



PathHunter[®] Pertuzumab Bioassay

Qualified with Perjeta[®]

[93-1042Y3-00095](#) (2-Plate Kit)

[93-1042Y3-00096](#) (10-Plate Kit)

OUR EXPERTISE
IN YOUR HANDS.
DISCOVER
CONFIDENTLY.

For Treatment of HER2+ breast cancer

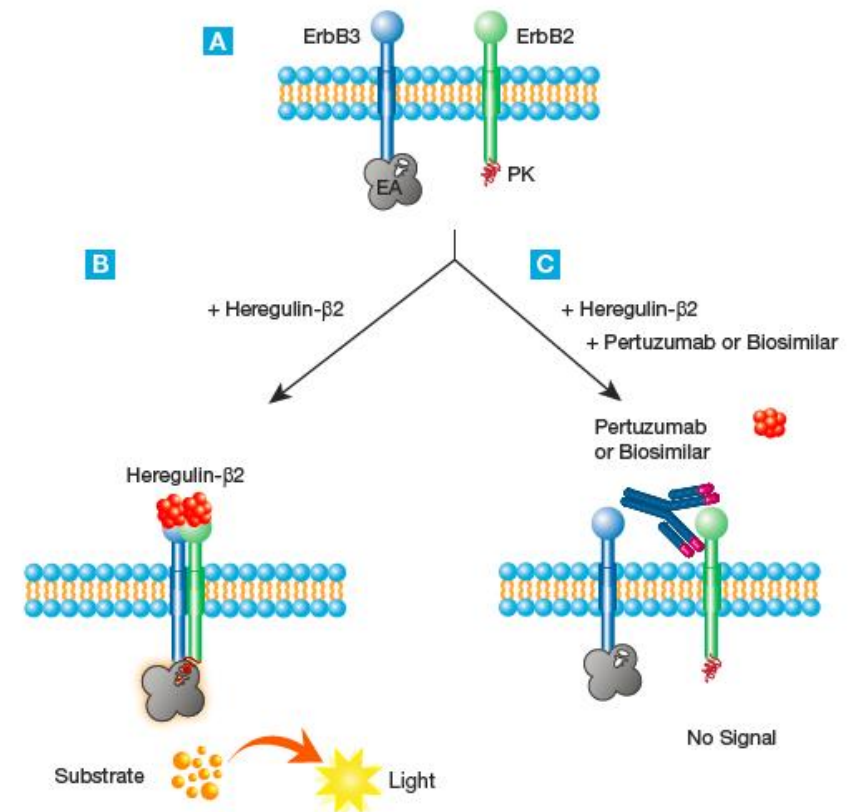
Mechanism of Action

HER2/ErbB2 is known to form homodimers as well as heterodimers with other ErbB family members (e.g. ErbB2, ErbB3) *in vivo*.

Heterodimerization with ErbB3 is the most potent combination for activating signaling pathways that lead to dysregulated growth / cell proliferation.

Pertuzumab is an IgG1 isotype mAb that acts by binding to the domain on ErbB2 that interacts with ErbB3, thus inhibiting the dimerization of ErbB2 with ErbB3, resulting in inhibition of downstream signaling.

Assay Principle



PathHunter[®] Pertuzumab Bioassay Kit

Kit Components

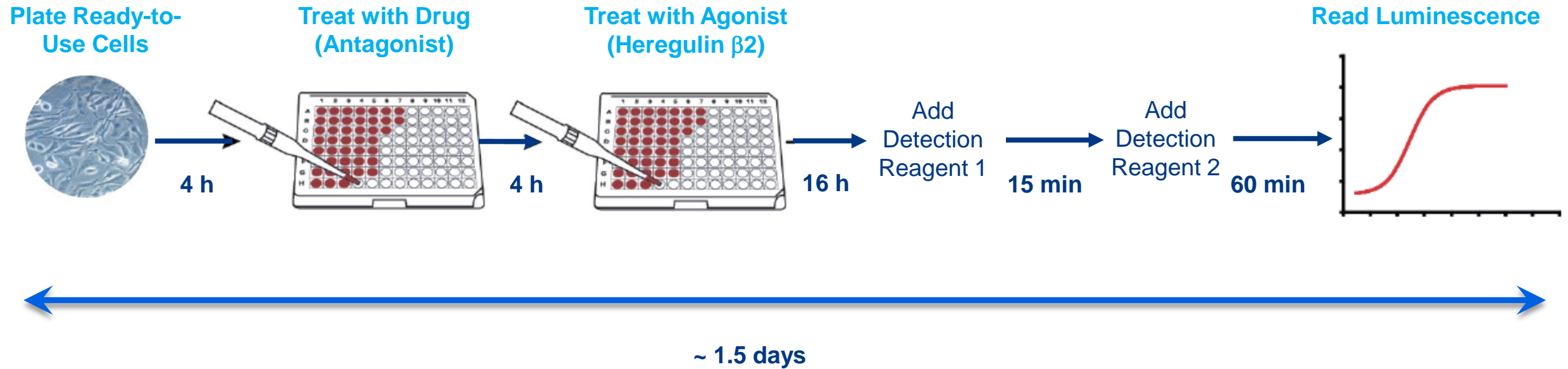
| List of Components | 93-1042Y3-00095 | 93-1042Y3-00096 |
|-------------------------------------------------------|-----------------|-----------------|
| PathHunter U2OS ErbB2/ErbB3 Bioassay Cells | 2 vials | 10 vials |
| PathHunter Bioassay Detection Kit | | |
| Detection Reagent 1 (mL) | 2 | 10 |
| Detection Reagent 2 (mL) | 8 | 40 |
| AssayComplete™ Cell Plating Reagent 5 | 1 X 100 mL | 3 X 100 mL |
| Protein Dilution Buffer | 1 X 50 mL | 2 X 50 mL |
| Control Agonist (Heregulin-β2) | 1 vial | 1 vial |
| 96-Well Clear-Bottom TC Treated, Sterile Plates w/Lid | 2 plates | 10 plates |

Sample data

Pertuzumab Bioassay Qualification

Bioassay Workflow

Simple, Homogenous and Rapid Protocol

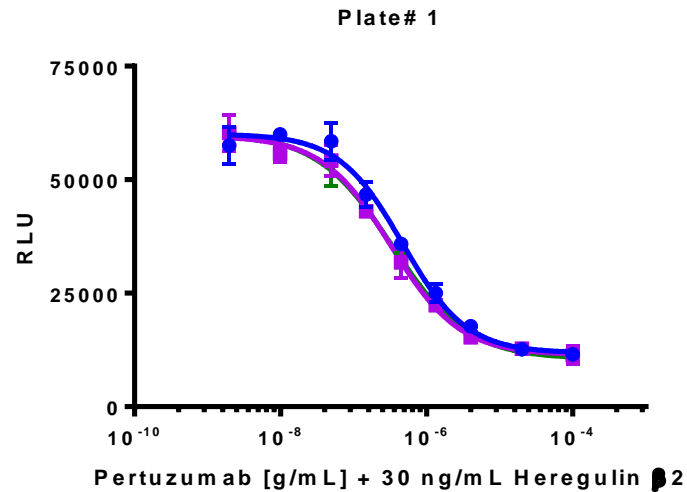


Assay Parameters Assessed

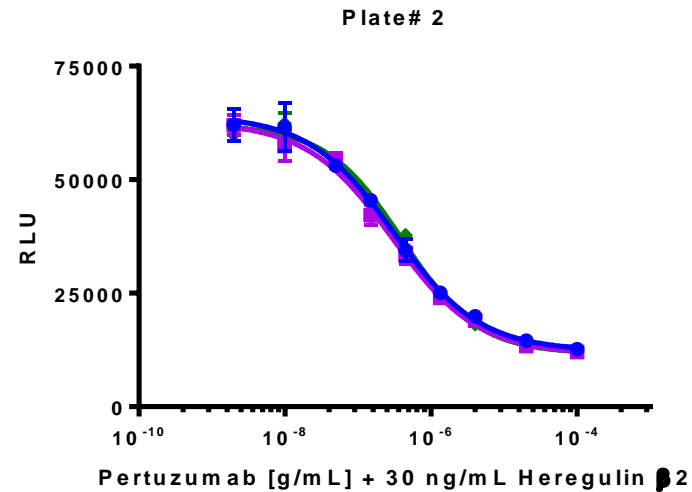
- % CV between 8 full plate DRCs
- Plate uniformity: EC_{80} and IC_{80} (of drug and stimulus) across entire plate
- Plate-to-Plate variability: 3 plates with full plate DRCs run on 3 days
- Slope consistency
- Relative potency across range of 50-150%
- Parallelism of relative potency data from two operators
 - Assay developer
 - Assay qualifier

Pertuzumab Bioassay Qualification with Perjeta®

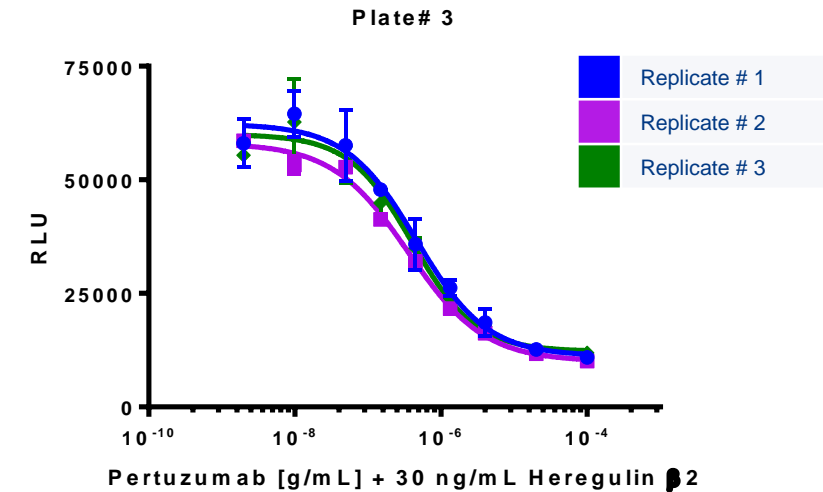
Plate to plate variability: 3 plates with full-plate DRC - Same Day



| Parameter | R1 | R2 | R3 |
|--------------------------|---------|---------|---------|
| S/B | 5.2 | 5.3 | 5.7 |
| Hill Slope | -0.9952 | -0.9197 | -0.8406 |
| IC ₅₀ (ng/mL) | 466.6 | 323.1 | 335.8 |



| Parameter | R1 | R2 | R3 |
|--------------------------|---------|---------|---------|
| S/B | 5.1 | 5.1 | 5.3 |
| Hill Slope | -0.7423 | -0.7404 | -0.8315 |
| IC ₅₀ (ng/mL) | 312 | 295.4 | 397.4 |



| Parameter | R1 | R2 | R3 |
|--------------------------|---------|---------|---------|
| S/B | 5.7 | 5.6 | 6.0 |
| Hill Slope | -0.9016 | -0.8358 | -0.9993 |
| IC ₅₀ (ng/mL) | 458.7 | 356.3 | 405.5 |

Pertuzumab Bioassay Qualification with Perjeta®

Assay Robustness

Intra-day Precision

| Plate # | Sample | S/B | %RSD, S/B | IC ₅₀ (ng/mL) | Mean IC ₅₀ (ng/mL) | %RSD IC ₅₀ |
|---------|--------|-----|-----------|--------------------------|-------------------------------|-----------------------|
| 1 | R1 | 5.2 | 5% | 466.6 | 375.2 | 21% |
| | R2 | 5.3 | | 323.1 | | |
| | R3 | 5.7 | | 335.8 | | |
| 2 | R1 | 5.1 | 2% | 312 | 334.9 | 16% |
| | R2 | 5.1 | | 295.4 | | |
| | R3 | 5.3 | | 397.4 | | |
| 3 | R1 | 5.7 | 4% | 458.7 | 406.9 | 13% |
| | R2 | 5.6 | | 356.3 | | |
| | R3 | 6 | | 405.6 | | |

Inter-day Precision

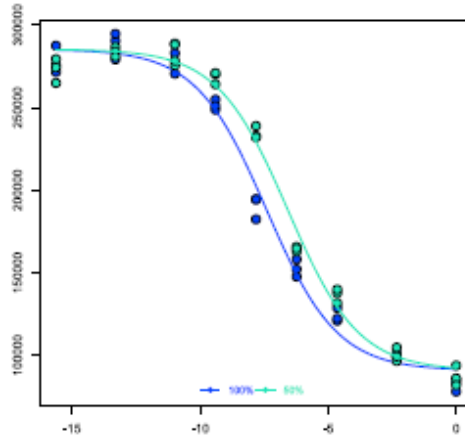
| Day | IC ₅₀ (ng/mL) | Mean IC ₅₀ (ng/mL) | %RSD |
|-----|--------------------------|-------------------------------|------|
| 1 | 453 | 474 | 24% |
| 2 | 597 | | |
| 3 | 372 | | |

Pertuzumab Bioassay Qualification with Perjeta®

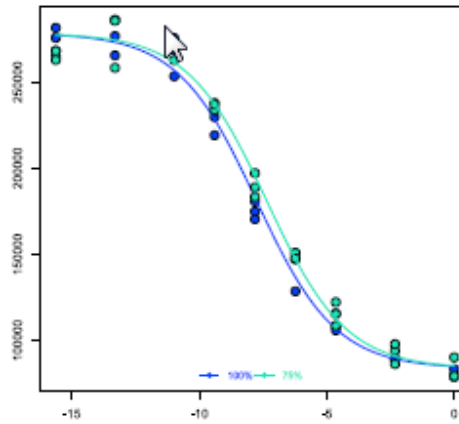
Parallelism and Potency Estimation (PLA)

Analyst 1

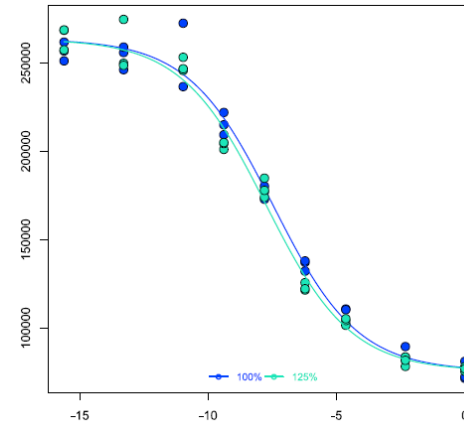
50%



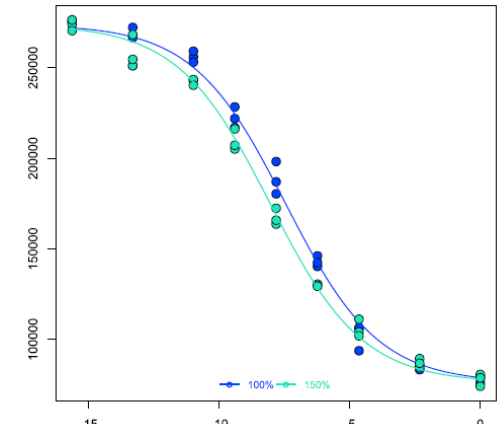
75%



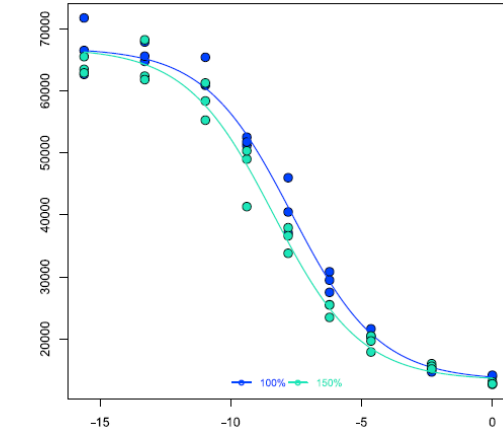
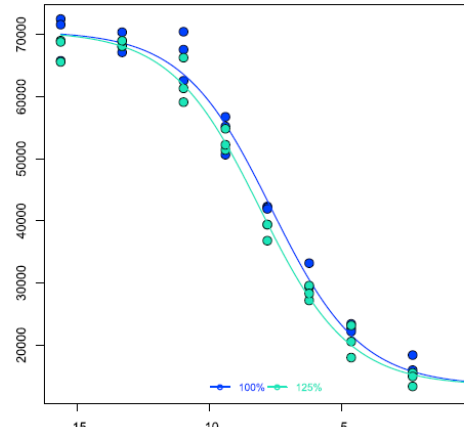
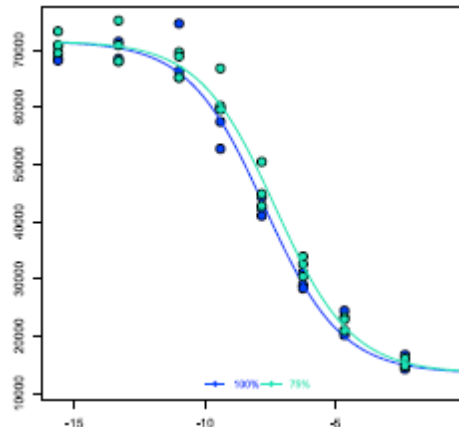
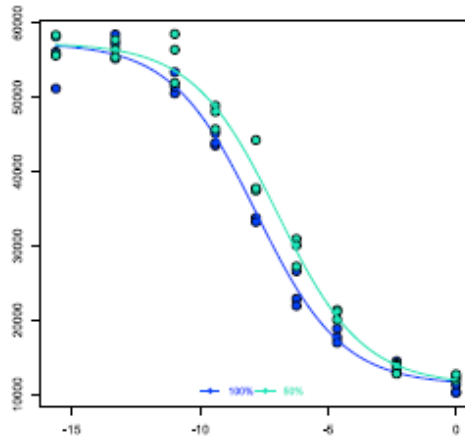
125%



150%



Analyst 2

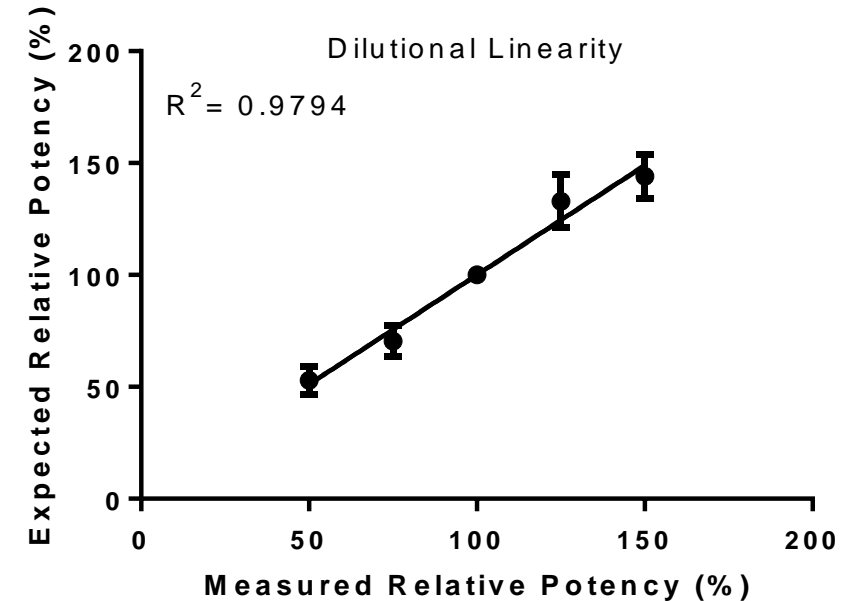


● Reference ● Sample

Pertuzumab Bioassay Qualification with Perjeta®

Summary: Accuracy, Precision and Dilutional Linearity

| Exp# | Expected RP (%) | Observed RP (%) | Average RP (%) | % RSD | % Recovery |
|------|-----------------|-----------------|----------------|-------|------------|
| 1 | 150 | 132.6 | 144 | 6.9 | 96 |
| 2 | | 148.9 | | | |
| 3 | | 150.6 | | | |
| 1 | 125 | 134.6 | 132 | 8.98 | 105.6 |
| 2 | | 120.3 | | | |
| 3 | | 144 | | | |
| 1 | 75 | 62.6 | 70.4 | 9.8 | 93.9 |
| 2 | | 75.8 | | | |
| 3 | | 72.8 | | | |
| 1 | 50 | 56.8 | 53 | 11.7 | 106 |
| 2 | | 45.8 | | | |
| 3 | | 56.3 | | | |



Accuracy: 100.4%

Precision: 9.35%

Benefits for “Ready-to-Use” Bioassay Kits

Functional response based on drug MOA

Verified and Qualified with innovator’s marketed drug

Simple protocol; Rapid results

Specific and Sensitive assay

Highly reproducible

Readily Implement with Optimized kit

- Frozen ready-to-assay cells
- Bioassay Detection Reagents
- Cell Plating Reagent
- Dilution Buffer
- Control Agonist
- Tissue Culture-Treated Plates

For More Info, Questions or Technical Support



Web:

[Cell-Based Bioassays for Biologics](#)

Technical Support

For NA:

DRX_SupportUS@eurofinsUS.com

For Europe, Africa & Middle East:

DRX_SupportEurope@eurofinsUS.com

For Asia-Pacific:

AsiaPacificSupport@eurofins.com