



LANCE *Ultra* STAT3 (Y705) Cellular Detection Kit

Product number: TRF4004 C/M

Research Use Only. Not for use in diagnostic procedures.

Product Information

Application:

This kit is designed for the detection of phosphorylated STAT3 (Y705) in cell lysates using a homogeneous LANCE *Ultra* assay (no wash steps).

Typical Performance

(Undiluted positive control lysate versus Buffer):

Signal/Background: 9.2

Z': 0.93

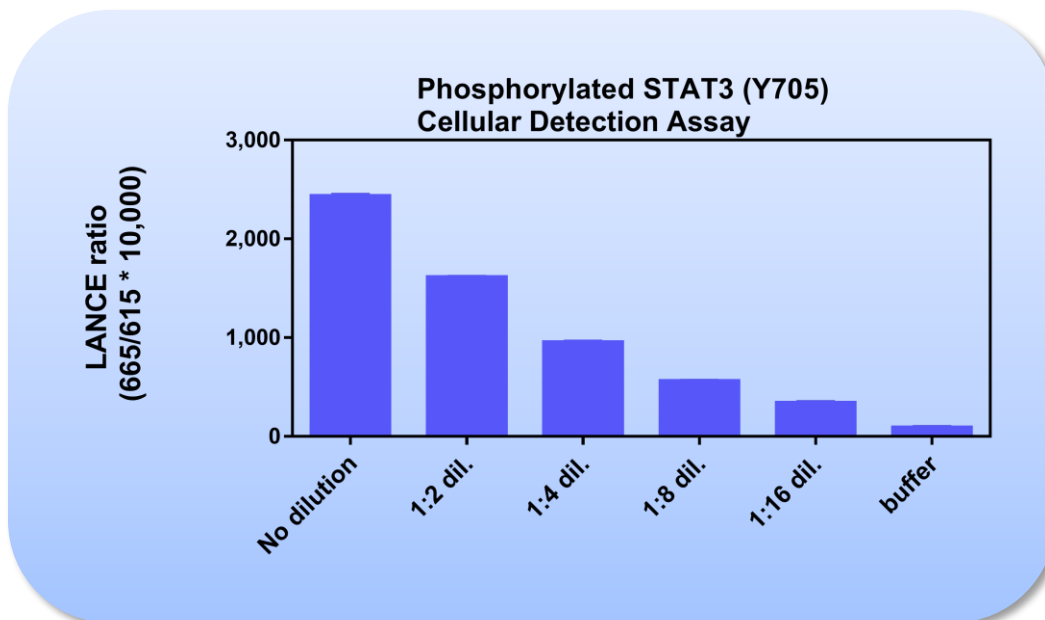


Figure 1. Typical positive control lysate (HeLa stimulated with IFN α) diluted in lysis buffer. The data was generated using a white Optiplate™-384 microplate and read on an EnVision™ Multilabel Plate Reader equipped with TR-FRET laser option. Total signal, signal/background window, and sensitivity may vary with other instruments. Positive control lysate is not supplied with the kit and must be purchased separately.

Storage:

Store kit in the dark at +4°C.

Stability:

This kit is stable for at least 6 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Analyte of Interest

STAT3 is an important transcriptional activator, and is involved in regulating cytokines and growth factor receptors. STAT3 is phosphorylated at Tyr-705 by the JAK pathway, upon which STAT3 will dimerize, translocate to the nucleus, and bind DNA. In some cancers STAT3 is shown to be constitutively activated. It has been implicated in both anti-apoptotic and oncogenic pathways.

Description of the LANCE *Ultra* Assay

LANCE® and LANCE® (Lanthanide chelate excite) *Ultra* are homogeneous (no wash) TR-FRET (time-resolved fluorescence resonance energy transfer) technologies. One antibody of interest is labeled with a donor fluorophore (a LANCE Europium chelate) and the second antibody is labeled with an acceptor fluorophore [ULight™ dye]. Upon excitation at 320 or 340 nm, energy can be transferred from the donor Europium chelate to the acceptor fluorophore if sufficiently close for FRET (~10 nm). This results in the emission of light at 665 nm. Data is represented as ratiometric (665/615 nm X 10,000).

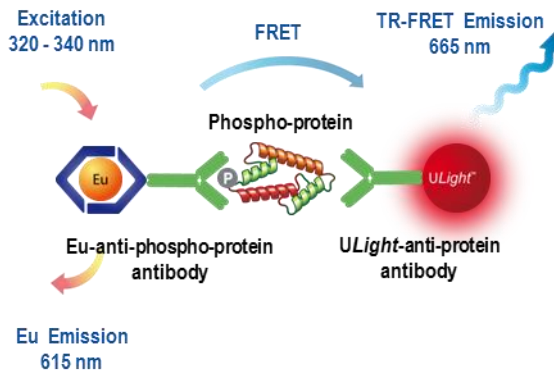


Figure 2. LANCE assay principle

Kit Content: Reagents and Materials

Kit components	TRF4004C (500 assay points**)	TRF4004M (10 000 assay points**)
LANCE <i>Ultra</i> Eu-labeled Anti-STAT3 (Y705) Antibody stored in TSA, 0.1% BSA	10 µL @ 500 nM (1 clear tube, yellow cap)	200 µL @ 500 nM (1 clear tube, yellow cap)
LANCE <i>Ultra ULight</i> -labeled Anti- STAT3 Antibody stored in TSA, 0.1% BSA	100 µL @ 500 nM (1 brown tube, blue cap)	2 X 1000 µL @ 500 nM (2 brown tubes, green caps)
LANCE Detection Buffer (10X) *	1.8 mL, 1 small bottle	250 mL, 1 large bottle
LANCE <i>Ultra</i> Lysis Buffer 1 (5X) *	10 mL, 1 small bottle	100 mL, 1 large bottle

* Extra detection buffer can be ordered separately (cat # CR97-100C: 1.8 mL or cat # CR97-100: 250 mL).
Extra Lysis Buffer can be ordered separately (cat # TRF001C: 10 mL or cat # TRF001F: 100 mL).

** The number of assay points is based on an assay volume of 20 µL in 384-well assay plates using the kit components at the recommended concentrations.

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the signal.

Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
VICTOR™ X, VICTOR Nivo™, ViewLux®, EnVision, EnSight™, EnSpire® Multilabel Plate Reader equipped with TR-FRET option	Revvity Inc.	Please consult our website-
TopSeal-A PLUS Adhesive Sealing Film	Revvity Inc.	6050185
Tissue culture treated clear SpectraPlates™, for culturing cells when using the 2-plate manual	Revvity Inc.	6005650
White OptiPlate-384, for LANCE <i>Ultra</i> detection assays when using the 2-plate manual	Revvity Inc.	6007290
White CulturPlate-384 when using the 1-plate manual	Revvity Inc.	6007680
Positive Control Cell Lysate (HeLa cells stimulated with IFN α)	Revvity Inc.	TRF4004S

Recommendations

General recommendations:

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube.
- Re-suspend all reagents by vortexing before use.
- Centrifuge all tubes before use to improve recovery of content (2000x g, 10-15 sec).
- Use Milli-Q® grade H₂O (18 M Ω ·cm) to dilute Detection and Lysis Buffers.
- When diluting the samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Films to reduce evaporation during incubation. LANCE *Ultra* TR-FRET assays cannot be read with the TopSeal-A Film attached. Please remove before reading.
- LANCE signal can be detected using a VICTOR X, ViewLux, EnVision, EnSpire, VICTOR NIVO, or EnSight Multilabel Reader equipped with TR-FRET. Use an excitation wavelength of 320 or 340 nm to excite the LANCE Europium chelate. We recommend you read this assay in dual emission mode, detecting both the emission from the Europium donor fluorophore at 615 nm, and the acceptor fluorophore (at 665 nm for *ULight* dye). The 665/615 nm x 10,000 calculation is used to process your data.
- Signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperatures should be used for each plate.
- The representative data shown in this technical data sheet are provided for information only.

Cell Handling and Lysis recommendations:

- Evaporation can be problematic with cells cultured in microtiter plates. For overnight incubation, it is recommended to add warm PBS or sterile water to unused wells. For longer incubation periods, a sterile breathable sealing membrane (Corning, cat. #3345) can be used to cover the plate. Alternatively, cells can be cultured in larger wells, and/or in a larger volume of culture medium.
- Phosphatase Inhibitors such as NaF and activated Na₃VO₄ can be added to lysis buffers to protect kinases without affecting LANCE detection.
- An incubation of 30 minutes is usually sufficient for cell lysis. However, the optimal lysis incubation time should be determined by each investigator using a time course study.
- A starving step with serum-free medium may be required depending on your target/cell line and should be evaluated in a separate experiment.
- For 2-plate manuals with adherent cells: cell seeding densities of 40K cells/well are usually sufficient for most cell lines. However, optimization of cell seeding density is recommended.
- For 1-plate manuals with suspension cells: Cell seeding densities of 100K cells/well are usually sufficient for most cell lines. However, optimization of cell seeding density is recommended.

Assay Procedure

Format	# of data points	Volume			Plate recommendation
		Final	Sample	Eu-Antibody/ <i>ULight</i> -Antibody MIX	
TRF4004C	250	40 µL	30 µL	10 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	20 µL	15 µL	5 µL	White OptiPlate-384 (cat # 6007290) White CulturPlate-384 (cat # 6007680)
	1 250	8 µL	6 µL	2 µL	ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290) White CulturPlate-384 (cat # 6007680)
	2 500	4 µL	3 µL	1 µL	White OptiPlate-1536 (cat # 6004290)
TRF4004M	10 000	20 µL	15 µL	5 µL	White OptiPlate-384 (cat # 6007290) White CulturPlate-384 (cat # 6007680)
	25 000	8 µL	6 µL	2 µL	ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290) White CulturPlate-384 (cat # 6007680)
	50 000	4 µL	3 µL	1 µL	White OptiPlate-1536 (cat # 6004290)

General Lysis Manual: Cells are lysed in 1X lysis buffer.

Each manual described below is designed for 500 assay points.

Lysate Preparation (2-plate manual for adherent cells):

Stimulation:

Plate 50 μ L of cells in 96-well tissue culture treated plate (allow adherence overnight in CO₂ atmosphere)

Add 50 μ L of 2X stimulator in cell culture media

Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Carefully remove cell supernatant and discard

Add 50 μ L of 1X lysis buffer (Recommended: LANCE *Ultra* Lysis Buffer 1)

Incubate while shaking plate at 350 rpm for 30 minutes at 23 °C

Transfer 15 μ L of lysate to 384-well plate

Inhibition:

Plate 50 μ L of cells in 96-well tissue culture treated plate (allow adherence overnight in CO₂ atmosphere)

Add 25 μ L of 3X inhibitor in cell culture media

Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Add 25 μ L of 4X stimulator in cell culture media

Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Carefully remove cell supernatant and discard

Add 50 μ L of 1X lysis buffer (Recommended: LANCE *Ultra* Lysis Buffer 1)

Incubate while shaking plate at 350 rpm for 30 minutes at 23 °C

Transfer 15 μ L of lysate to 384-well plate

Lysate Preparation (1-plate manual for suspension cells):

Stimulation:

Plate 8 μ L of cells in 384-well OptiPlate

Add 4 μ L of 3X stimulator in cell culture media

Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Add 3 μ L of 5X lysis buffer
(Recommended: LANCE *Ultra* Lysis Buffer 1)

Incubate while shaking plate at 350 rpm for 30 minutes at 23 °C

Inhibition:

Plate 8 μ L of cells in 384-well Optiplate

Add 2 μ L of 5X inhibitor in cell culture media

Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Add 2 μ L of 6X stimulator in cell culture media

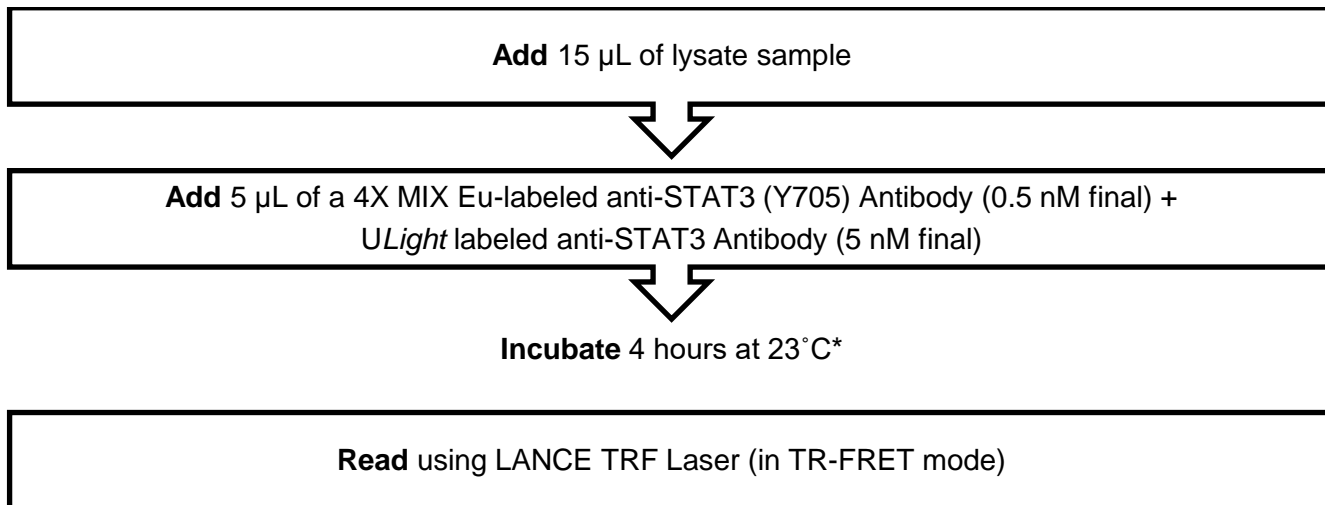
Incubate while shaking at 350 rpm for 30 minutes at 23 °C

Add 3 μ L of 5X lysis buffer
(Recommended: LANCE *Ultra* Lysis Buffer 1)

Incubate while shaking plate at 350 rpm for 30 minutes at 23 °C

Reagent Preparation:

- 1) Preparation of 1X LANCE Detection Buffer:
 - a. Add 1 mL of 10X LANCE Detection Buffer to 9 mL H₂O.
- 2) Preparation of 4X MIX Eu-labeled anti-STAT3 (Y705) Antibody (2 nM) + ULight labeled anti-STAT3 Antibody (20 nM):
 - a. Prepare just before use.
 - b. Add 10 μ L of 500 nM Eu-labeled anti-STAT3 (Y705) Antibody and 100 μ L of 500 nM ULight-labeled anti-STAT3 Antibody to 2390 μ L of LANCE Detection Buffer
- 3) In a white Optiplate (384 wells):



***In order to reduce evaporation, we recommend covering the OptiPlate with TopSeal-A PLUS during the incubation. Longer incubation times can be used and in some cases may improve assay signal/background.**

Important: LANCE signal is detected using an EnVision Multilabel Reader equipped with a TR-FRET laser. Use an excitation wavelength of 320 or 340 nm to excite the LANCE Europium chelate. We recommend you read this assay in dual emission mode, detecting both the emission from the Europium donor fluorophore at 615 nm, and the acceptor fluorophore (at 665 nm for ULight dye). Data is calculated and presented ratiometrically by dividing the signal at 665 nm by the signal at 615 nm and multiplying by 10,000.

Data Analysis

- Data is represented ratiometrically. Divide the signal at 665 nm by the signal at 615 nm and multiply by 10,000.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a 1/Y² data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).

Assay Performance Characteristics

LANCE Ultra assay performance described below was determined using the 2-plate manual.

Dose Response Curve:

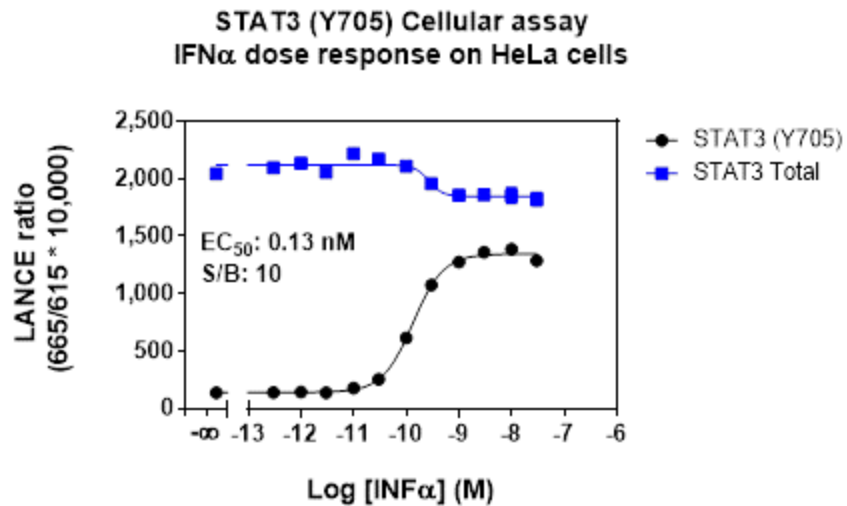


Figure 3: Dose Response Curve HeLa cells stimulated with IFN α . HeLa cells (40K/well) were treated with increasing concentrations of IFN α for 30 minutes prior to lysis with LANCE Ultra Lysis Buffer 1 for 30 minutes at room temperature. TRF4005 was used to detect Total STAT3.

Inhibition Dose Response:

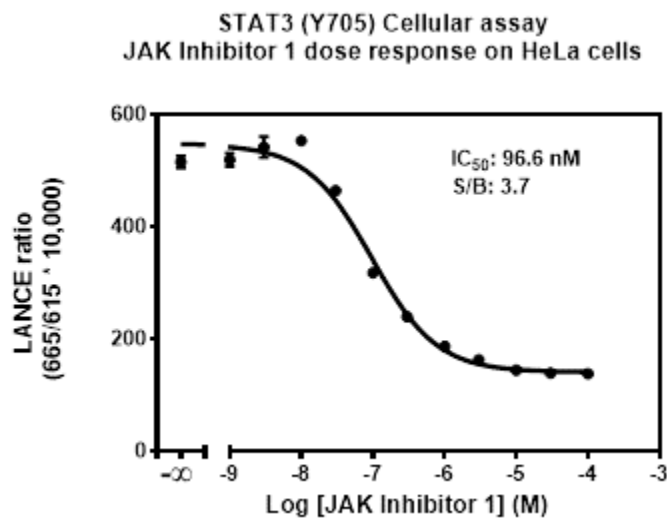


Figure 4: Inhibition Dose Response Curve HeLa cells treated with JAK1 and stimulated with IFN α . HeLa cells (40K/well) were treated with increasing concentrations of JAK1 and incubated for 30 minutes. The cells were then stimulated with 1 nM IFN α for 30 minutes prior to lysis with LANCE Ultra Lysis Buffer 1 for 30 minutes at room temperature.

Assay Robustness (Z')

The assay was tested with Stimulated and Unstimulated cell lysates in multiple replicates (n=36) and a Z' score was determined.

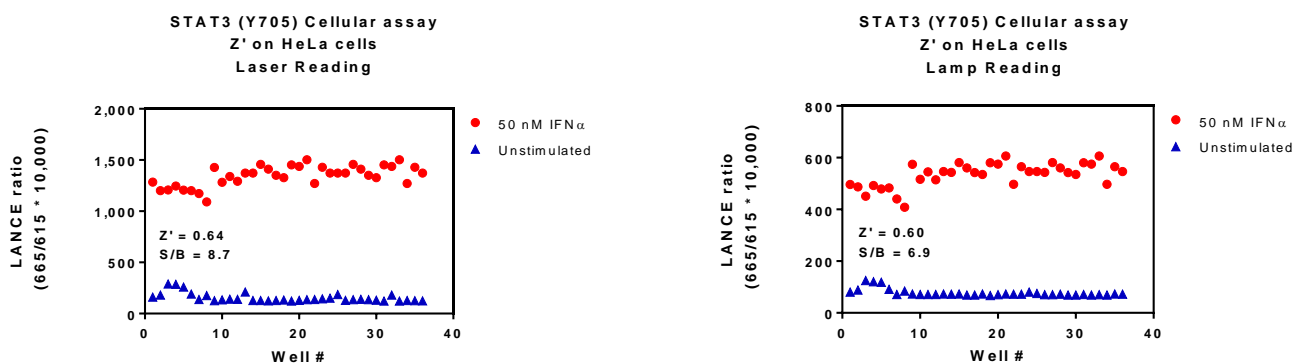


Figure 5: Z' determination in Stimulated versus Unstimulated cell lysates read using laser or lamp configuration. HeLa cells (40K/well) were treated with 50 nM IFN α for 30 minutes prior to lysis with LANCE *Ultra* Lysis Buffer 1 for 30 minutes at room temperature. The same plate was read using both the TR-FRET laser or lamp options.

Additional Resources

For more information regarding Lance Ultra Assays follow the link below:

www.revivity.com

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.