumab (ng/mL) | Measured Pertuzumab (ng/mL) | CV% | Accuracy% |
|-----------------|-----------------------------|-----------------------------|-----|-----------|
| 1000 | 20 | 19.04 | 7.4 | 95.2 |
| 2000 | 10 | 10.56 | 4.6 | 105.6 |
| 100 | 20 | 18.74 | 4.9 | 93.7 |
| 50 | 20 | 19.60 | 8.1 | 98.0 |

- **Specificity**

Pertuzumab QC samples at two concentrations (ULOQ of 400 ng/mL and LLOQ of 6.25 ng/mL) were spiked with different amounts of human IgG1 (800 and 8000 ng/mL). The test result demonstrated that the high concentration of human IgG1 did not interfere with the detection of Pertuzumab (Table 10).

Table 10. Specificity analysis of the kit

| Pertuzumab (ng/mL) | Human IgG1 (ng/mL) | Measured Pertuzumab (ng/mL) | CV% | Accuracy% |
|--------------------|--------------------|-----------------------------|-----|-----------|
| 400 | 4000 | 380.78 | 5.6 | 95.2 |
| 400 | 400 | 379.27 | 1.2 | 94.8 |
| 6.25 | 4000 | 7.80 | 2.0 | 124.8 |
| 6.25 | 400 | 7.74 | 3.4 | 123.8 |

XII. TROUBLESHOOTING

| Problem | Probable Cause | Solution |
|------------------------|---|---|
| Poor Precision | Wells are not washed or aspirated properly | Make sure the washing apparatus works properly and wells are dry after aspiration |
| | Wells are scratched with pipette tips or washing needles | Dispense and aspirate solution into and out of wells with caution |
| | Particulates are found in the samples | Remove any particulates by centrifugation prior to the assay |
| Weak/No Signal | Substrate is not added or added at the wrong time | Follow the manual to add the substrate properly |
| | Components are used from other lots or sources | Use only lot-specific components |
| | Substrate is contaminated | Use a new substrate from the same Lot |
| | Volumes of reagents are not correct | Repeat the assay with the required volumes as noted in the manual |
| | The plate is not incubated for proper time or temperature | Follow the manual to repeat the assay |
| | The plate is not read within the specified time range | Read the plate within 5 minutes |
| High Background | Plate is not washed properly | Make sure the washing apparatus works properly |
| | Substrate is contaminated | Use new substrate from the same Lot |
| | Evaporation of wells during incubations | Perform incubation steps with a plate sealer in a repeat assay |
| | Incorrect incubation times and/or temperatures | Follow the manual to repeat the assay |

XIII. REFERENCES

1. International Council for Harmonisation of Technical Requirements for Pharmaceuticals for Human Use. ICH Harmonised Guideline, Bioanalytical Method Validation M10 (2019).
www.ema.europa.eu/en/documents/scientific-guideline/draft-ich-guideline-m10-bioanalytical-method-validation-step-2b_en.pdf
2. US FDA. Bioanalytical Method Validation Guidance for Industry (2018).
www.fda.gov/files/drugs/published/Bioanalytical-Method-Validation-Guidance-for-Industry.pdf
3. European Medicines Agency. Guideline on Bioanalytical Method Validation (2011).
www.ema.europa.eu/en/documents/scientific-guideline/guideline-bioanalytical-method-validation_en.pdf
4. DeSilva B, Smith W, Weiner R et al. Recommendations for the bioanalytical method validation of ligand-binding assays to support pharmacokinetic assessments of macromolecules. *Pharm. Res.* 20(11), 1885–1900 (2003).

For research use only. Not intended for human and animal therapeutic or diagnostic use.

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China