



# PathHunter<sup>®</sup> Adalimumab Bioassay

Qualified with Humira<sup>®</sup> (Adalimumab)

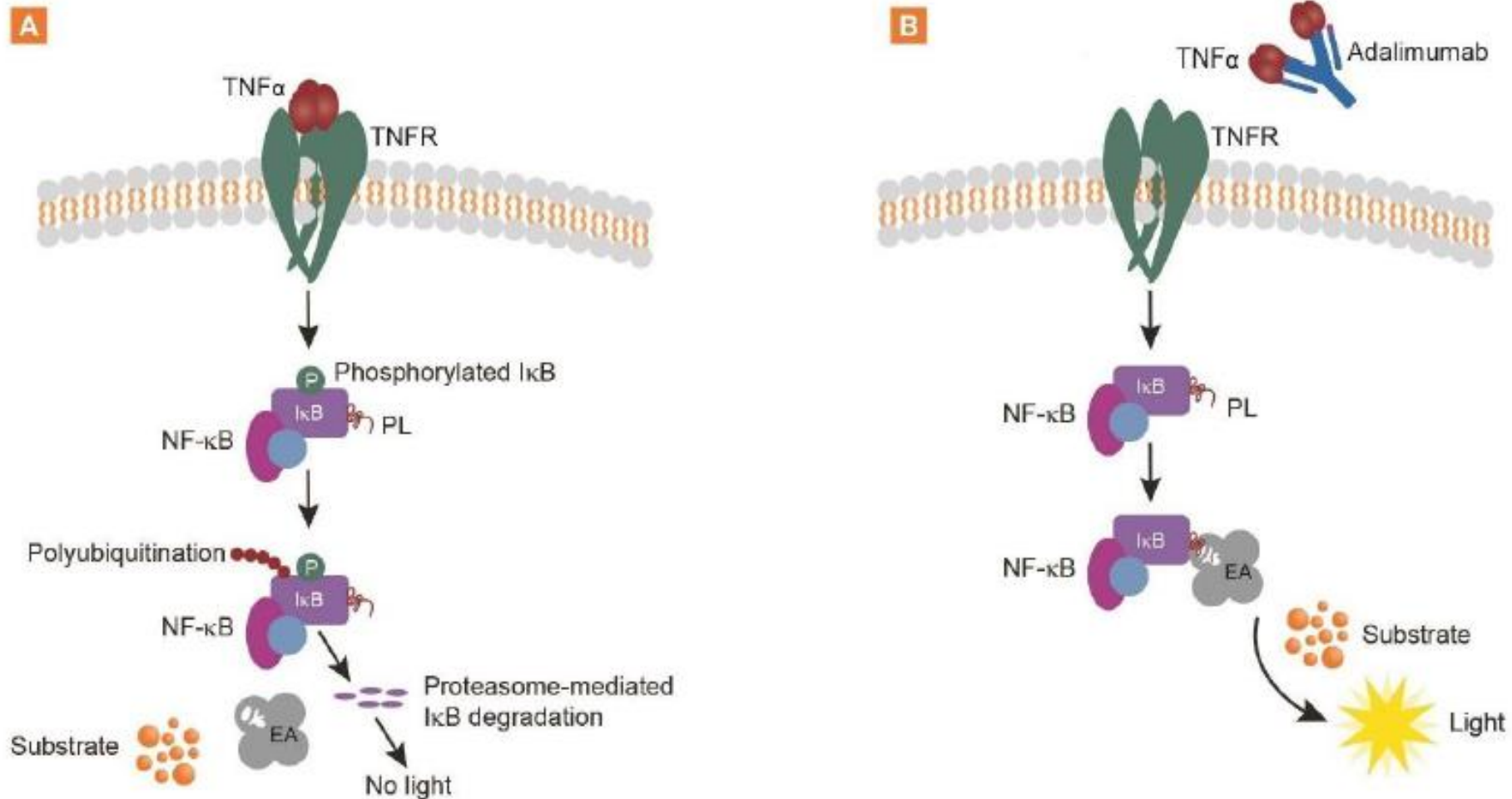
[93-0538B15-00131](#) (2-Plate Kit)

[93-0538B15-00132](#) (10-Plate Kit)

OUR EXPERTISE  
IN YOUR HANDS.  
DISCOVER  
CONFIDENTLY.

# Adalimumab-mediated I $\kappa$ B Degradation

## Assay Design



# PathHunter<sup>®</sup> Adalimumab Bioassay Kit

## Kit Components

### Materials Provided

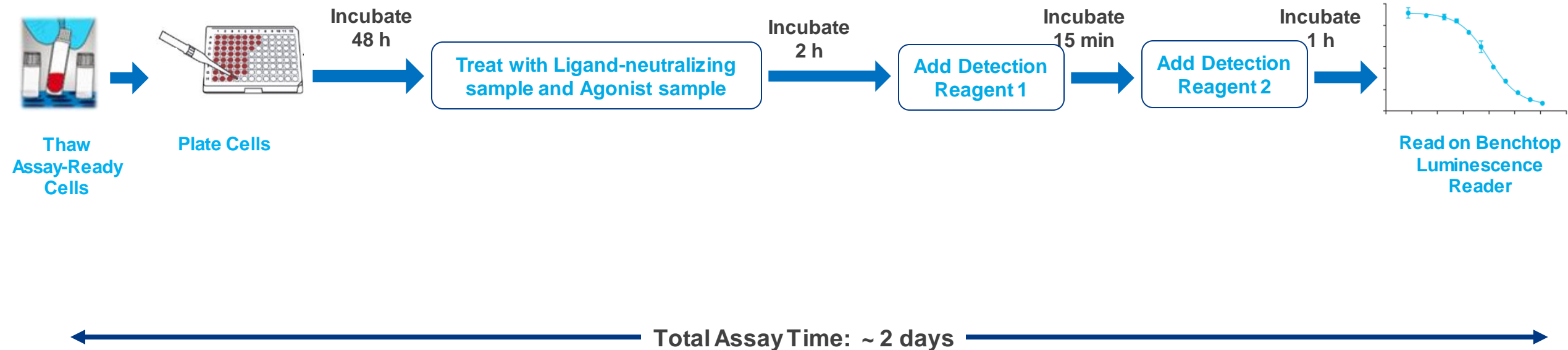
List of Components	93-0538B15-00131	93-0538B15-00132
PathHunter A549 IκB Bioassay Cells (1.2 x 10 <sup>6</sup> cells in 0.1 mL per vial)	2	10
AssayComplete™ Cell Plating 38 Reagent (100 mL per bottle)	1	2
Recombinant Human TNFα (10 μg per vial)	1	2
AssayComplete Protein Dilution Buffer (50 mL per bottle)	1	2
PathHunter Bioassay ED Detection Kit		
Detection Reagent 1 (Bottle)	1 x 6 mL	1 x 30 mL
Detection Reagent 2 (Bottle)	1 x 20 mL	1 x 100 mL
Detection Reagent 3 (Bottle)	1 x 5 mL	1 x 25 mL
96-Well White, Clear Flat-Bottom, TC-Treated, Sterile Plates with Lid	2	10

Sample Data

# Adalimumab Bioassay Kit Qualification

# Adalimumab Bioassay Workflow

*Simple, Homogenous and Rapid Protocol*



## *Assay Parameters Assessed*

- % CV between eight 11-pt DRCs
- Plate uniformity:  $EC_{80}$  and  $IC_{80}$  (of drug and stimulus) across entire plate
- Plate-to-Plate variability: 3 plates with 11-pt DRCs run on 3 days
- Slope consistency
- Accuracy, precision and linearity of the assay over a range of 50-150% from two operators
  - Assay developer
  - Assay qualifier
- Parallel line analysis

# Adalimumab Bioassay Qualification with Humira® Reproducibility Study (Over 3 Days)

## Full Plate

Experiment	Mean S/B	% RSD	Interday Mean S/B	%RSD	Mean IC <sub>50</sub> , ng/mL	% RSD	Interday Mean IC <sub>50</sub> , ng/mL	%RSD
Day 1	5.8	4.22	5.9	11.96	62.72	10.84	54.9	12.38
Day 2	5.3	3.03			50.52	6.09		
Day 3	6.7	8.10			51.44	4.21		

## Minus Rows A and H

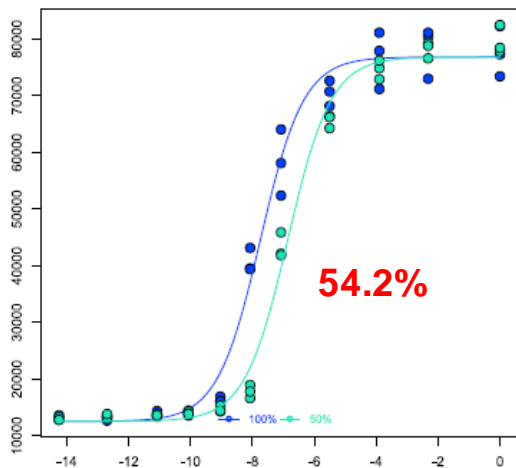
Experiment	Mean S/B	% RSD	Interday Mean S/B	%RSD	Mean IC <sub>50</sub> , ng/mL	% RSD	Interday Mean IC <sub>50</sub> , ng/mL	%RSD
Day 1	5.8	4.68	5.9	11.11	64.3	11.31	55.7	13.3
Day 2	5.3	3.01			51.2	3.77		
Day 3	6.6	3.39			51.7	2.81		

# Adalimumab Bioassay Qualification with Humira<sup>®</sup>

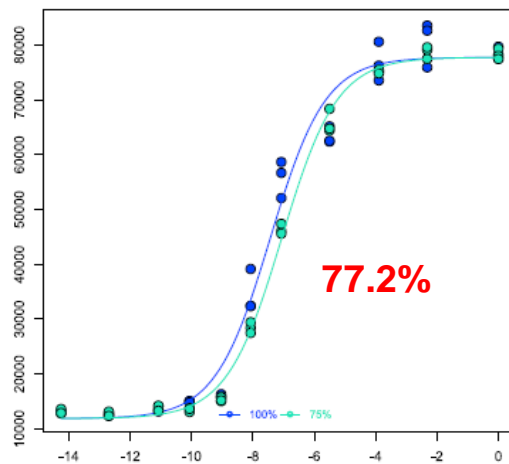
## Representative Relative Potency Data (50%-150%)

### Analyst 1

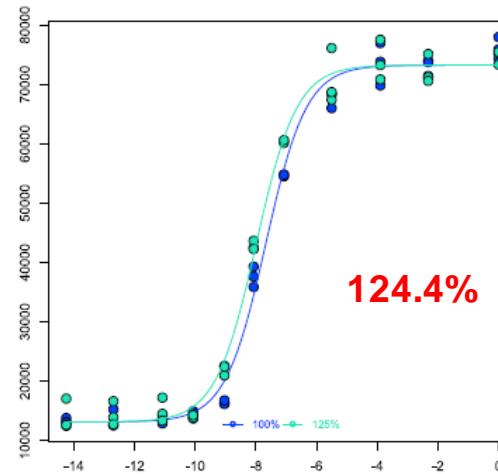
#### 100% vs 50%



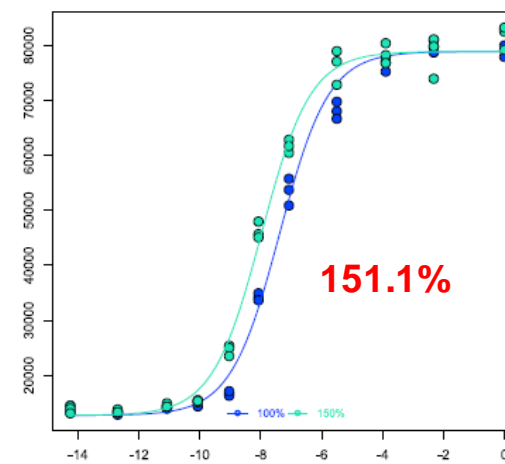
#### 100% vs 75%



#### 100% vs 125%

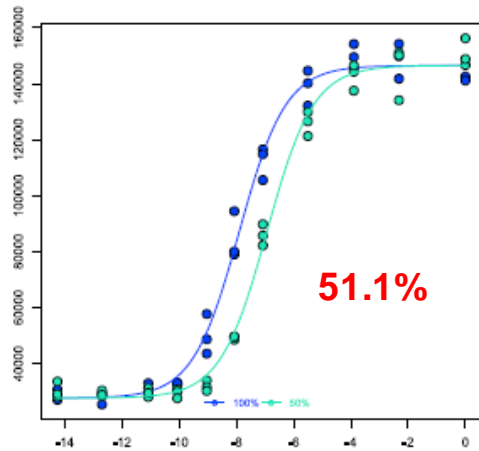


#### 100% vs 150%

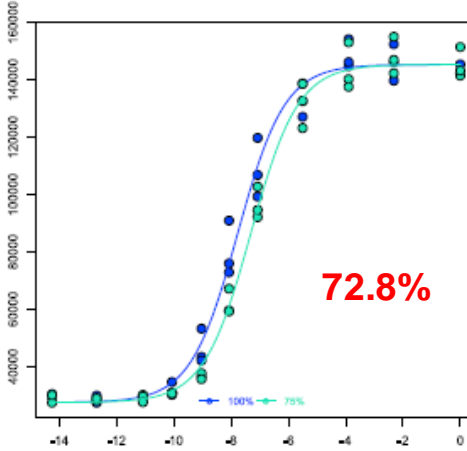


### Analyst 2

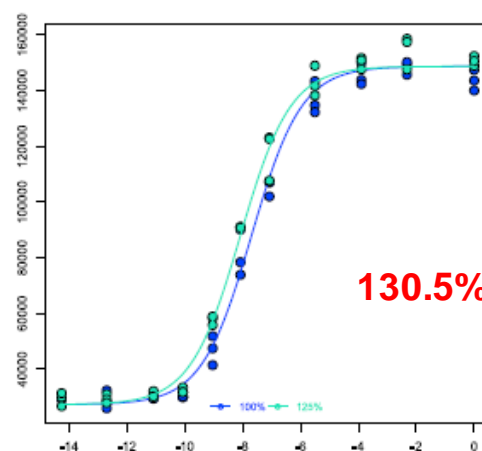
#### 100% vs 50%



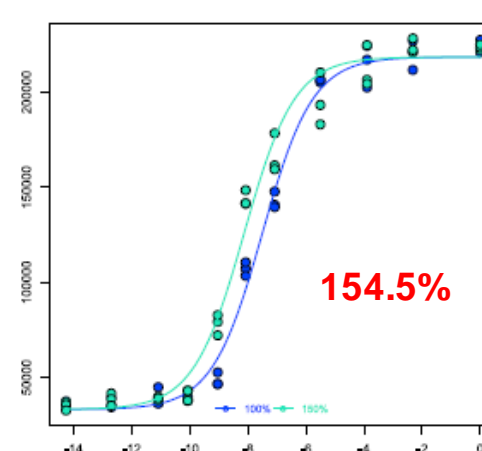
#### 100% vs 75%



#### 100% vs 125%



#### 100% vs 150%





# Adalimumab Bioassay Qualification with Humira®: Repeatability (Relative Potency)

The ICH Q2B, Section 5.1 recommends a minimum of 6 determinations at 100% nominal potency be used to assess repeatability. Repeatability was assessed using six determinations at 100% nominal relative potency by a single Analyst during the assessment of accuracy in the Adalimumab Bioassay.

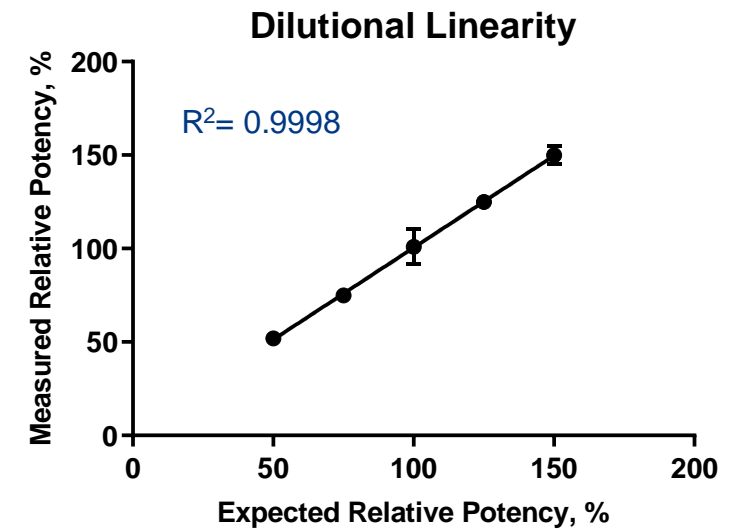
Analyst	Observed Potency at 100% Nominal Potency	Average Observed Potency	% RSD
A	88.3%	100.4%	8.44%
	94%		
	107%		
	104.8%		
	110.4%		
	97.9%		
<b>Repeatability Acceptance Criterion:</b>			≤ 20%

**Repeatability= 8.4%**

# Adalimumab Bioassay Qualification with Humira®: Accuracy, Intermediate Precision and Dilutional Linearity

Expected RP (%)	Exp #	Analyst #	Measured RP (%)	Average RP (%)	% RSD	% Recovery
150	1	1	150.6	149.9	3.26	99.9
	2	1	141.6			
	3	1	151.7			
	4	1	151.1			
	5	2	154.5			
125	1	1	126.4	124.9	3.06	99.9
	2	1	122.3			
	3	1	120.7			
	4	1	124.4			
	5	2	130.5			
75	1	1	80	75	4.71	100
	2	1	71.4			
	3	1	73.5			
	4	1	77.2			
	5	2	72.8			
50	1	1	55.5	51.9	5.65	103.8
	2	1	48.4			
	3	1	50.1			
	4	1	54.2			
	5	2	51.1			

*Intermediate precision addressed multiple analysts and multiple days*



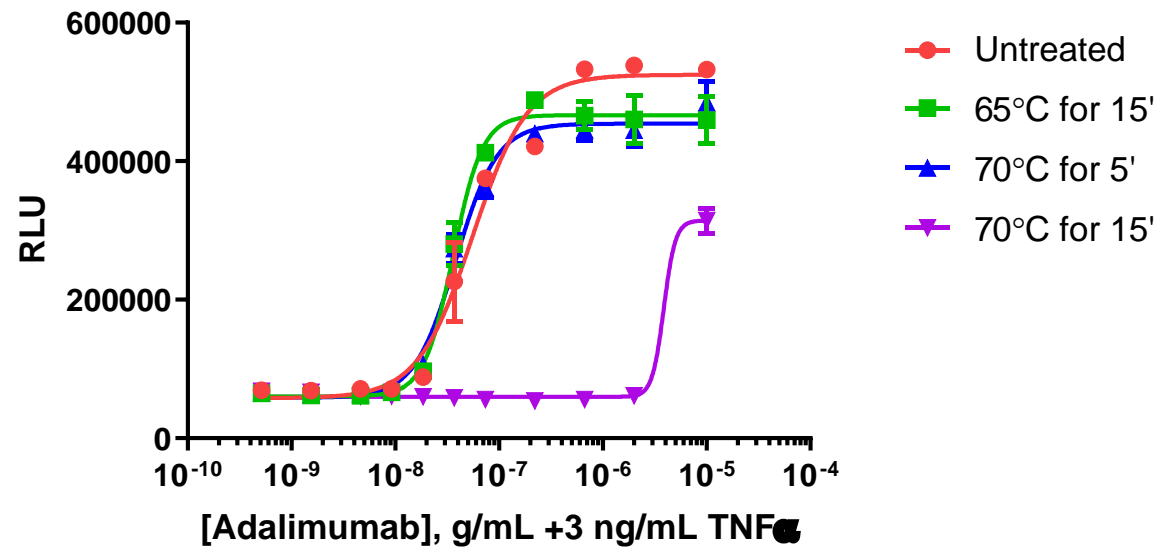
Accuracy = 100.9%  
 Intermediate precision  $\leq 5.65\%$   
 Linearity ( $R^2$ ) = 0.9998

# Adalimumab Bioassay Qualification with Humira® Evaluation of Forced Degradation Samples

## A549 $\alpha$ B Bioassay

P/N: 93-0538C15; L/N: 19B1807

48h recovery; 15' pre-incubation of drug +TNF $\alpha$  @ R/T;  
2h @ R/T drug incubation



*Treatment of Adalimumab at 70°C for 15 minutes produced a change in potency of the drug outside the range of the assay (right-shifted by two orders of magnitude).*

	Untreated	65C for 15'	70C for 5'	70C for 15'
Bottom	57786	60377	58409	59813
Top	524711	466425	454184	~ 314258
LogEC50	-7.256	-7.443	-7.419	~ -5.413
HillSlope	1.647	3.066	2.178	~ 7.765
EC50	5.550e-008	3.606e-008	3.810e-008	~ 3.866e-006

S/B            8.0            7.0            7.5            4.5

# For More Info, Questions or Technical Support



## **Web:**

[Cell-Based Bioassays for Biologics](#)

## **Technical Support**

For NA:

[DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

For Europe, Africa & Middle East:

[DRX\\_SupportEurope@eurofinsUS.com](mailto:DRX_SupportEurope@eurofinsUS.com)

For Asia-Pacific:

[AsiaPacificSupport@eurofins.com](mailto:AsiaPacificSupport@eurofins.com)