



# PROGRANULIN KITS

## PROTOCOL

### Part # 62PROPEG & 62PROPEH

**Test size#:** 500 tests (62PROPEG) and 10,000 tests (62PROPEH) - assay volume: 20  $\mu$ L

**Revision:** 03-May 2020

**Store at:** -60°C or below (62PROPEG); -60°C or below (62PROPEH)

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

### ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of Progranulin in cell/tissue culture supernatants and offers a fast alternative to ELISA.

The detection principle of this kit is based on HTRF® technology (Homogeneous Time-Resolved Fluorescence). As shown in Figure 1, Progranulin is detected in a sandwich assay by using anti Progranulin antibody labeled with Europium cryptate (donor), and anti Progranulin antibody labeled with d2 (acceptor).

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). Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the progranulin concentration.

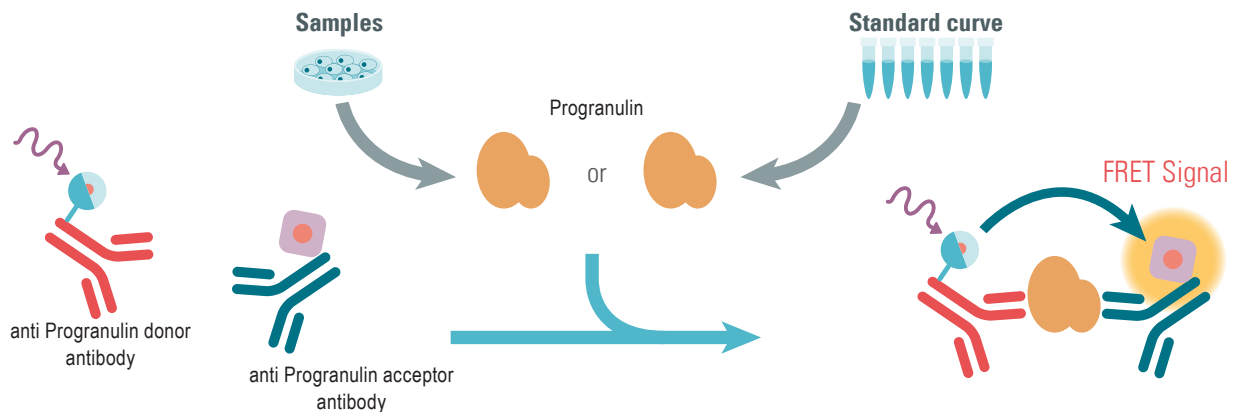
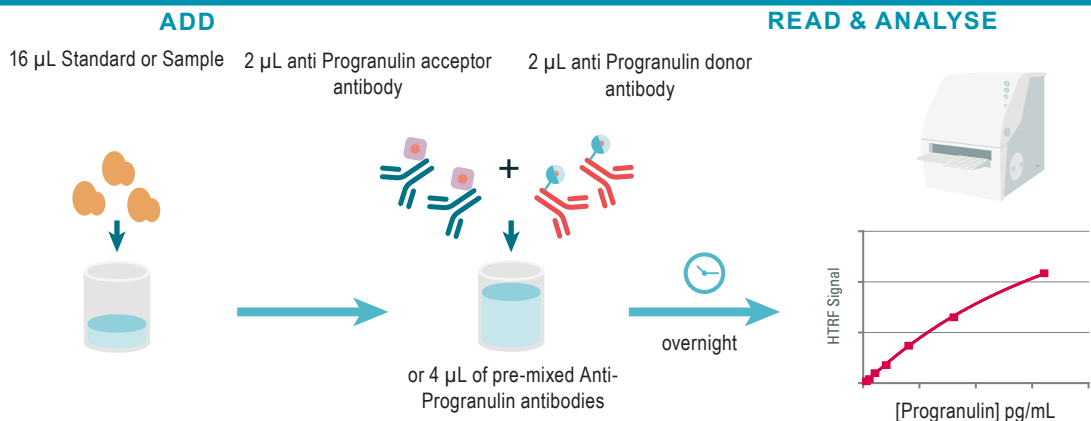


Figure 1: Principle of HTRF Progranulin sandwich assay.

### PROTOCOL AT A GLANCE



Make sure to use the set-up for Eu Cryptate.

**MATERIALS PROVIDED:**

KIT COMPONENTS	500 TESTS * CAT # 62PROPEG	10,000 TESTS * CAT # 62PROPEH
Progranulin Standard Frozen	1 vial - 50 µL 1 µg/mL	1 vial - 50 µL 1 µg/mL
Progranulin Eu Cryptate Antibody	1 vial - 20 µL Frozen - 50X	1 vial - 0.4 mL Frozen - 50X
Progranulin d2 Antibody	1 vial - 20 µL Frozen - 50X	1 vial - 0.4 mL Frozen - 50X
Diluent #5 ** 5X	1 vial 2 mL	1 vial 10 mL
Detection buffer *** ready-to-use	2 vials 1.5 mL Detection Buffer #3	1 vial 50 mL Detection Buffer #3

\* When used as advised, the two available kit sizes will provide sufficient reagents for 500 tests and 10,000 tests respectively in 20 µL final volume..

Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

\*\* Medium like cell culture medium can be an alternative to the diluent.

\*\*\* 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.cisbio.com/compatible-readers](http://www.cisbio.com/compatible-readers)

- Small volume (SV) detection microplates - Use white plate only.

For more information about microplate recommendations, please visit our website at: [cisbio.com/microplates-recommendations](http://cisbio.com/microplates-recommendations)

**STORAGE AND STABILITY**

Store the kit at -60°C or below.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label. Diluent and detection buffer are shipped frozen, but can be stored at 2-8°C in your premises.



Reagents

If lyophilized, reconstituted reagents, antibodies, and standard stock solutions may be frozen and thawed only once. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below .







Volume of Progranulin standard aliquots should not be under 10 µL.

**REAGENT PREPARATION****BEFORE YOU BEGIN:**

- It is very important to prepare reagents in the specified buffers. The use of an incorrect diluent may affect reagent stability and assay results.
- Thaw the frozen reagents at room temperature, allow them to warm up to room temperature for at least 30 mins before use
- Before use, allow Diluent and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- Progranulin standards (for standard curve) must be prepared in diluent or in the same medium as the samples.

**TAKE CARE TO PREPARE STOCK AND WORKING SOLUTIONS ACCORDING TO THE DIRECTIONS FOR THE KIT SIZE YOU HAVE PURCHASED.**

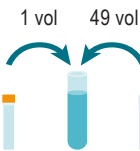
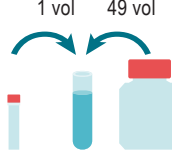
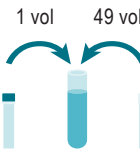
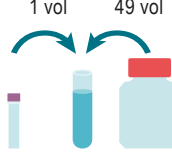

## TO PREPARE REAGENT STOCK SOLUTIONS:

500 TESTS KIT - 62PROPEG		10,000 TESTS KIT - 62PROPEH	
Anti-Progranulin Eu Cryptate antibody			
Thaw the Progranulin Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.			Thaw the Progranulin Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.
Anti-Progranulin d2 antibody			
Thaw the Progranulin d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.			Thaw the Progranulin d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below.
Progranulin Standard			
Thaw the Progranulin standard solution in order to obtain a 1 µg/mL stock solution. Mix gently.			Thaw the Progranulin standard solution in order to obtain a 1 µg/mL stock solution. Mix gently.
Diluent			
Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 1 mL of diluent + 4 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below.	4 vol 	1 vol	Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 10 mL of diluent + 40 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below.
Detection buffer			
The Detection buffer is ready-to-use.			The Detection buffer is ready-to-use.

## TO PREPARE ANTIBODY WORKING SOLUTIONS:

Each well requires 2 µL of Progranulin-Eu Cryptate Antibody and 2 µL of Progranulin-d2 Antibody.

Prepare the two antibody solutions in separate vials.

500 TESTS KIT - 62PROPEG		10,000 TESTS KIT - 62PROPEH	
Progranulin Eu Cryptate antibody			
Dilute 50-fold the 50X stock solution (thawed reagent) of Progranulin Eu Cryptate antibody stock solution with the Detection buffer #3 : add 1 volume of Eu Cryptate antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 20 µL of Eu Cryptate antibody stock solution + 980 µL of Detection Buffer #3).			Dilute 50-fold the 50X stock solution (thawed reagent) of Progranulin Eu Cryptate antibody stock solution with the Detection buffer #3 : add 1 volume of Eu Cryptate antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 0.4 mL of Eu Cryptate antibody stock solution + 19.6 mL of Detection Buffer #3).
Progranulin d2 antibody			
Dilute 50-fold the 50X stock solution (thawed reagent) of Progranulin d2 antibody stock solution with the Detection buffer #3 : add 1 volume of d2-antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 20 µL of d2-antibody stock solution + 980 µL of Detection Buffer #3).			Dilute 50-fold the 50X stock solution (thawed reagent) of Progranulin d2 antibody stock solution with the Detection buffer #3 : add 1 volume of d2 antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 0.4 mL of d2 antibody stock solution + 19.6 mL of Detection Buffer #3).
Antibody mix			
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 (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).			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 (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).

## TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 16  $\mu\text{L}$  of standard.
- Dilute the standard stock solution serially with diluent #5 (1X)
- In order to check for a potential interference effect from your own assay buffer when using the assay for the first time, we highly recommend the parallel preparation of a standard curve in your own supplemented cell culture medium and in diluent #5 (1X).
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

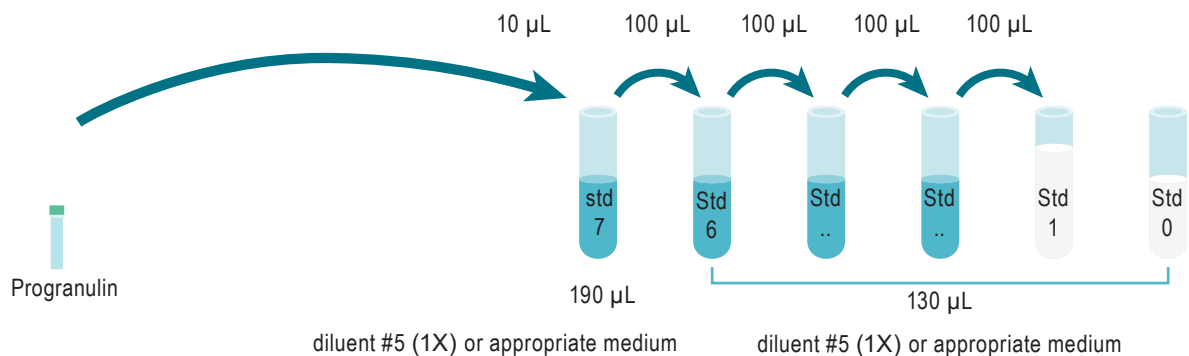
A recommended standard dilution procedure is listed and illustrated below:

Dilute the standard stock solution 20-fold with diluent #5 (1X) to prepare high standard (std 7): e.g. take 10  $\mu\text{L}$  of standard stock solution and add it to 190  $\mu\text{L}$  of diluent #5 (1X). Mix gently.

Use the high standard (std 7) to prepare the standard curve using 1/2.3 serial dilutions as follows:

- Dispense 130  $\mu\text{L}$  of diluent #5 (1X) in each vial from Std 6 to Std 0.
- Add 100  $\mu\text{L}$  of standard to 130  $\mu\text{L}$  of diluent #5 (1X), mix gently and repeat the 1/2.3 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 7 standards for the analyte. Std 0 (Negative control) is diluent #5 (1X) or appropriate culture medium alone.








STANDARD	SERIAL DILUTIONS	PROGRANULIN WORKING SOLUTIONS (PG/ML)
Standard Stock solution	Thawed stock solution	1,000,000
Standard 7	10 $\mu\text{L}$ stock solution + 190 $\mu\text{L}$ Diluent #5 (1X)	50,000
Standard 6	100 $\mu\text{L}$ standard 7 + 130 $\mu\text{L}$ Diluent #5 (1X)	21,739
Standard 5	100 $\mu\text{L}$ standard 6 + 130 $\mu\text{L}$ Diluent #5 (1X)	9,452
Standard 4	100 $\mu\text{L}$ standard 5 + 130 $\mu\text{L}$ Diluent #5 (1X)	4,109
Standard 3	100 $\mu\text{L}$ standard 4 + 130 $\mu\text{L}$ Diluent #5 (1X)	1,787
Standard 2	100 $\mu\text{L}$ standard 3 + 130 $\mu\text{L}$ Diluent #5 (1X)	777
Standard 1	100 $\mu\text{L}$ standard 2 + 130 $\mu\text{L}$ Diluent #5 (1X)	338
Standard 0	100 $\mu\text{L}$ Diluent #5 (1X)	0

## TO PREPARE SAMPLES:

- Each well requires 16  $\mu$ L of sample.
- Just after their collection, put the samples at 4°C and test them immediately. For later use, samples should be dispensed into disposable plastic vials and stored at -60°C or below. Avoid multiple freeze/thaw cycles.
- Samples with a concentration above the highest standard (std 7) must be diluted diluent #5 (1X) or in your appropriate sample medium.
- In order to measure human Progranulin in cell lysates, cells must be lysed with Lysis Buffer #4 (1X) for 30 min at RT under gentle shaking. Please note that the 4X stock solution of Lysis Buffer #4 must be ordered separately (Ref# 64KL4FDF, 130 mL) and 4-fold diluted with distilled water before use.

## ASSAY PROTOCOL

		Standard (Std 0 - std 7)	Samples
Step 1		Dispense 16 $\mu$ L of each Progranulin standard (Std 0 - std 7) into each standard well	Dispense 16 $\mu$ L of each sample into each sample well
Step 2		Add 2 $\mu$ L of Progranulin d2 antibody working solution to all wells	
Step 3		Add 2 $\mu$ L of Progranulin Eu Cryptate antibody working solution to all wells	
Step 4		Seal the plate and incubate overnight	
Step 5		Remove the plate sealer and read on an HTRF® compatible reader	



## DATA REDUCTION

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 ratio of the acceptor and donor emission signals for each individual well. The Standard 0 (Negative control) plays the role of an internal assay control.

$$\text{delta Ratio} = \text{Ratio Standard or sample} - \text{Ratio Standard 0}$$

For more information about data reduction, please visit <http://www.cisbio.com/htrf-ratio-and-data-reduction>

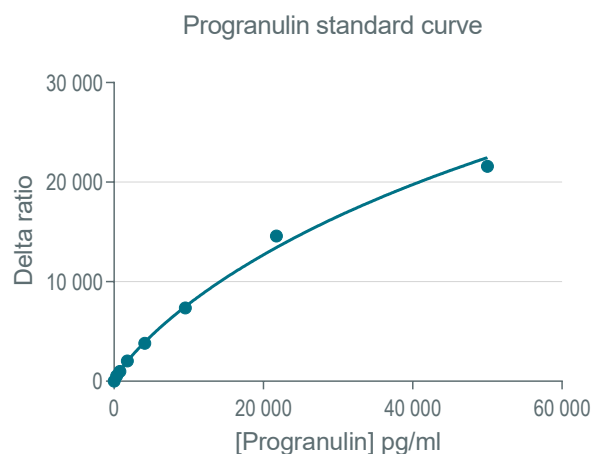
## RESULTS

This data must not be substituted for the data obtained in the laboratory and should be considered only as an example.

Results may vary from one HTRF® compatible reader to another.

Standard curve fitting with the 4 Parameter Logistic (4PL) model (with 1/Y<sup>2</sup> weighting):

	Ratio (1)	CV (2)	Delta Ratio
Standard 0 - Negative control	687	2%	0
Standard 1 - 338 pg/mL	1222	4%	535
Standard 2 - 777 pg/mL	1673	1%	986
Standard 3 - 1,787 pg/mL	2718	2%	2031
Standard 4 - 4,109 pg/mL	4489	1%	3802
Standard 5 - 9,452 pg/mL	8038	2%	7351
Standard 6 - 21,739 pg/mL	15262	1%	14575
Standard 7 - 50,000 pg/mL	22282	1%	21595



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.

Copyright 2020 Cisbio Bioassays. All rights reserved. HTRF and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays.

### FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880

Japan +81-(0)43-306-8712 Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio