

# Turn on the light with HTRF

## Technology

### Principle

Based on FRET technique :  
Light emission is caused by the proximity of 2 fluorophores.

NO interaction  
↓  
NO light

Interaction  
↓  
LIGHT

Energy transfer

⚡ : excitation

### Advantages

GET RID OF THE NOISE.  
Thanks to HTRF® stability, background fluorescence is reduced.



### Process

Simple Procedure:

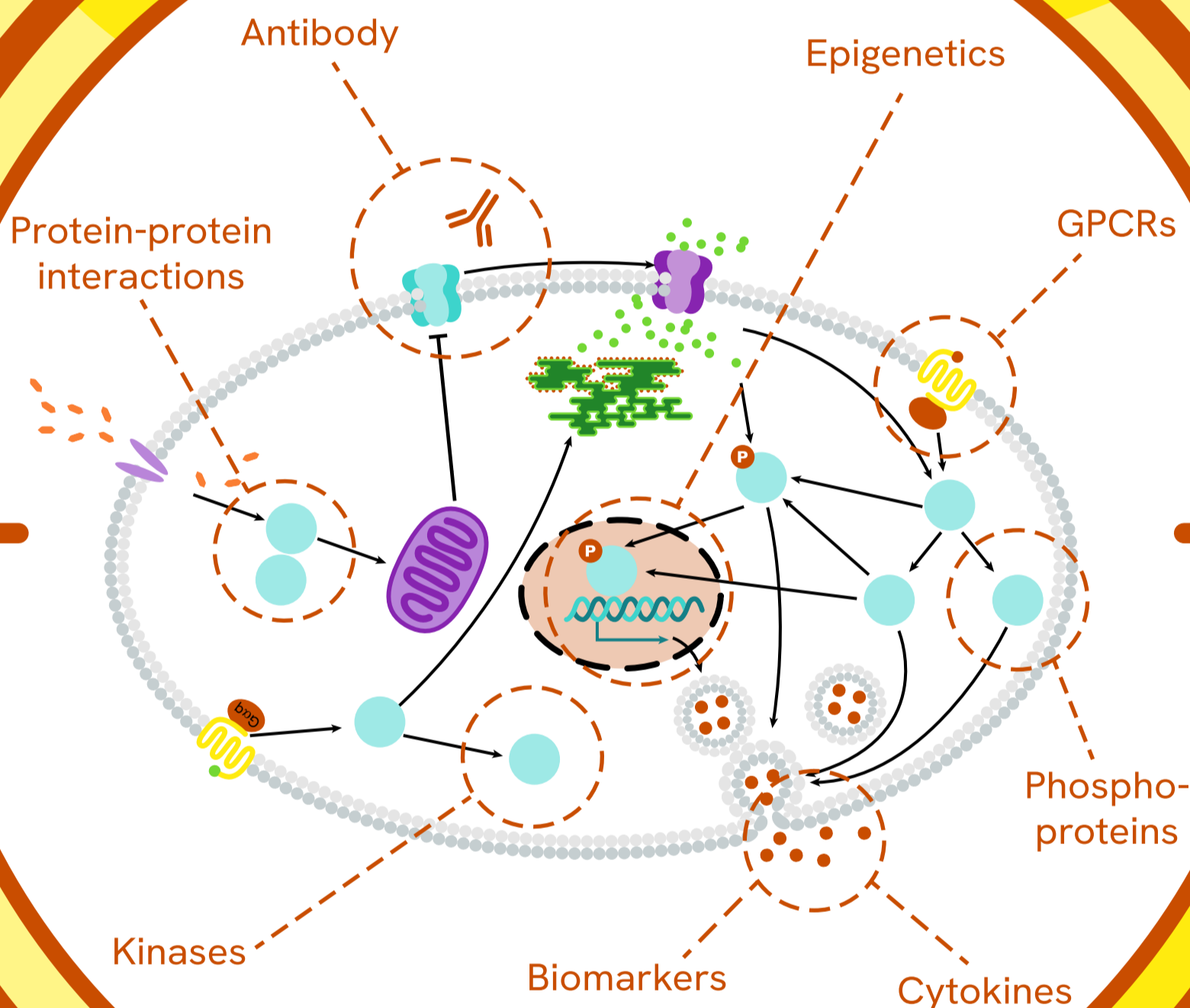
Add HTRF reagent  
↓  
Incubate  
↓  
Read

### About

Developed with  
Jean-Marie Lehn,  
Nobel prize  
in Chemistry  
(1987).

Over 25 years of experience  
Over 1,000 product references  
Over 9,000 scientific publications mentioning HTRF products

## Targets



## Applications

With HTRF, Revvity provides advanced products to boost your research on many therapeutic areas.



Cardiology



Diabetes



Inflammation



Immuno-oncology



NASH / fibrosis



Oncology



Neurosciences



Rare diseases



Custom services



### References

HTS-compatible FRET-based conformational sensors clarify membrane receptor activation.  
Scholler P, Moreno-Delgado D, Lecat-Guillet N, Doumazane E, Monnier C, Charrier-Savourin F, Fabre L, Chouvet C, Soldevila S, Lamarque L, Donsimoni G, Roux T, Zwier JM, Trinquet E, Rondard P, Pin JP. Nat Chem Biol. 2017 Apr;13(4):372-380

Time-resolved FRET strategy to screen GPCR ligand library.  
Oueslati N, Hounsou C, Belhocine A, Rodriguez T, Dupuis E, Zwier JM, Trinquet E, Pin JP, Durroux T. Methods Mol Biol. 2015;1272:23-36

Energy Transfer Luminescence of Europium(III) and Terbium(III) Cryptates of Macrobicyclic Polypyridine Ligands.  
Alpha B, Lehn JM, Mathis G. Angew Chem Int Ed Engl. 1987;26:266-7