



## MDA5 TOTAL KITS

**Part # 63ADK112PEG & 63ADK112PEH**

**Test size#:** 500 tests (63ADK112PEG), 10,000 tests (63ADK112PEH) - assay volume: 20  $\mu$ L

**Revision:** #04 of March 2024

**Store at:**  $\leq$ - 60°C (63ADK112PEG);  $\leq$ - 60°C (63ADK112PEH)

**For research use only. Not for use in diagnostic procedures.**

### ASSAY PRINCIPLE

Revvity MDA5 total assay is only intended for quantitative measurement of MDA5 total in cells using HTRF<sup>®</sup> technology.

MDA5 total is detected in a sandwich assay format using 2 different specific antibodies, labeled with Europium Cryptate (donor) and with d2 (acceptor).

The principle of detection is based on HTRF<sup>®</sup> technology. When the dyes are in close proximity, the excitation of the donor with a light source (laser or flash lamp) triggers a Fluorescence Resonance Energy Transfer (FRET) towards the acceptor, which in turn fluoresces at a specific wavelength (665 nm). The donor & acceptor labeled antibodies bind to the MDA5 present in the sample, thereby generating FRET. Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the MDA5 total concentration (Fig. 1).

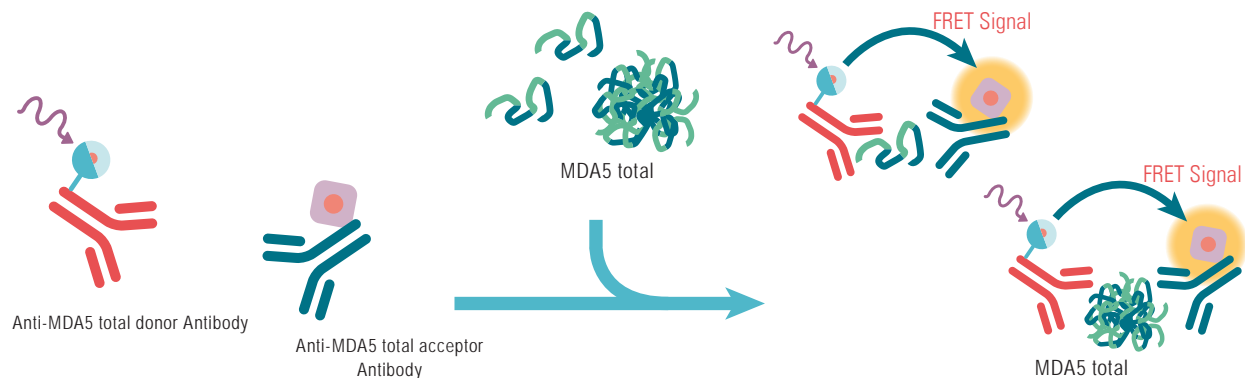
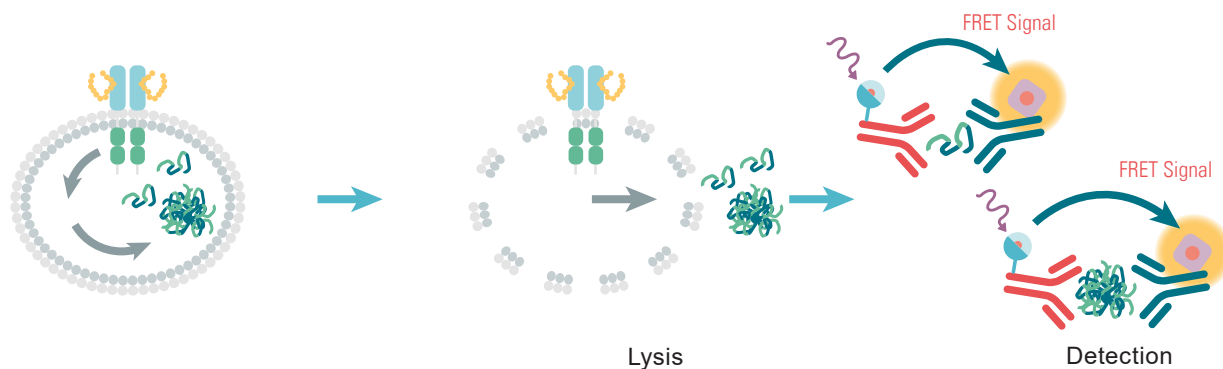


Figure 1: Principle of HTRF MDA5 total sandwich assay.

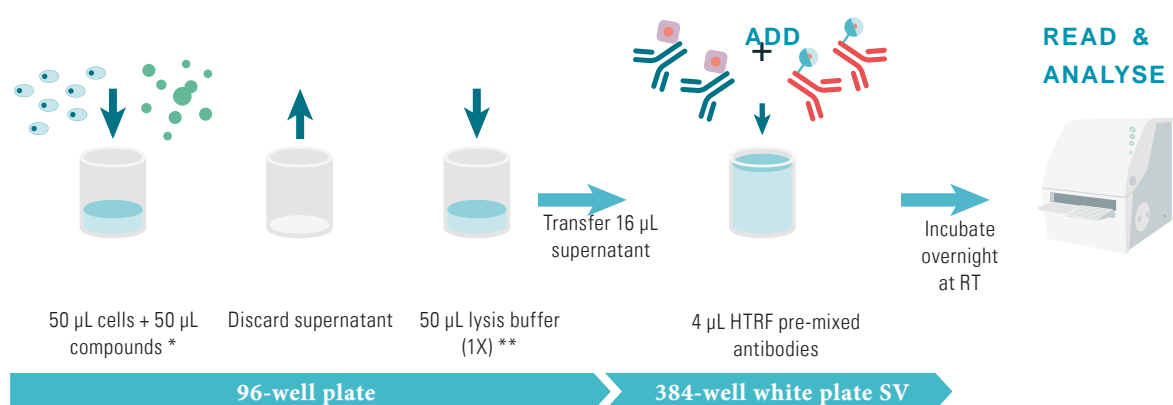
The assay is run under a two-plate assay manual, where cells are plated, stimulated and lysed in the same culture plate. Lysates are then transferred to the assay plate for the detection of MDA5 total by HTRF<sup>®</sup> reagents. This manual gives the cells viability and confluence to be monitored.

Technical support team can help you to set-up this manual or another one.  
Please contact us at [www.revvity.com](http://www.revvity.com)

## MANUAL AT A GLANCE



### TWO-PLATE ASSAY MANUAL (FOR ADHERENT CELLS):



\* Note that concentration above 0.5% DMSO will impair assay performances.

\*\* Depending on cell lines used, volume of lysis should be optimized, it can also be necessary to dilute the cell lysate to ensure samples are within the assay linear range.

### MATERIALS PROVIDED:

Kit components	500 tests Cat # 63ADK112PEG	10,000 tests Cat # 63ADK112PEH
Control lysate Frozen/ready-to-use	1 vial - 150 µL	2 vials - 150µL
Anti-MDA5 total-Eu Cryptate Antibody	1 vial - 20 µL Frozen - 50 X	1 vial - 400µL Frozen - 50 X
Anti-MDA5 total-d2 Antibody	1 vial - 20 µL Frozen - 50 X	1 vial - 400 µL Frozen - 50 X
Lysis buffer #3 * stock solution 4X	4 vials - 2 mL Frozen	1 vial - 130 mL Frozen
Detection Buffer #8 ** ready-to-use	1 vial - 2 mL Frozen	1 vial - 50 mL Frozen

\* Amounts of reagents provided are sufficient for generating 50 µL of cell lysate per well.

\*\* The Detection Buffer is used to prepare working solutions of acceptor and donor reagents.

### PURCHASE SEPARATELY:

- HTRF®-Certified Reader\*\*. Make sure the setup for Eu Cryptate is used
- For a list of HTRF-compatible readers and set-up recommendations, please visit [www.revvy.com](http://www.revvy.com)
- Small volume (SV) detection microplates - Use white plate only.
- For more information about microplate recommendations, please visit our website at: [www.revvy.com](http://www.revvy.com)

## STORAGE AND STABILITY

Antibodies, control lysate and buffers should be stored frozen until use.

Thawed detection buffer can be stored at 2-8°C in your premises. Thawed antibodies are stable 48 hours at 2-8°C; they can be refrozen (at -20°C or below) and thawed at least one more time. Control lysate must be stored frozen at -60°C or below. Thawed control lysate can be refrozen (at -60°C or below) and thawed one more time.

## REAGENT PREPARATION

Allow all reagents to thaw before use.

We recommend centrifuging the vials gently after thawing, before pipeting the stock solutions.

Prepare the working solutions from stock solutions by following the instructions below.

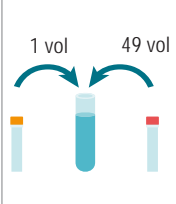
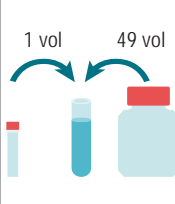
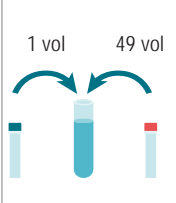
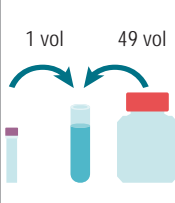
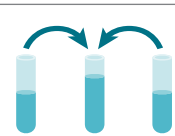
### POSITIVE CONTROL SOLUTION: READY-TO-USE

The control cell lysate is only provided as an internal assay control to check the quality of the results obtained. The window between control lysate and negative control should be greater than 2.

### TO PREPARE WORKING ANTIBODY SOLUTIONS:

HTRF® reagent concentrations have been set for optimal assay performances. Note that any dilution or improper use of the d2 and Cryptate-antibodies will impair the assay's quality. Be careful, as working solution preparation for antibodies may differ between the 500 and 10,000 tests data point kit.

Antibody working solutions are stable for 2 days at 4°C. Dilute the antibodies with detection buffer #8.

500 TESTS KIT - 63ADK112PEG			10,000 TESTS KIT - 63ADK112PEH
<b>Anti-MDA5 total- Cryptate antibody</b>			
Dilute 50-fold the frozen stock solution with Detection buffer #8: e.g. add 2.45 mL of detection buffer to the 0.05 mL of Cryptate-antibody stock solution.			Dilute 50-fold the frozen stock solution with Detection buffer #8: e.g. add 49 mL of detection buffer to the 1 mL of Cryptate antibody stock solution.
<b>Anti-MDA5 total-d2 antibody</b>			
Dilute 50-fold the frozen stock solution with Detection buffer #8: e.g. add 2.45 mL of detection buffer to the 0.05 mL of d2- antibody stock solution.			Dilute 50-fold the frozen stock solution with Detection buffer #8: e.g. add 49 mL of detection buffer to the 1 mL of d2- antibody stock solution.
<b>Antibody mix</b>			
It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2-antibody solution to 1 volume of Cryptate-antibody solution.			It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2-antibody solution to 1 volume of Cryptate-antibody solution.

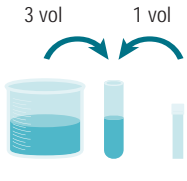

## TO PREPARE LYSIS BUFFER:

Make sure that the lysate has been generated by using the kit reagents.

Prepare the required amount of lysis buffer before running the assay, working solutions are stable for 2 days at 2-8°C.

Lysis buffer 1X:

Determine the amount of lysis buffer needed for the experiment. Each well requires generally 50 µL of lysis buffer. Prepare a lysis buffer solution 1X by diluting 4-fold the lysis buffer 4X with distilled water.

500 TESTS KIT - 63ADK112PEG	Preparation of lysis buffer 1X		10,000 TESTS KIT - 63ADK112PEH
Dilute the "lysis buffer 4X" 4-fold with distilled water to prepare lysis buffer 1X. e.g. take 1.25 mL of lysis buffer 4X and add it to 3.75 mL of distilled water. Mix gently.			Dilute the "lysis buffer 4X" 4-fold with distilled water to prepare lysis buffer 1X. e.g. take 1.25 mL of lysis buffer 4X and add it to 3.75 mL of distilled water. Mix gently.

## TWO PLATE ASSAY MANUAL

### FOR ADHERENT CELLS

#### GENERAL LAB WORK PRIOR USING REVVITY KIT: CELLS PREPARATION

Plate 50 µL of cells in 96-well tissue-culture treated plate in appropriate growth medium and incubate overnight, at 37°C in CO<sub>2</sub> atmosphere.

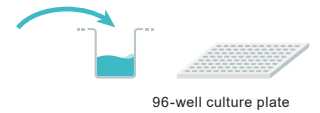
- Cell seeding densities of 200 K cells/well are generally sufficient for most cell lines, but optimization of cell seeding densities is recommended. Depending on receptor a starving step with serum-free medium could be essential.

Dispense 50 µL of compounds (2X) diluted in cell culture medium

- For most compound, incubation time should be above 24 hours at 37°C. We recommend a time course study to determine the optimal stimulation time. Note that concentration above 0.5% DMSO will impair assay performances. Same final concentration of DMSO must be used for each compound dilutions.



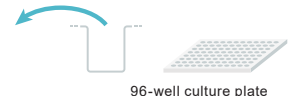
96-well culture plate



96-well culture plate

- Remove carefully cell supernatant either by aspirating supernatant or by flicking the plate.

Discard supernatant (for adherent cells)

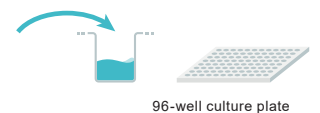


96-well culture plate

#### MDA5 TOTAL DETECTION USING REVVITY KIT

Immediately add 50 µL of lysis buffer #3 (1X) and incubate for at least 30 minutes at room temperature under shaking.

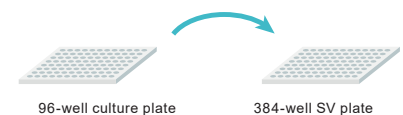
- Use the appropriate lysis buffer and incubate at room temperature with shaking. We recommend a time course study to determine the optimal lysis incubation time. Lysis volume can be decreased down to 25 µL.



96-well culture plate

- After homogenization by pipeting up and down, transfer 16 µL of cell lysate from the 96-well cell-culture plate to a 384-well small volume white plate.

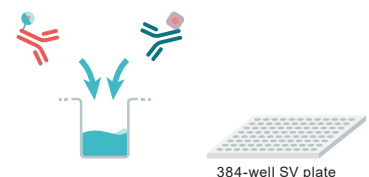
Depending on cell lines used, it can be necessary to dilute the cell lysate to ensure samples are within the assay linear range



96-well culture plate

384-well SV plate






- Add 4 µL of premixed antibody solutions (vol/vol) prepared in the detection buffer #8. Cover the plate with a plate sealer. Incubate overnight at RT. Set up your reader for Eu Cryptate and read the fluorescence emission at two different wavelengths (665nm and 620nm) on a compatible HTRF® reader\*.



384-well SV plate

\* For more information about HTRF® compatible readers and for set-up recommendations, please visit our website at: [www.revvity.com](http://www.revvity.com)

Standard manual for two-plate assay manual in 20 µL final volume (after lysis step)

	Non treated cell lysate	Treated cell lysate	Positive control	Negative control	Blank control
<b>Step 1</b> 	Dispense 16 µL of non treated cell lysate	Dispense 16 µL of treated cell lysate	Dispense 16 µL of control lysate	Dispense 16 µL of lysis buffer 1X	Dispense 16 µL of non treated cell lysate
<b>Step 2</b> 	Add 2 µL of Anti-MDA5 total-d2 Antibody working solution to all wells				Add 2 µL of detection buffer
<b>Step 3</b> 	Add 2 µL of Anti MDA5 total-Eu Cryptate Antibody working solution to all wells				
<b>Step 4</b> 	Cover the plate with a plate sealer. Incubate overnight at room temperature.				
<b>Step 5</b> 	Remove the plate sealer and read on an HTRF® compatible reader				

The blank control is used to check the Cryptate signal at 620 nm.

The Negative control is used to check the non-specific signal. The ratio between control lysate signal / non-specific signal should be greater than 2.

## DATA REDUCTION & INTERPRETATION

1. Calculate the ratio of the acceptor and donor emission signals for each individual well.

$$\text{Ratio} = \frac{\text{Signal 665 nm}}{\text{Signal 620 nm}} \times 10^4$$

2. Calculate the % CVs. The mean and standard deviation can then be worked out from ratio replicates.

$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

3. Calculate the % delta F which reflects the signal to background of the assay. The negative control plays the role of an internal assay control. Delta F is used for the comparison of day to day runs of the same assay.

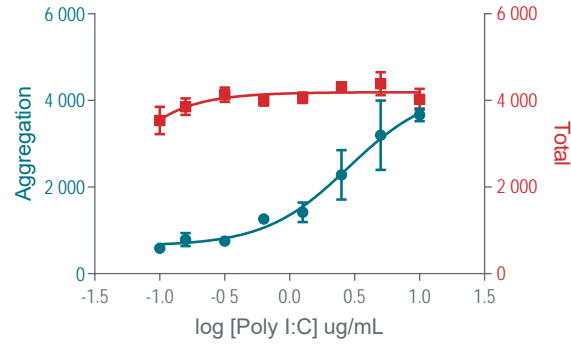
$$\text{delta F (\%)} = \frac{\text{Ratio Standard or sample} - \text{Ratio Negative Control}}{\text{Ratio Negative Control}} \times 100$$

For more information about data reduction, please visit [www.revivity.com](http://www.revivity.com)

## RESULTS

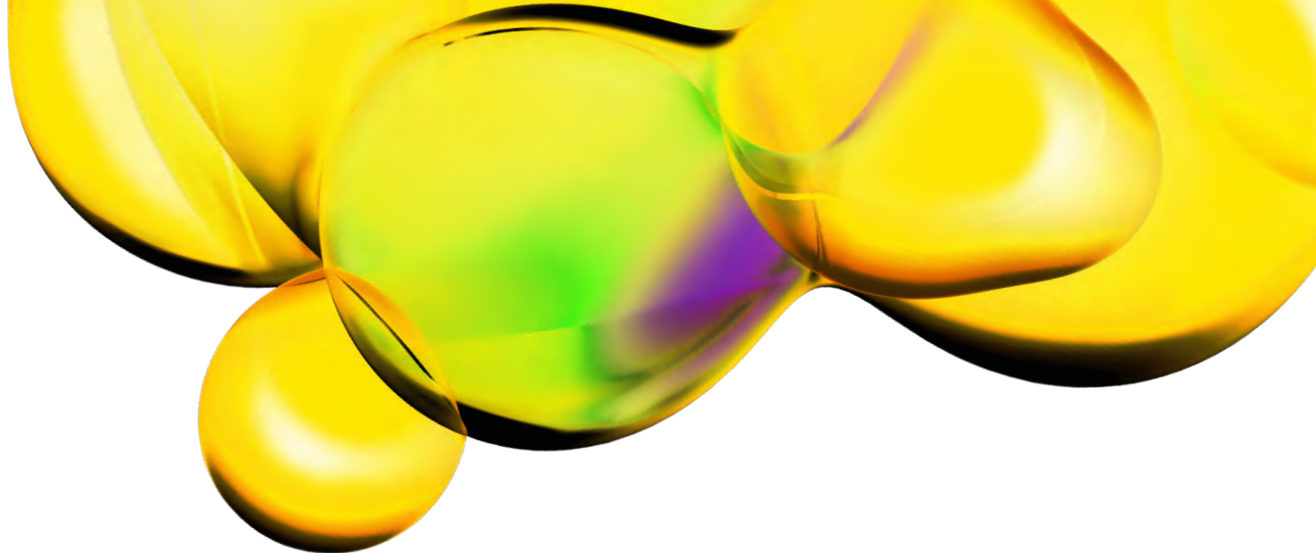
These data should be considered only as an example. Results may vary from one HTRF® compatible reader to another. PHERAstarFS with flash lamp (BMG) was used for reading.

MDA5 Aggregation and Total MDA5 assays



REACH European regulations and compliance This product and/or some of its components include a Triton concentration of 0.1% or more and as such, it is concerned by the REACH European regulations. We recommend researchers using this product to act in compliance with REACH and in particular: to only use the product for in vitro research in appropriate and controlled premises by qualified researchers, ii) to ensure the collection and the treatment of subsequent waste, and iii) to make sure that the total amount of Triton handled does not exceed 1 ton per year.

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage. The use of the cell line will be done with appropriate safety and handling precautions to minimize health and environmental impact. Remaining disclaimer.



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](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.