

# ***Bioluminescent Cellular Reporter Technologies Featuring NanoLuc Luciferase***

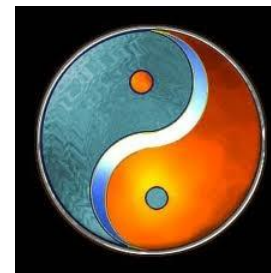
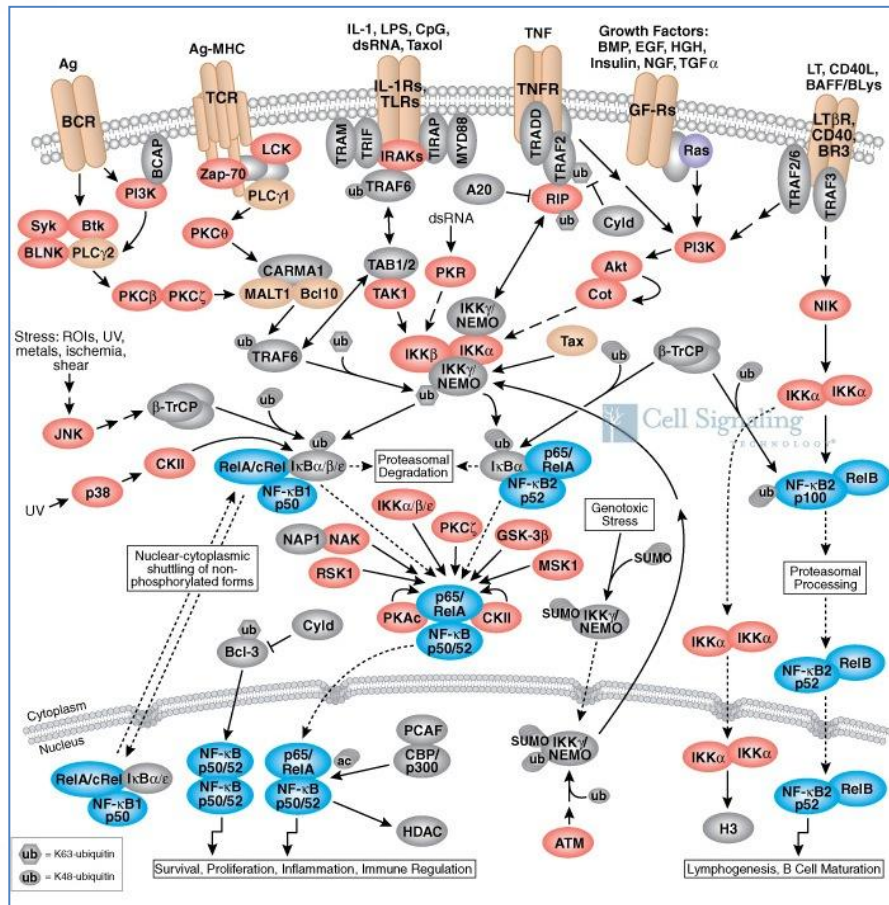
**Frank Fan, Ph.D.  
Director of Research**

**March, 2014**

# *Outline*

- Bioluminescence
- NanoLuc luciferase
- Genetic reporter
  - Dual luciferase assays
  - Animal imaging
- Protein reporter
  - Location
  - Stability
  - Interactions (NanoBRET, NanoBiT)

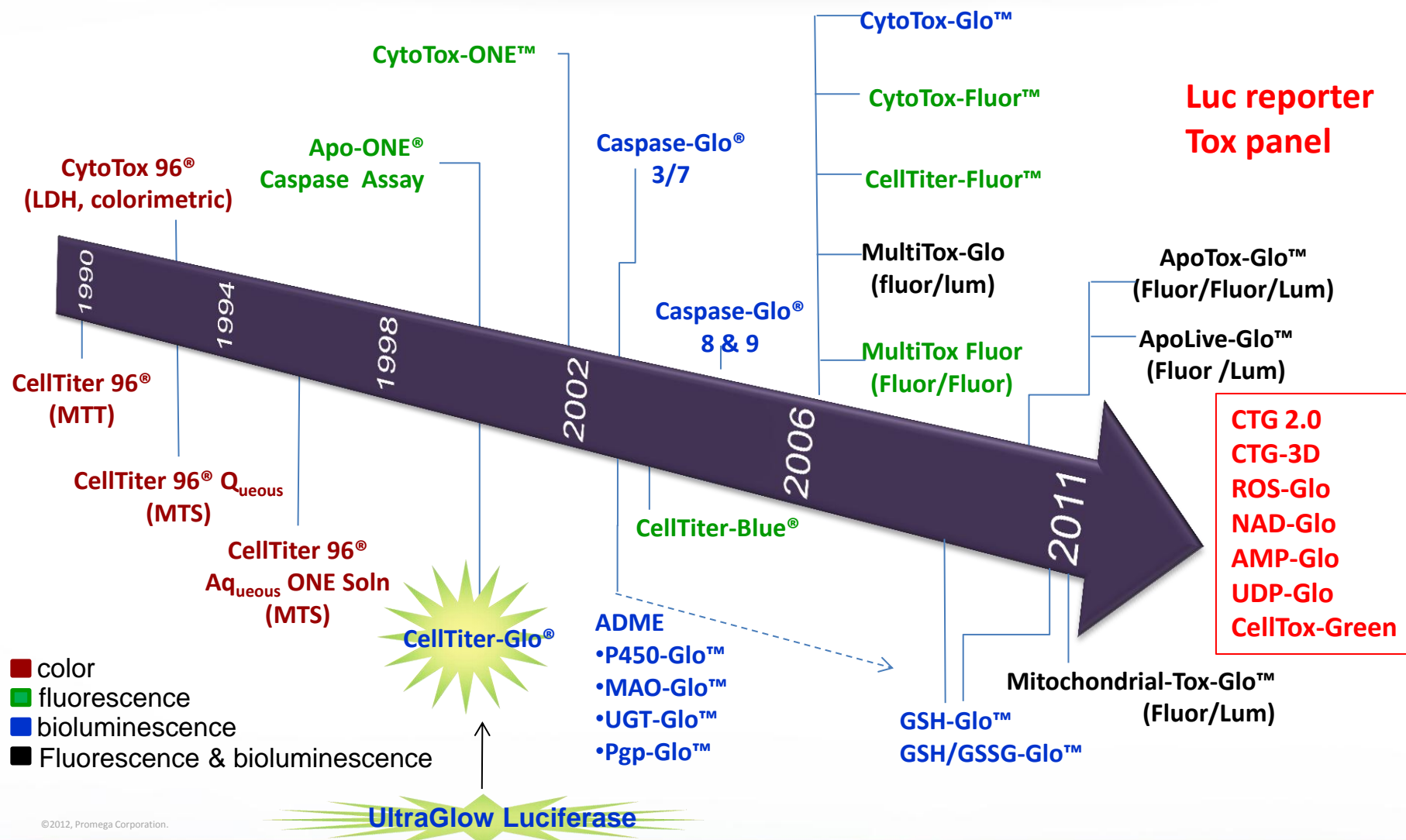
# Complicated Biology ---- Simple Assay



## Add-Mix-Read



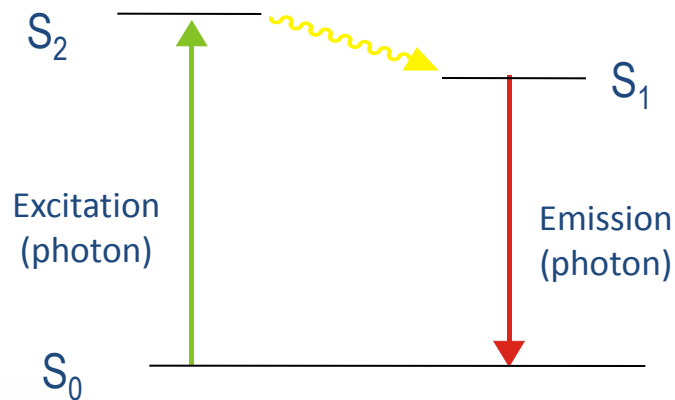
# Evolution of assay design



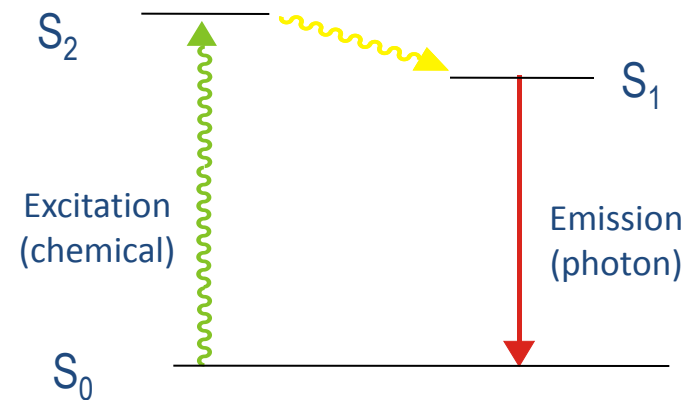
# Fluorescence vs Luminescence

- Fluorescence and luminescence differ primarily by the method for creating the excited-state of the photon emitter.
- The difference in how the excited state is created dictates practical differences in how these technologies perform.

Fluorescence

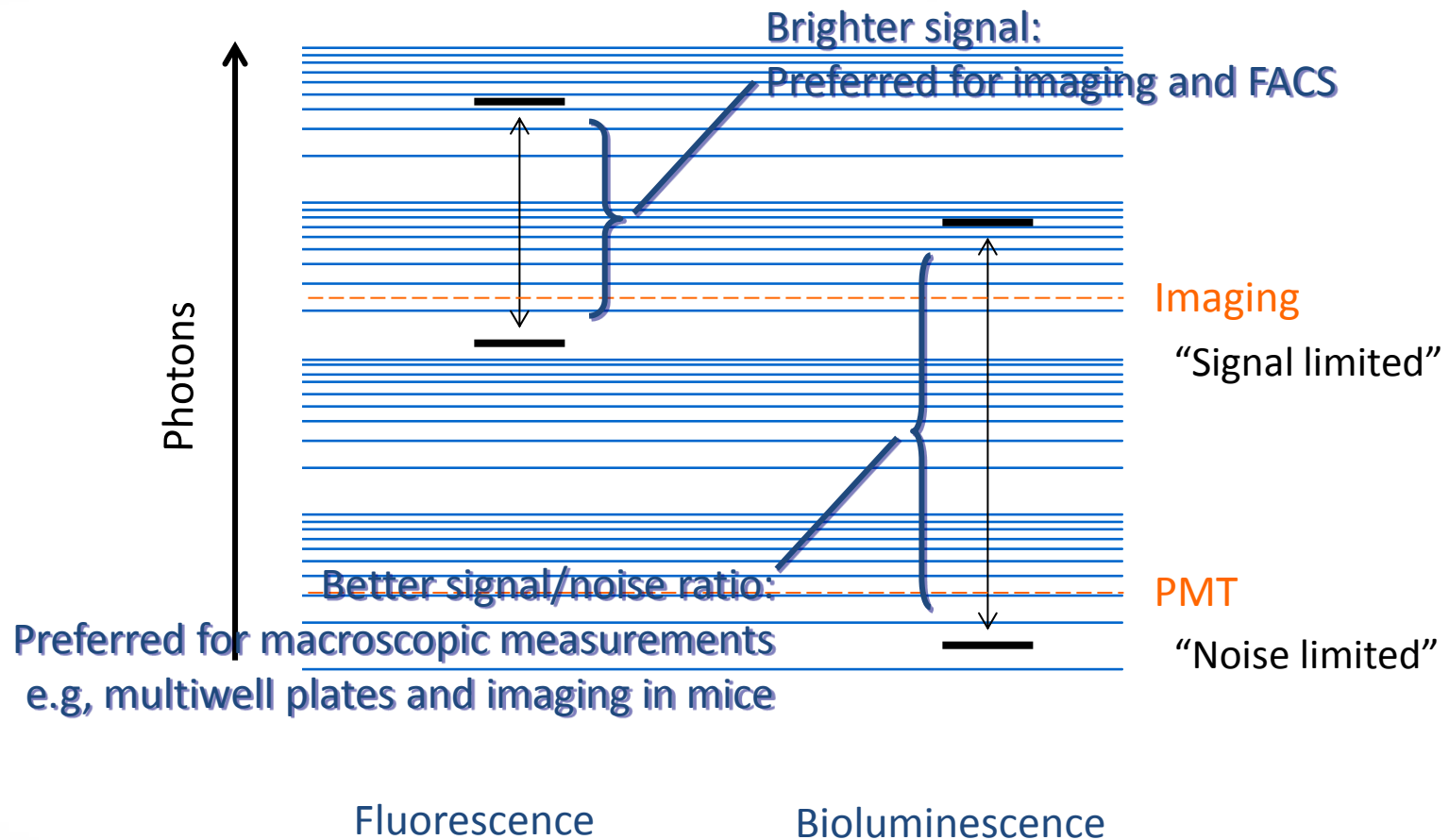


Luminescence

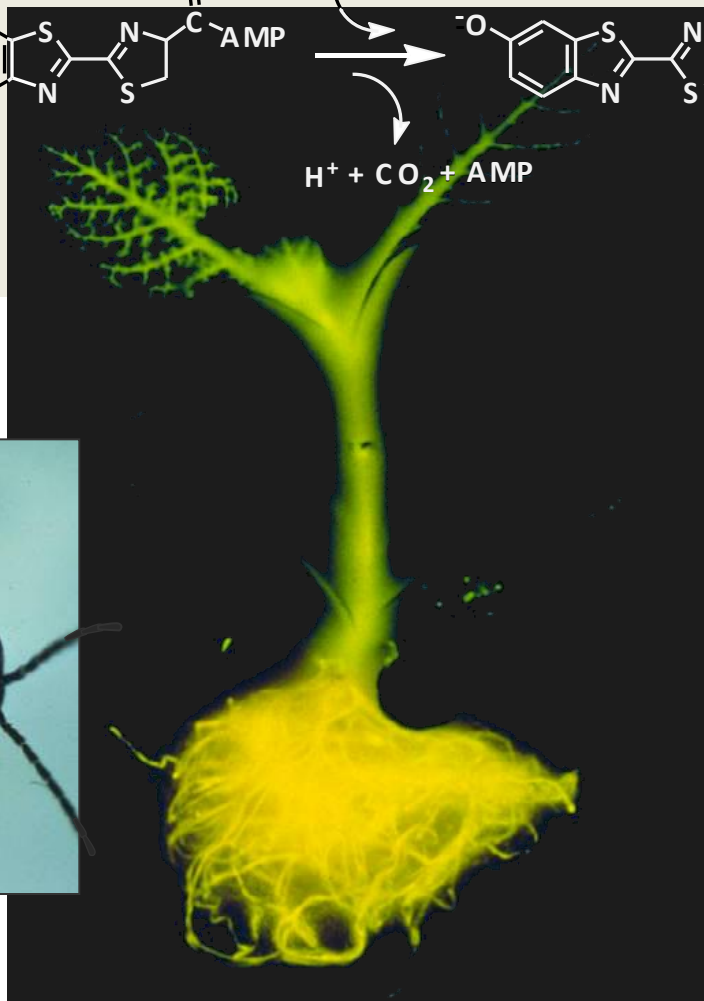
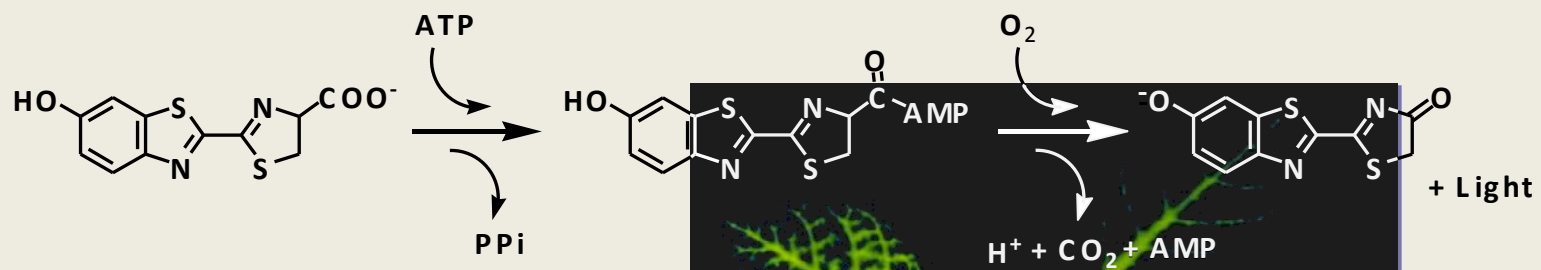




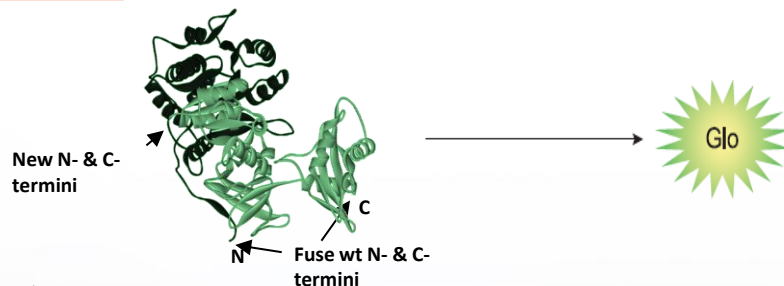
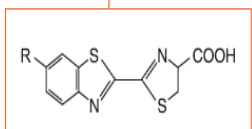
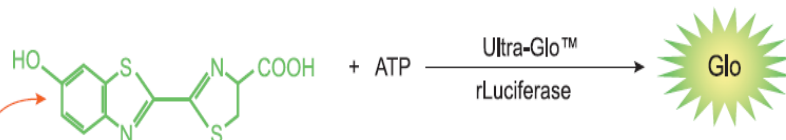
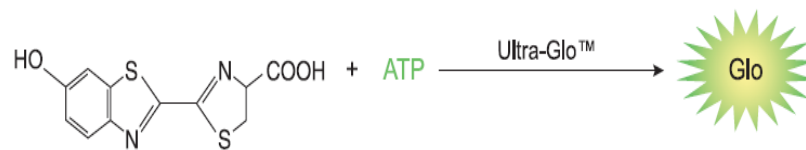
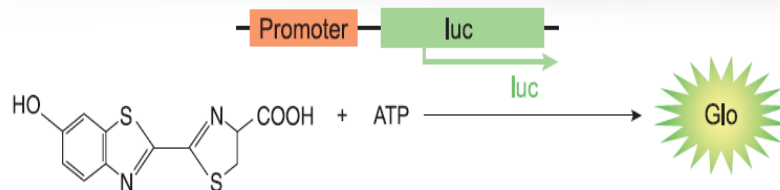
# Luminescent assay has superior sensitivity and dynamic range



Great for HTS, why?



# Bioluminescent Assays Developed From Firefly Luciferase Chemistry



- Reporter gene assays  
GPCR & Nuclear Receptor

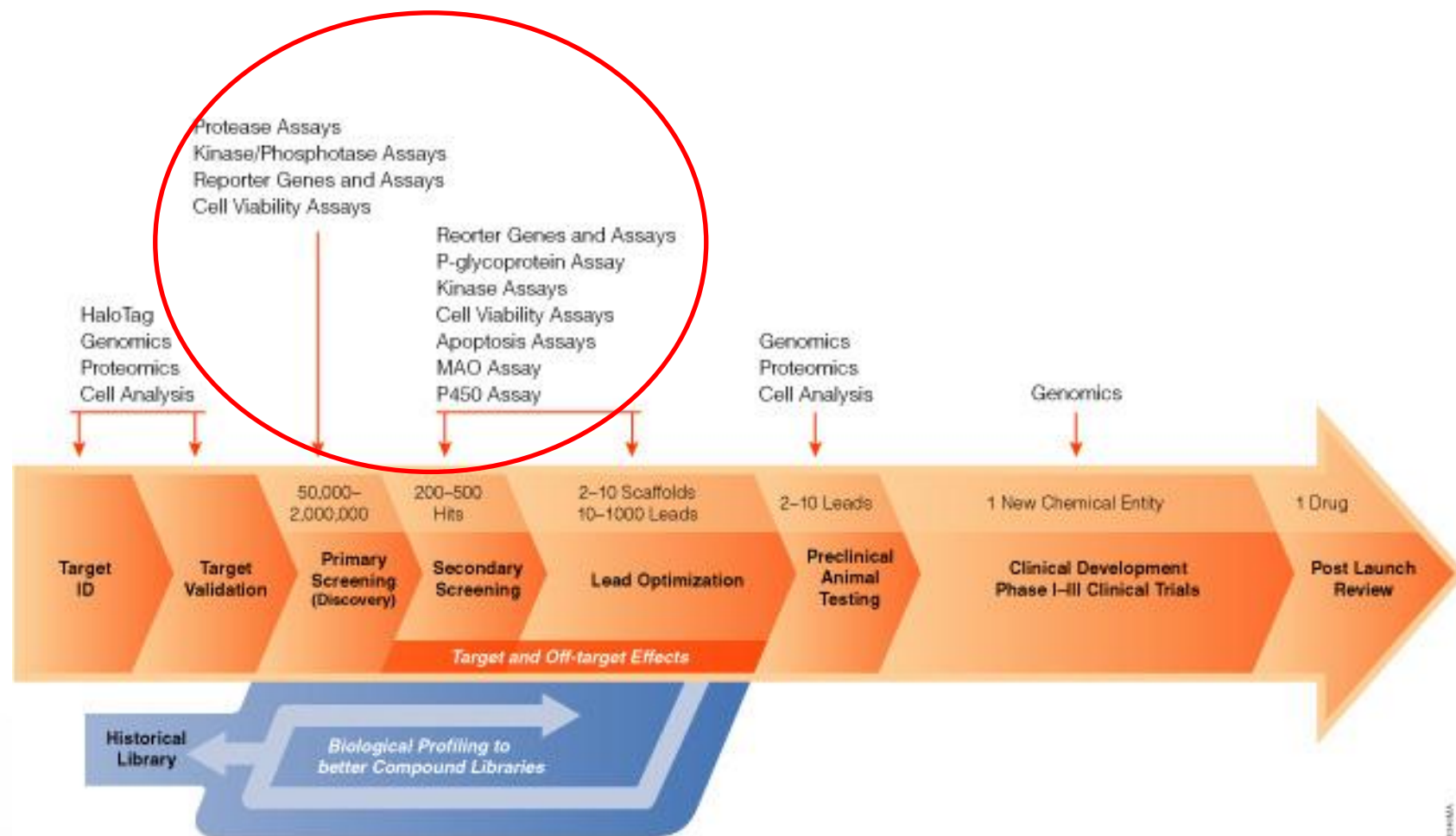
- Cell Viability
- Kinase assays
- cAMP & PDE assays
- P-glycoprotein assay

- Caspases/proteases
- CYP450 assays
- MAO assays
- GSH assays
- HDAC assays

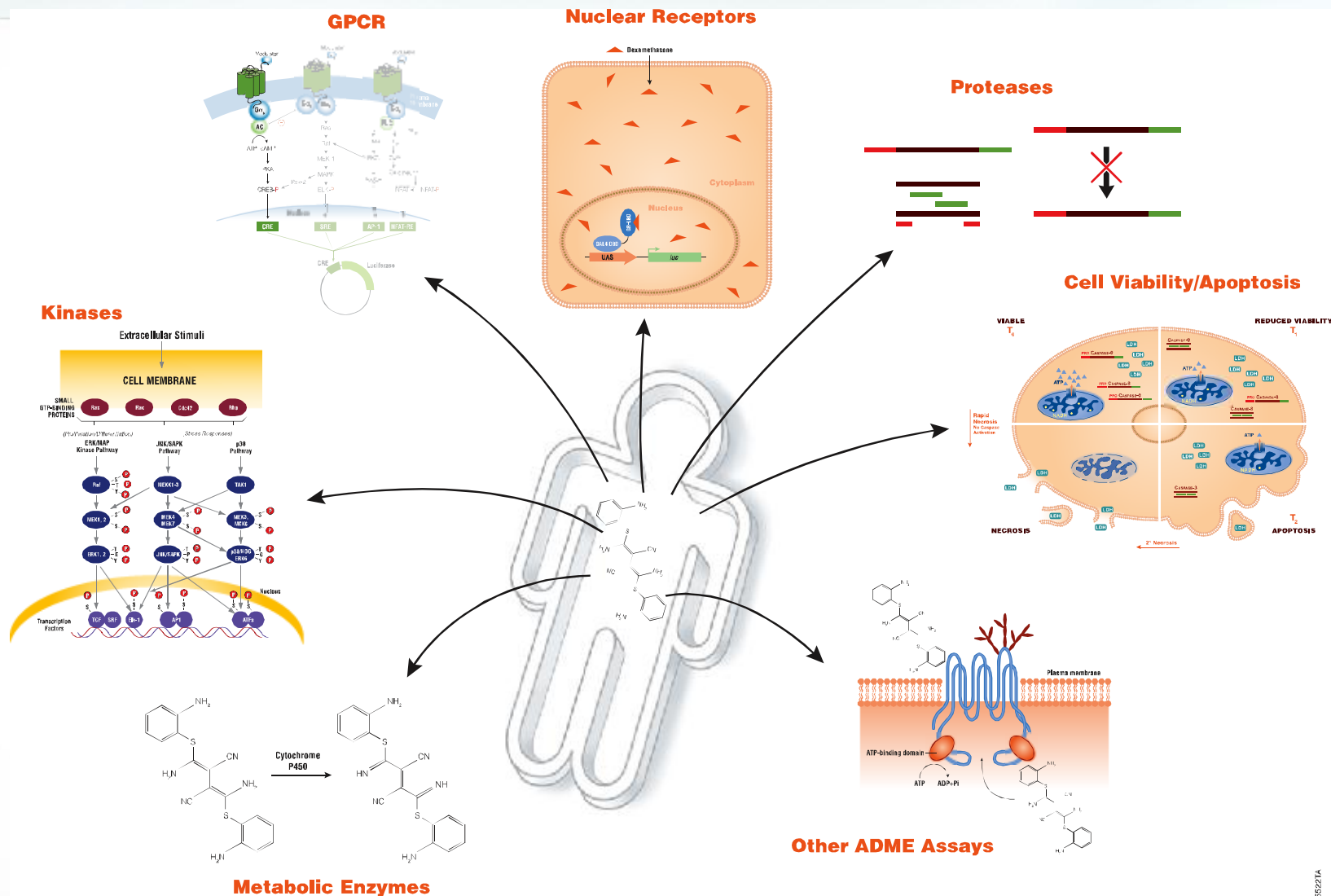
- GloSensor cAMP
- GloSensor caspase



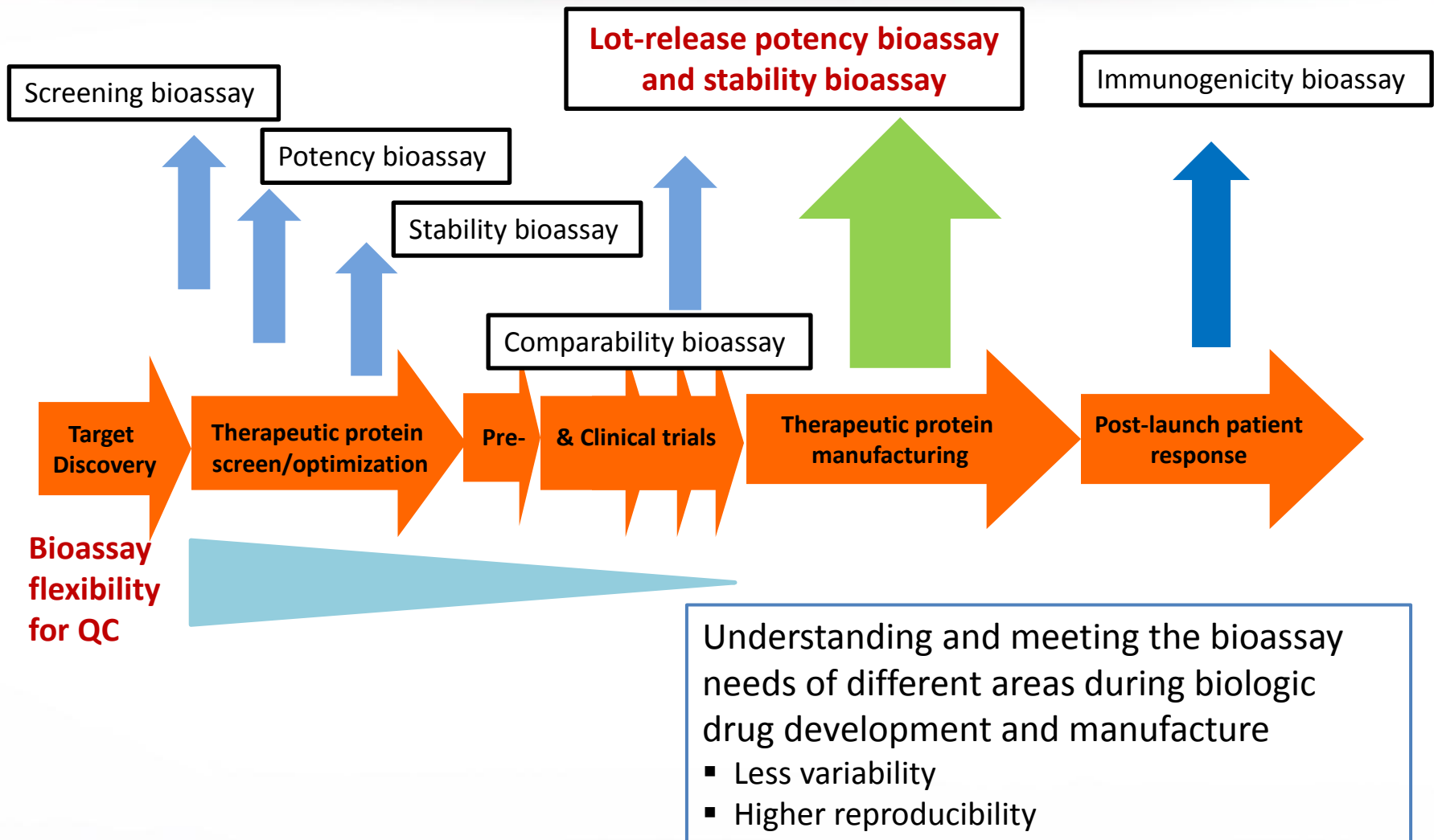
# ***Bioluminescent assays for chemical drug discovery***



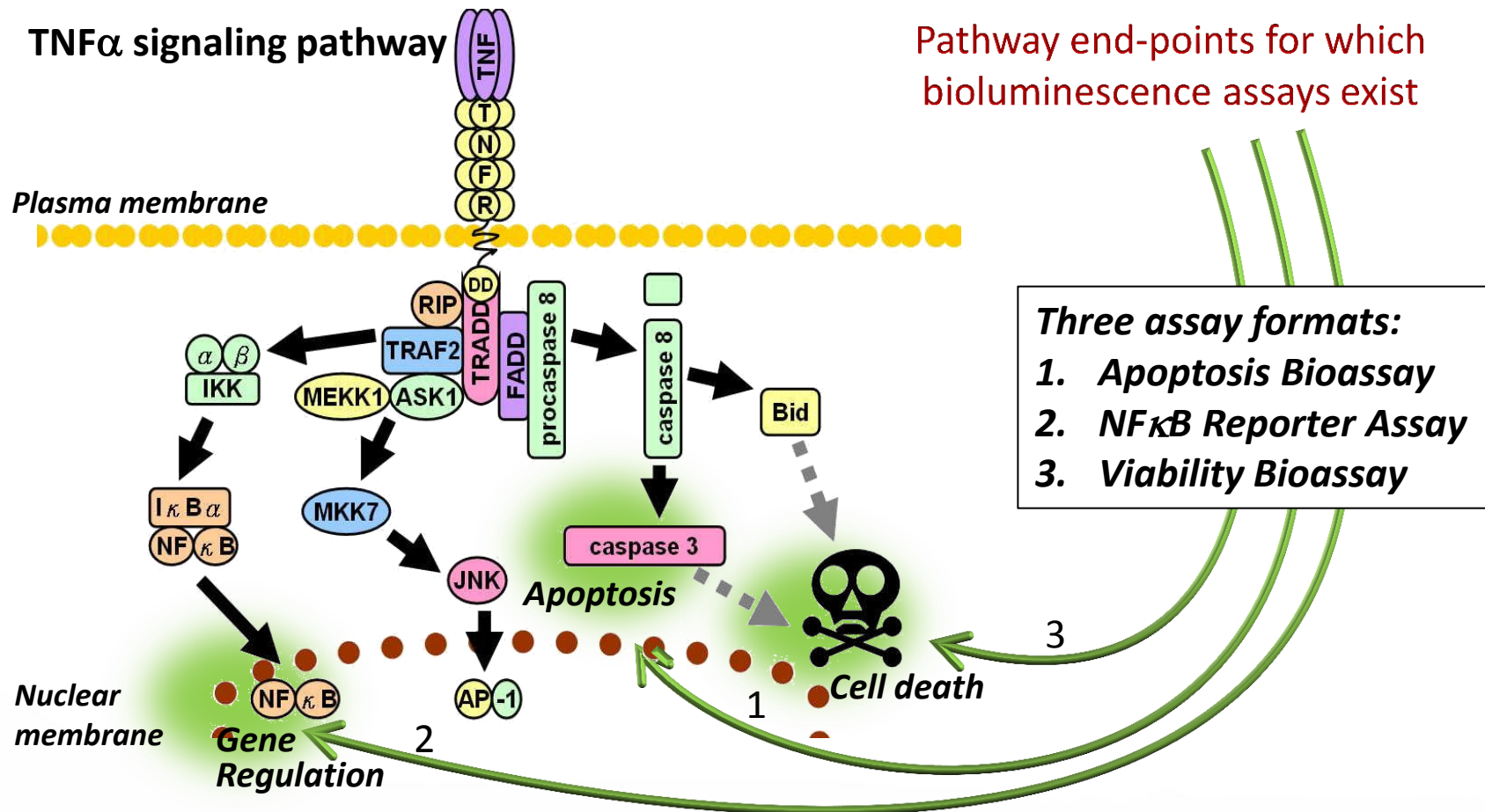
# Bioluminescent assays for chemical drug discovery



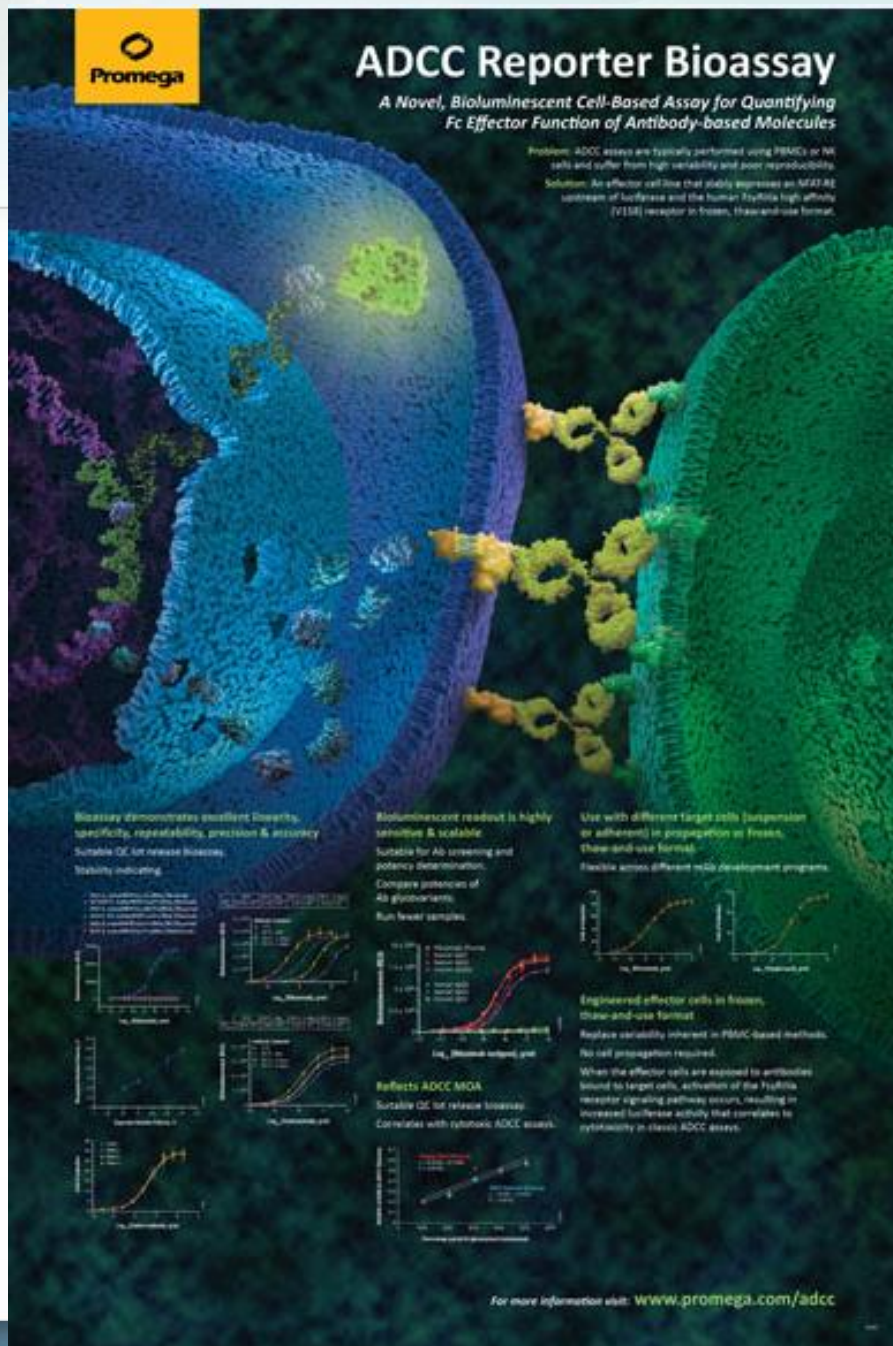
# ***Bioluminescent bioassays for antibody drug development***



# Bioluminescent Cell-based Bioassays







**ADCC Reporter Bioassay**  
A Novel, Bioluminescent Cell-Based Assay for Quantifying Fc Effector Function of Antibody-based Molecules

**Problem:** ADCC assays are typically performed using PBMCs or NK cells and suffer from high variability and poor reproducibility.

**Solution:** An effector cell line that stably expresses an NFAT2.B2 construct of luciferase and the human FcγRIIIa high affinity (V158) receptor in frozen, thaw-and-use format.

**Bioassay demonstrates excellent sensitivity, specificity, repeatability, precision & accuracy.**  
Suitable QC for release bioassays.  
Stability indicating.

**Bioluminescent readout is highly sensitive & scalable.**  
Suitable for Ab screening and potency determination.  
Compare potencies of Ab glycoforms.  
Run fewer samples.

**Use with different target cells (suspension or adherent) in propagation or frozen, thaw-and-use format.**  
Flexible across different cell development programs.

**Engineered effector cells in frozen, thaw-and-use format.**  
Replace variability inherent in PBMC-based methods.  
No cell propagation required.  
When the effector cells are exposed to antibodies bound to target cells, activation of the FcγRIIIa receptor signaling pathway occurs, resulting in increased luciferase activity that correlates to cytotoxicity in classic ADCC assays.

**Reflects ADCC MOA.**  
Suitable QC for release bioassays.  
Correlates with cytotoxic ADCC assays.

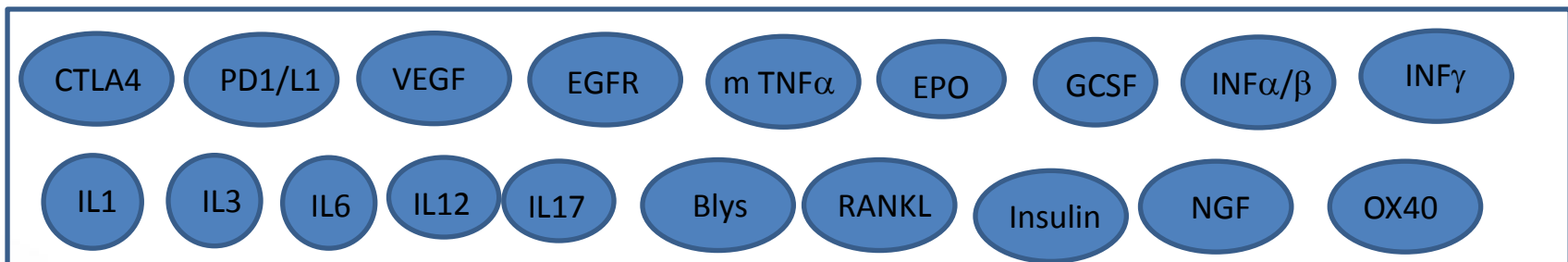
For more information visit: [www.promega.com/adcc](http://www.promega.com/adcc)

## Top10 Innovations 2013 The Scientist Magazine

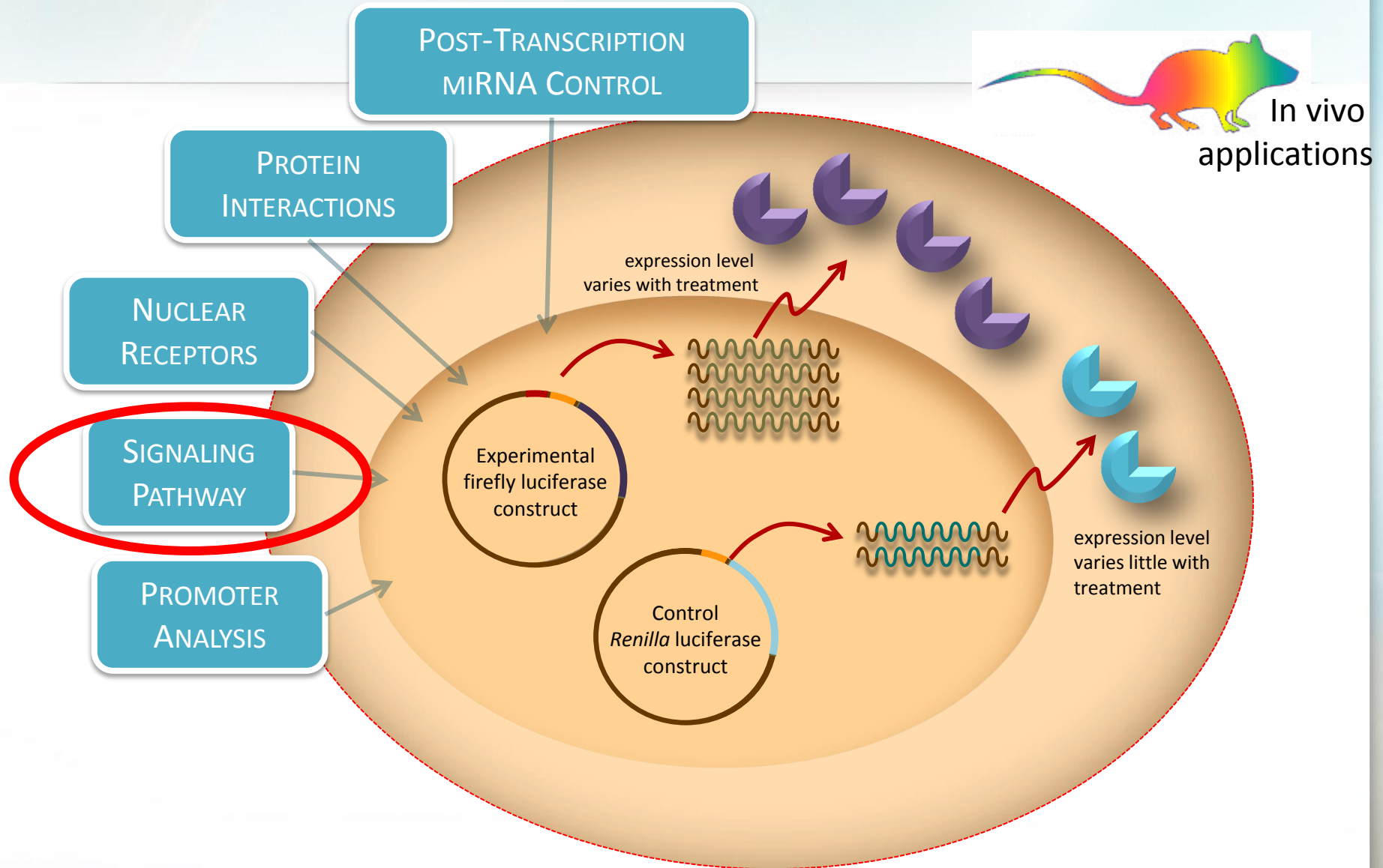


## ***Bioluminescent bioassays***

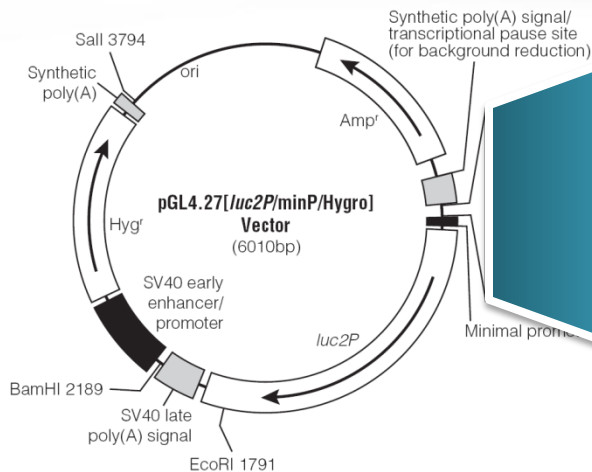
- ADCC Reporter Assay F version
- Mouse ADCC Assay
- ADCP Reporter Assay
- T cell Activation Assay



# Application Overview



# Promega offers complete solutions



Pre-designed,  
ready-to-use  
response  
element pGL4  
Vectors

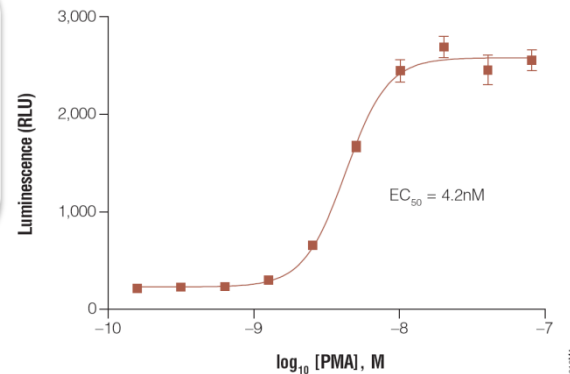
- Genes
- Vectors
- Cell lines
- Complete assays
- Catalog
- CAS
- CAM

HEK 293 Hyg<sup>R</sup> cell line  
made with the pGL4 vectors



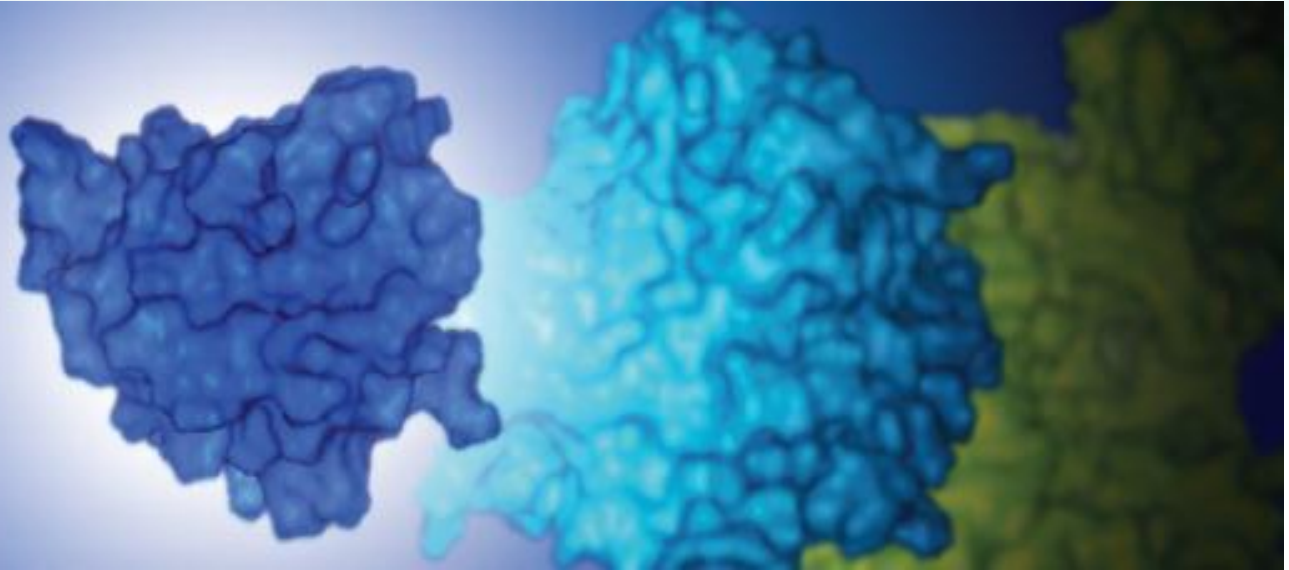
GloResponse<sup>™</sup>  
HEK 293 cell lines:

- cAMP RE
- NFAT RE
- NF- $\kappa$ B RE



Dose response curve  
GloResponse NFAT-RE-luc2P  
HEK293 Cell Line

# NanoLuc™ Luciferase Technology



## NanoLuc Luciferase as a genetic reporter

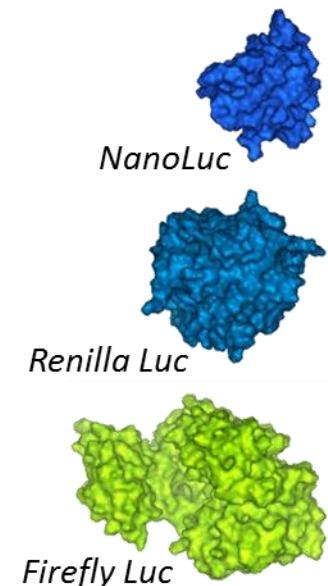
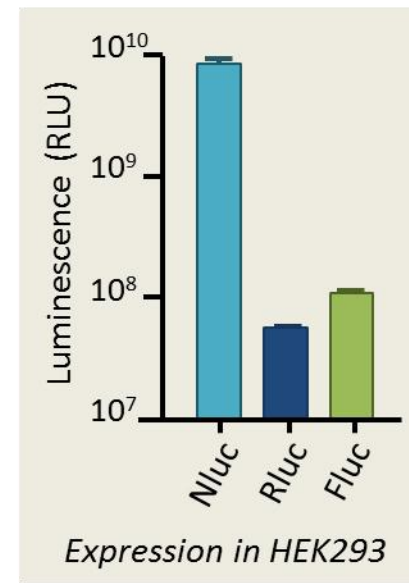
- Brighter, more responsive
- Smaller
- More versatile

## NanoLuc Luciferase as a protein function reporter

- Protein Translocation
- Protein Stability
- Protein:Protein Interactions

# ***NanoLuc: a new generation of luciferase technology***

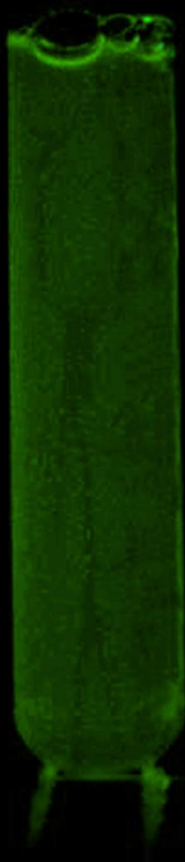
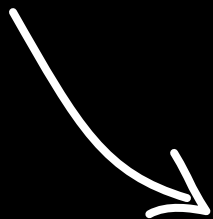
- **Extremely bright**
  - 100-fold greater light intensity
  - Much brighter in mammalian cells
  - Glow kinetics
- **Very small (19kD)**
  - Much smaller than Fluc or Rluc
  - Smaller than GFP
  - Ideal as fusion tag
- **High physical stability**
  - Compatible with fusion partners
  - Robust to experimental conditions
- **Narrow emission spectrum**
  - Less overlap with energy acceptor
  - Greater detection sensitivity



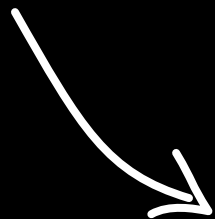
*Deep sea shrimp secretes extremely bright cloud of luminescence as a survival mechanism*



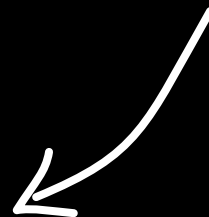
firefly



*firefly*



*NanoLuc*

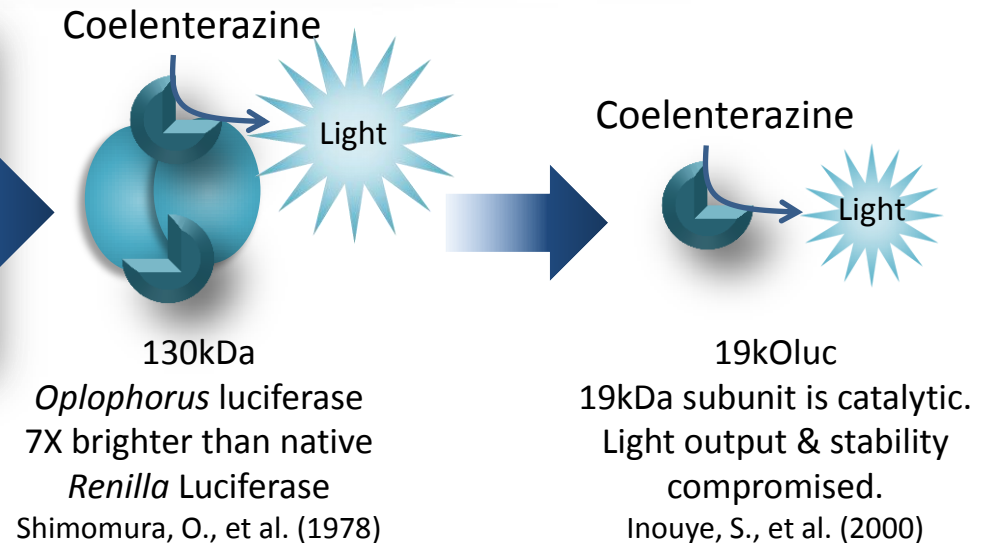


# Building a Better Luciferase

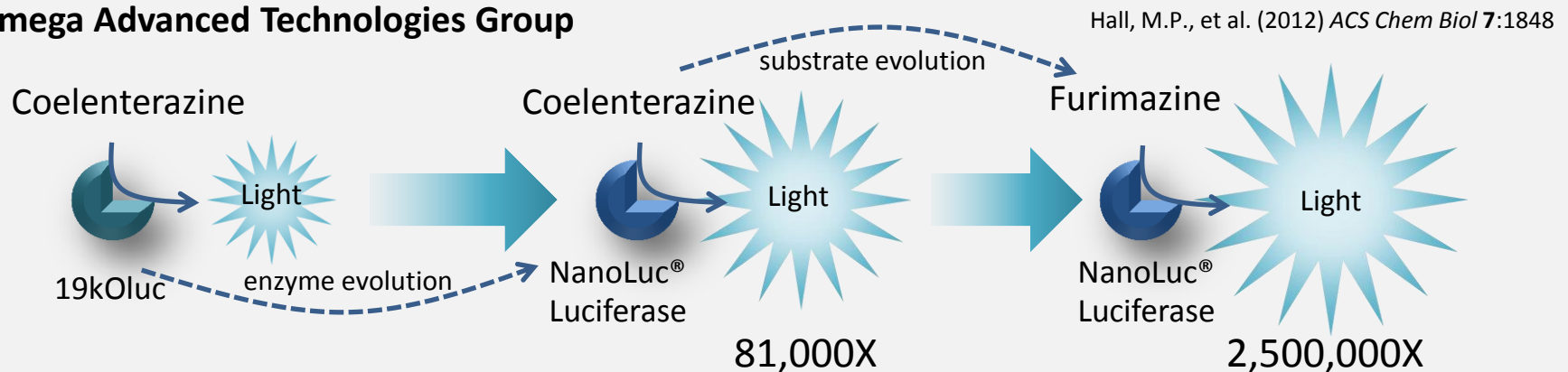
## NanoLuc<sup>®</sup> Evolution: From Ocean to Lab Bench



*Oplophorus gracilirostris* first cataloged  
in 1881

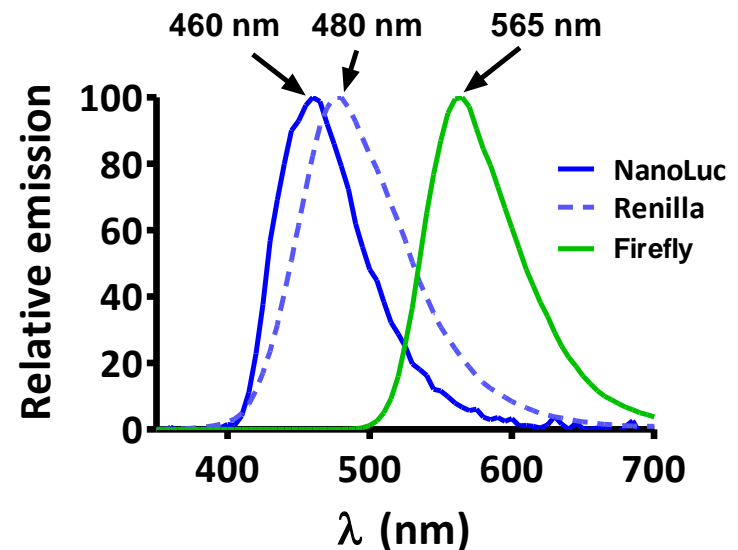


### Promega Advanced Technologies Group

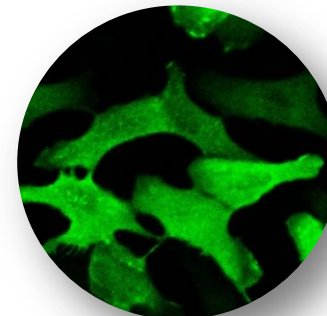


# Excellent Physical Properties Make NanoLuc® a Powerful Research Tool

- ✓ **Thermal stable**
  - Retains activity after 30 min @ 55°C
  - Melting temps: Nluc, 58°C; Fluc, 31°C
- ✓ **Active over broad pH range**
  - Fully active between pH 7-9
  - Retains significant activity at pH 5-7
  - Fluc: sharp decrease in activity below pH 8
- ✓ **Monomeric enzyme**
  - Facilitates use as transcriptional reporter or fusion partner
- ✓ **No mammalian post-translational modifications detected**
- ✓ **No disulfide bonds**
  - Supports high levels of activity in living cells
- ✓ **Uniform distribution in cells**
  - No apparent compartmental bias in the absence of targeting sequences

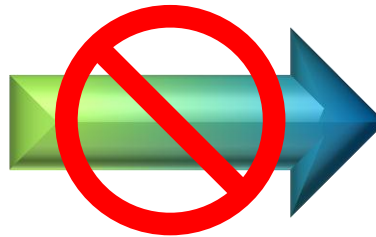
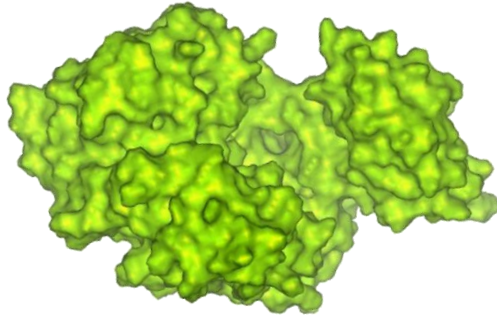


Unfused Nluc Immunofluorescence



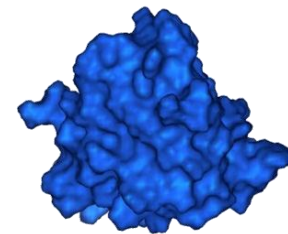
# Should I switch from Firefly to NanoLuc™ Luciferase?

*Firefly (Fluc)*



***Not necessarily***

*NanoLuc™ (Nluc)*



**Does it allow you to do your work?**  
**Do you plan to do work in vivo?**

**Firefly is a great reporter**

- ✓ Excellent signal:background
- ✓ Excellent dynamic range

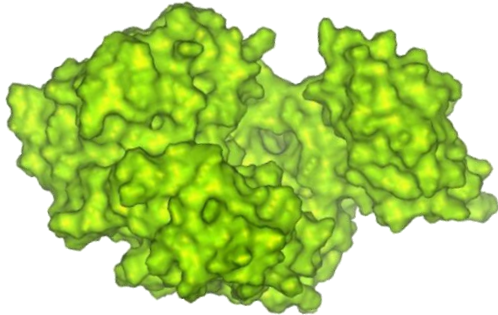
We just released new response element signaling pathway detection pGL4 vectors:

ARE	HSE	ISRE	STAT5	SRE
p53	HRE	SIE	NFAT	SRF
ATF6	XRE	SBE	CRE	
MRE	AP1	TCF-LEF	NF-κB	

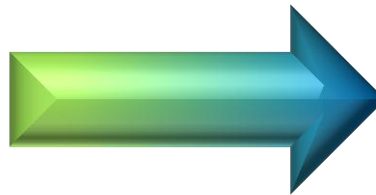


## *Should I switch from Firefly to NanoLuc™ Luciferase?*

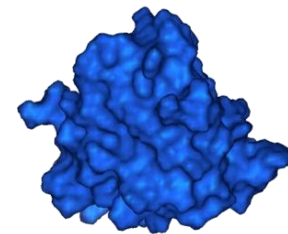
*Firefly (Fluc)*



*Yes, if ...*

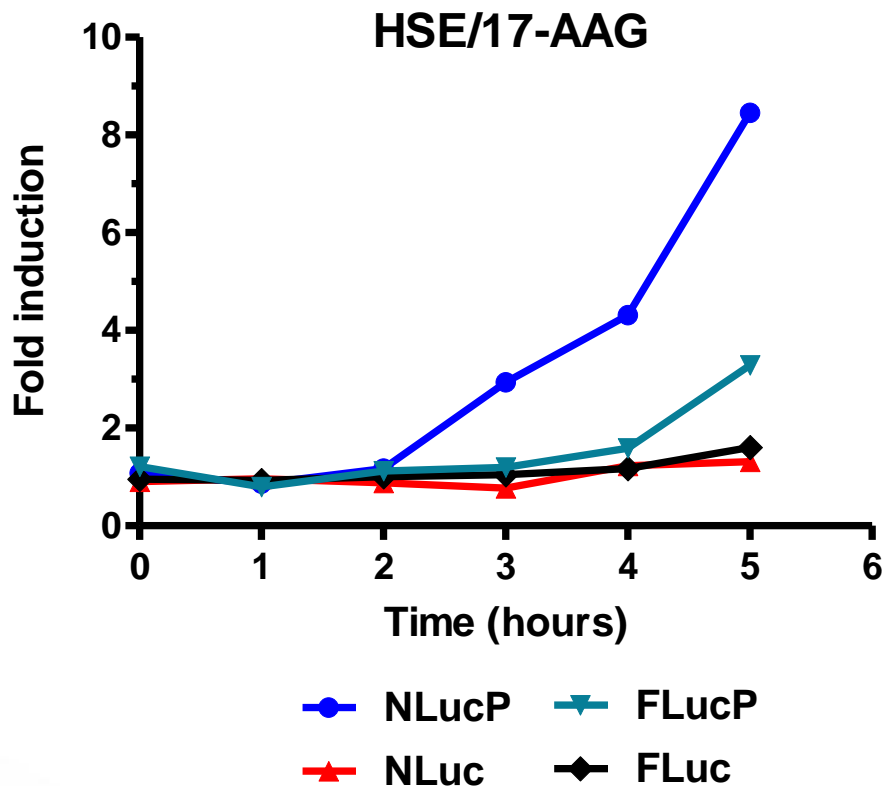


*NanoLuc™ (Nluc)*



- ✓ **Transfection efficiency limits you to easy-to-transfect cell lines**
- ✓ **Signals are too weak to move to 96-well plates**
- ✓ **FLuc is just too big**
- The increased brightness could allow a subtle signal become a reliable signal.
- The small size could allow gene replacement with minimal impact, especially in viral constructs

## Relative response of reporter genes



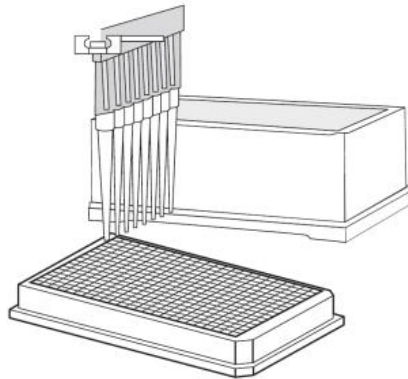
### Relative Response

**NLucP** > **FLucP** > **Fluc, NLuc**

Experimental details: transient transfection of HeLa cells w/ Hsf1 inducible constructs; addition of 500 nM 17-AAG at time zero.

# ***NanoDLR: Fluc then Nluc multiplex -in development***

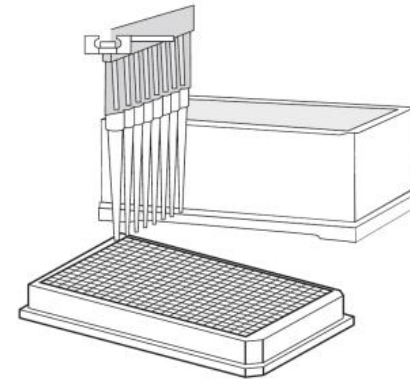
*Read #1: Fluc*



+ Fluc inhibitor  
+ Furimazine



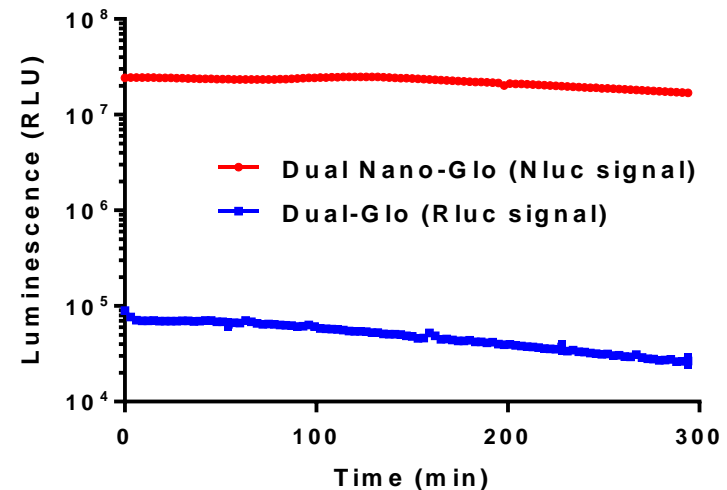
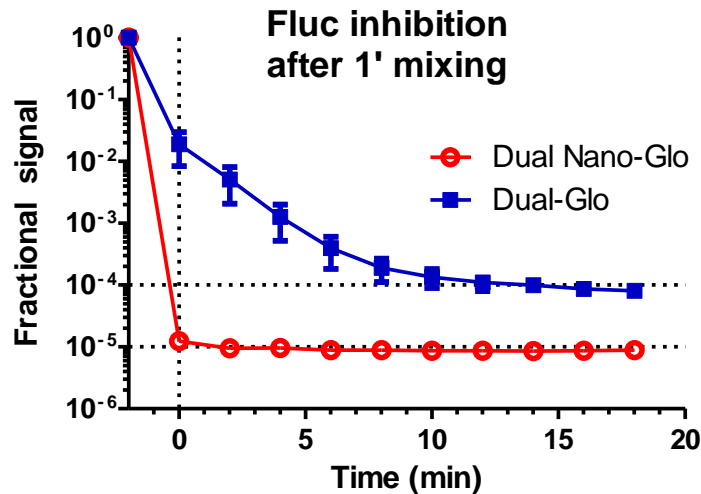
*Read #2: Nluc*



Possible formats (+/- PEST in all cases):

- Fluc dynamic/Nluc control
- Fluc control/Nluc dynamic
- Fluc dynamic/Nluc dynamic

# NanoDLR Stop & Glo reagent

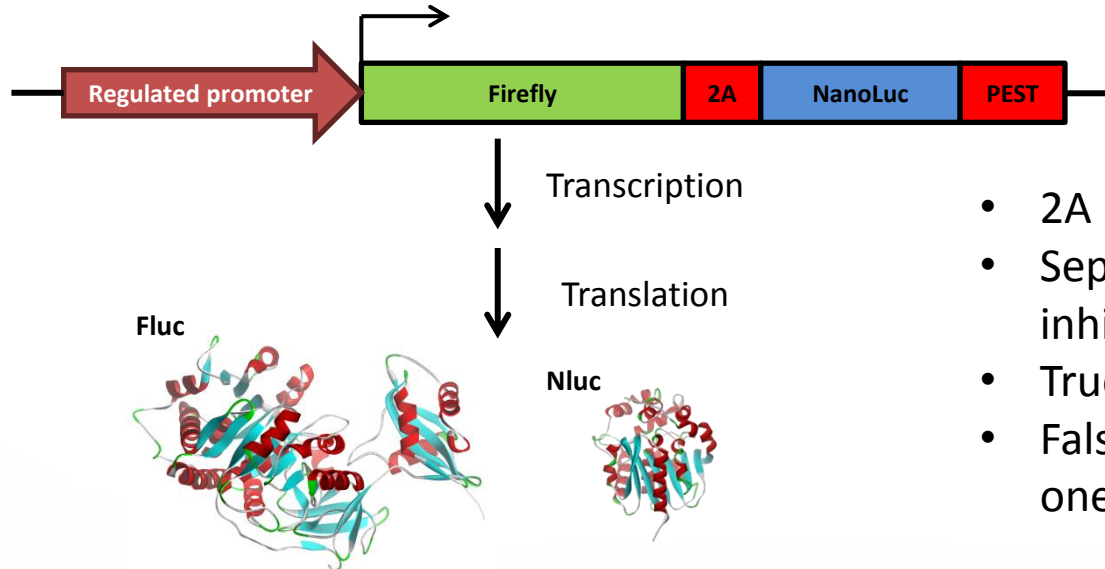


- Provides faster and more potent Fluc inhibition
  - Quickly take both reads for a single plate to facilitate HTS work-flow
  - Minimizes background for Nluc detection
- Large increase in sensitivity for detection of Nluc vs. Rluc
  - Lower residual Fluc signal
  - Nluc **much** brighter than Rluc

# NanoLuc and firefly Luc in HTS

- Compounds that inhibit/stabilize luciferase give too many false positives
  - Huge resource commitment to find the real hits
- Nanoluc less susceptible than other luciferases, but NlucP more susceptible than Nluc

## Co-incidence reporter developed at NIH NCGC\*

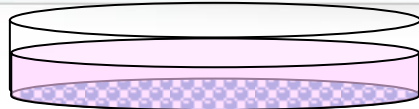


- 2A promotes 'ribosome skip'
- Separate Luc enzymes w/ own inhibitor profiles
- True hits show induction of both
- False positives show induction of one, not both

\* Cheng & Inglesse *Nature Methods* vol. 9, p. 937.



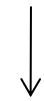
# Insert a luciferase biomarker in the test cells



tumor cells



stably transfect  
with luciferase  
construct



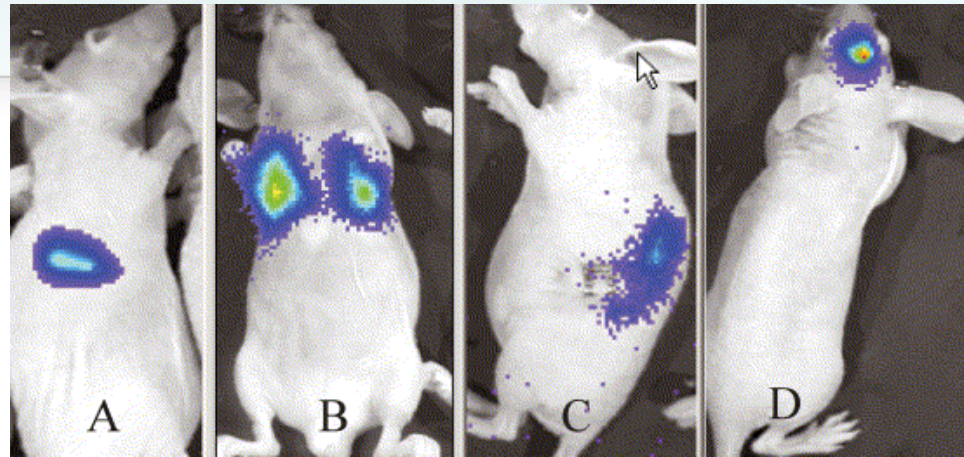
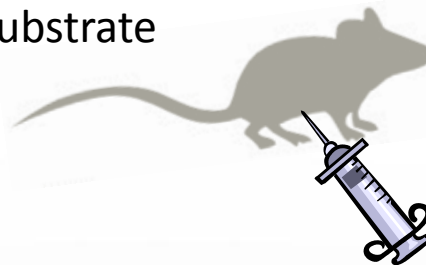
Inject cells



treat



Inject  
Luciferase  
substrate



Bioluminescent imaging with CCD device  
(overlaid on mouse photos)

In this case, luciferase is acting as a viability marker.  
This model can be used to evaluate treatments in vivo.

# **NanoLuc<sup>®</sup> for Viral Applications**

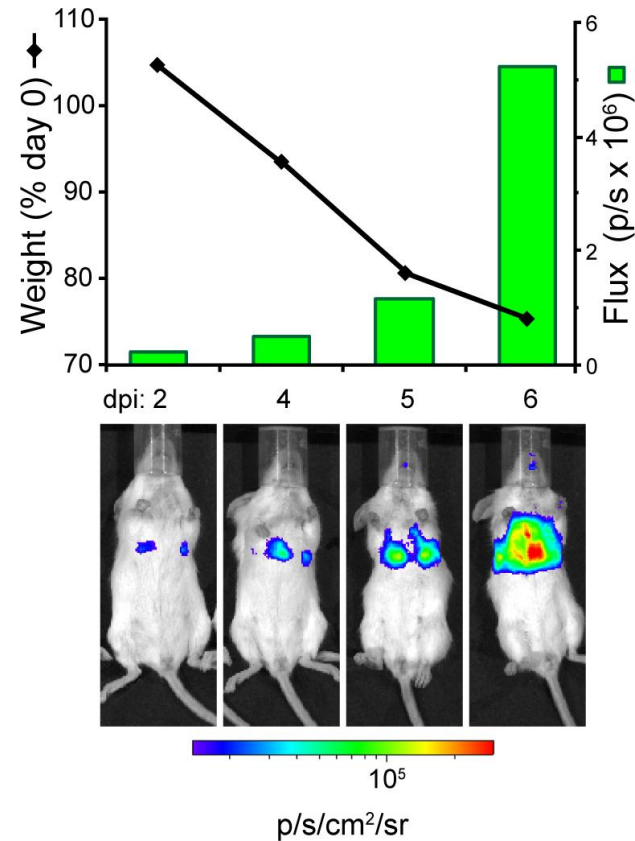
## *Small Size is Highly Advantageous*



- Data presented is from **Prof. Andrew Mehle's Lab** at the University of Wisconsin
- Historically, it's been difficult to stably integrate reporter genes into influenza
- Created influenza recombinant virus stably expressing NanoLuc<sup>®</sup>
  - No details on construction provided here
  - If you're interested in working with Andy or would like more info, please contact him directly:  
[amehle@wisc.edu](mailto:amehle@wisc.edu)

# Influenza - NanoLuc<sup>®</sup> Reporter Virus

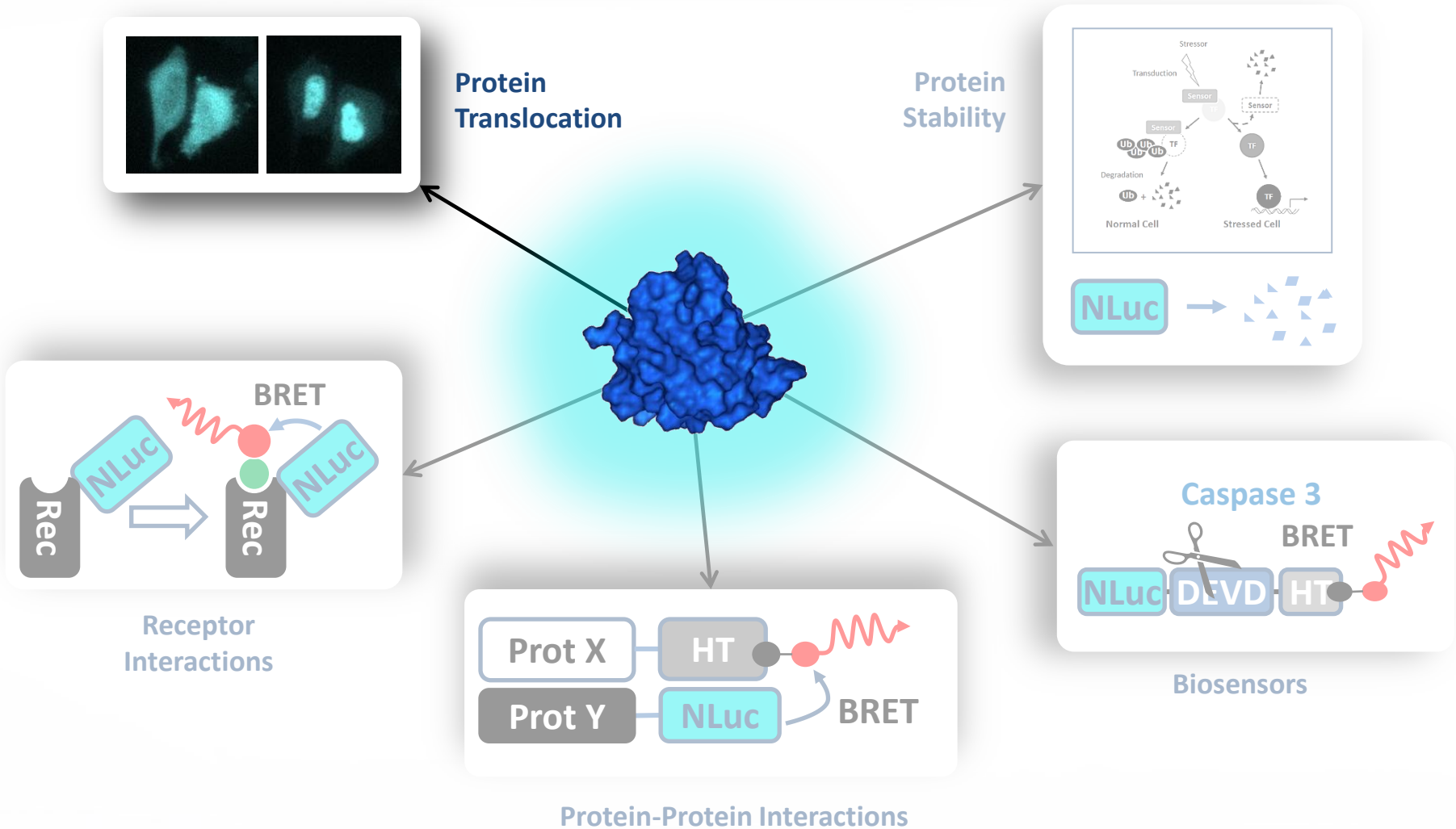
## Real-time In Vivo Imaging of Infection



# 20687

[http://en.wikipedia.org/wiki/ENCODE\\_\(ENCyclopedia\\_Of\\_DNA\\_Elements\)](http://en.wikipedia.org/wiki/ENCODE_(ENCyclopedia_Of_DNA_Elements))

# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments





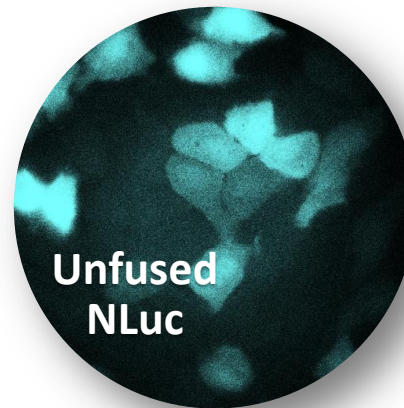
# *NanoLuc™ Luciferase excels in bioluminescent imaging applications*

Nluc brightness leads to short exposure times:

- Fluc/Rluc: 1-5min/exposure
- Nluc: 1-5sec/exposure

Why bother? Fluorescence works.

- Fluors are susceptible to photobleaching.
- Excitation can cause autofluorescence of other fluors
- Luciferases will generate light as long as substrate is available



*NanoLuc &  
LV200 featured*

*@ASCB 2012*

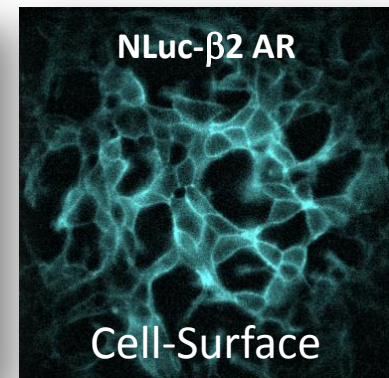
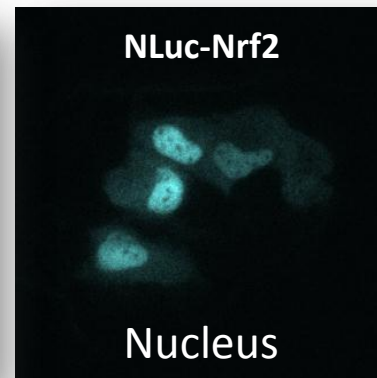
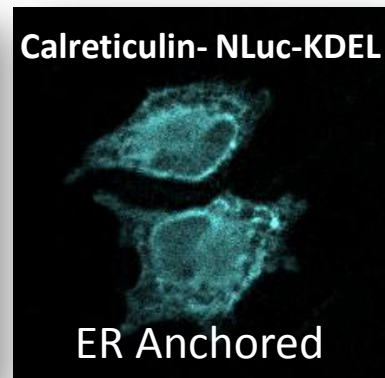
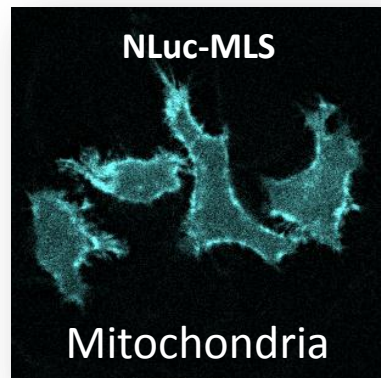
*Olympus  
Product Showcase*



Olympus LV200  
Bioluminescence Imager



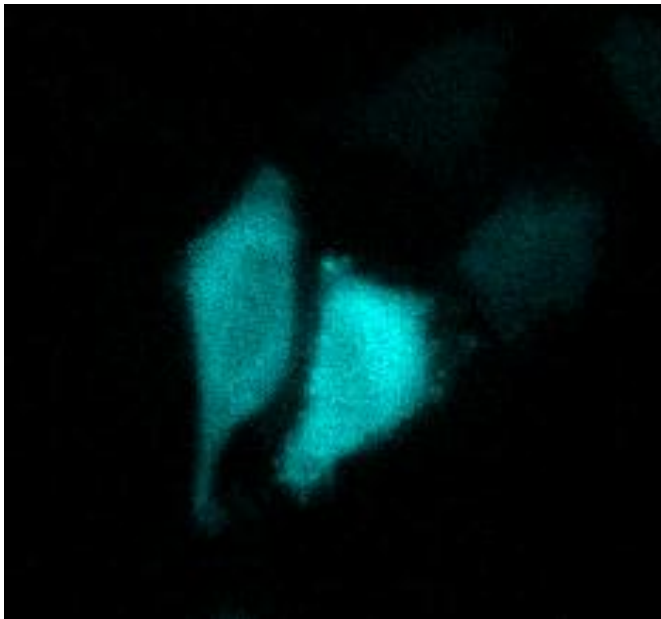
## *NanoLuc™ Fusions can go anywhere...*



NanoLuc™ Luciferase fusions could be a useful tool to investigate cell biology

## ***Bioluminescence imaging of NanoLuc fusion proteins***

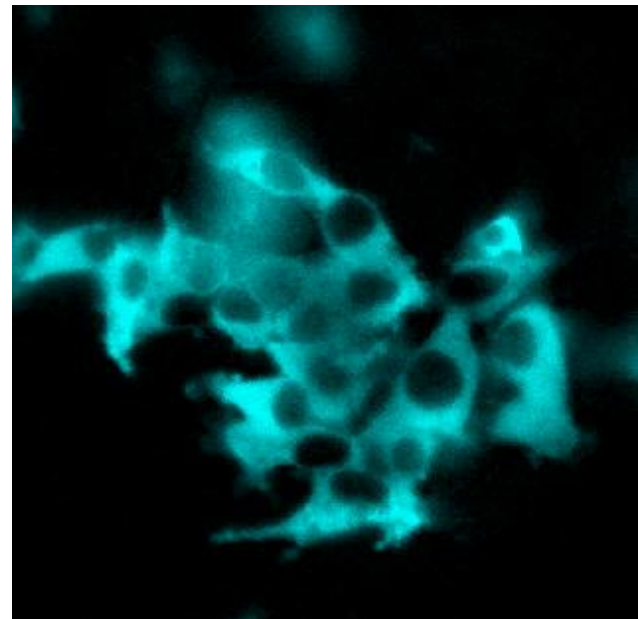
### Glucocorticoid receptor (GR)



*Time lapse: 13 minutes*

Upon treatment with dexamethasone, the receptor dissociates from the Hsp90 complex and translocates to the nucleus.

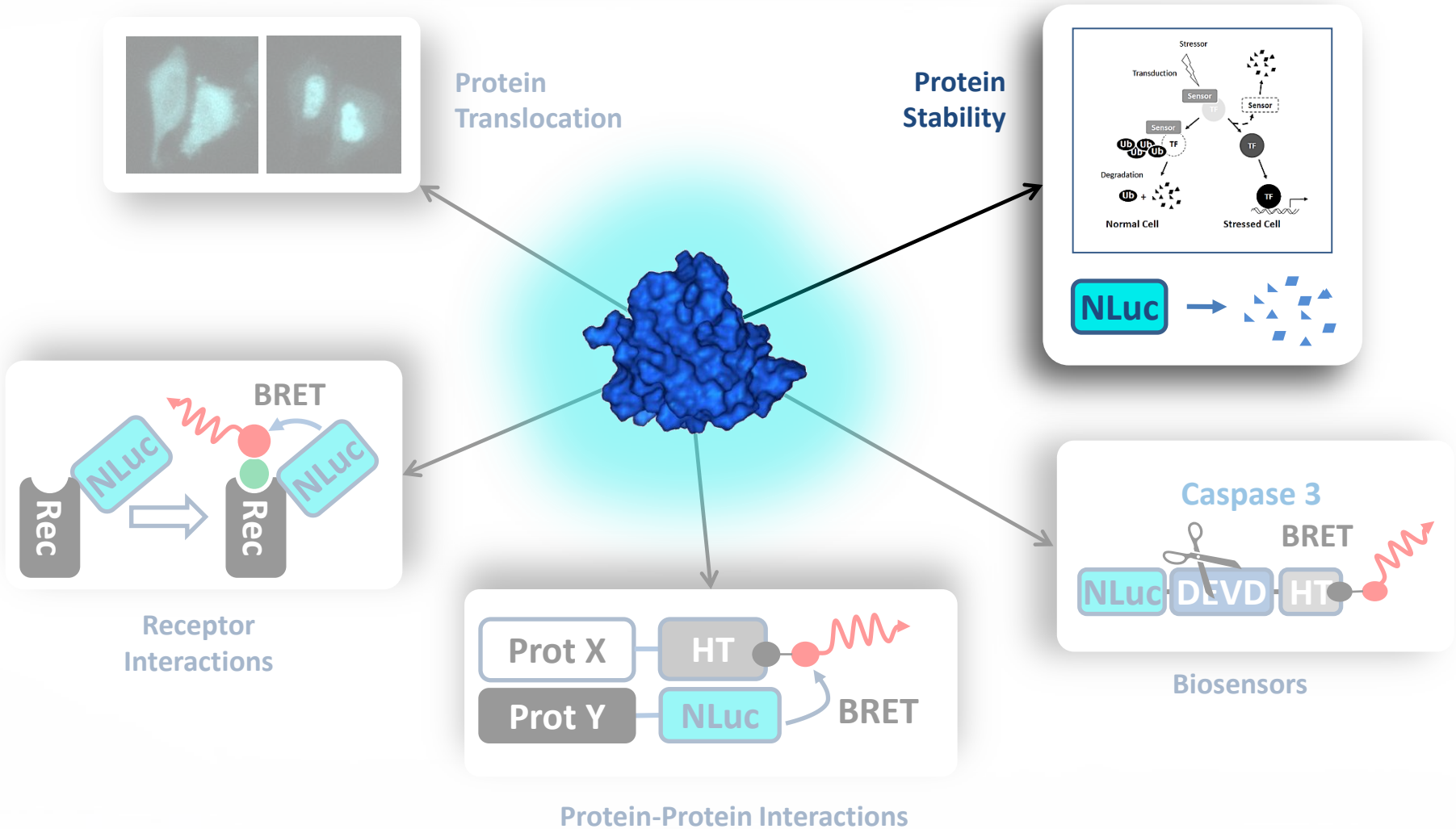
### Protein Kinase C alpha (PKC $\alpha$ )



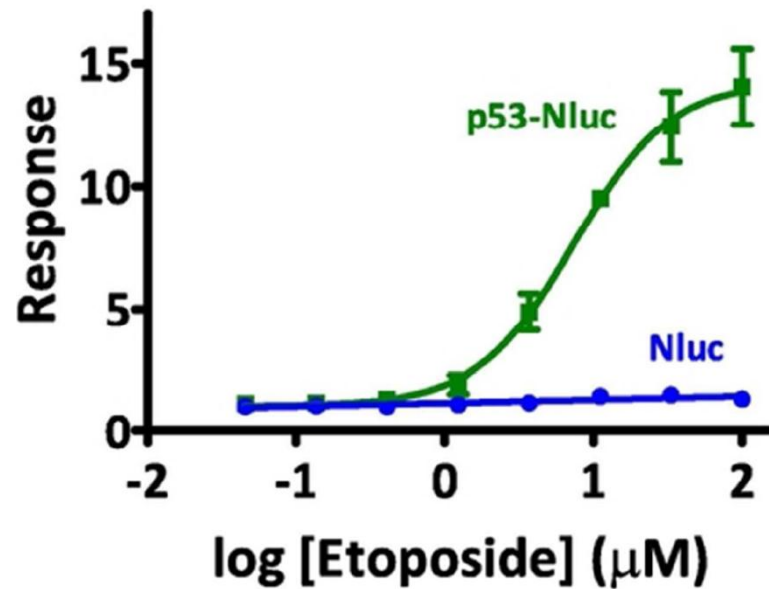
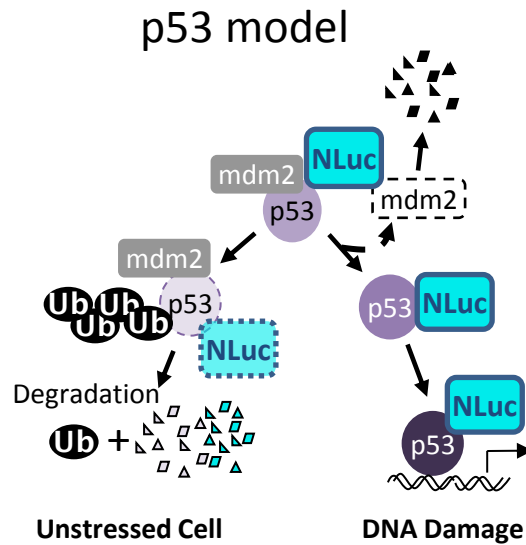
*Time lapse: 20 minutes*

Upon treatment with phorbol esters, PKC $\alpha$  is recruited to the membrane.

# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments

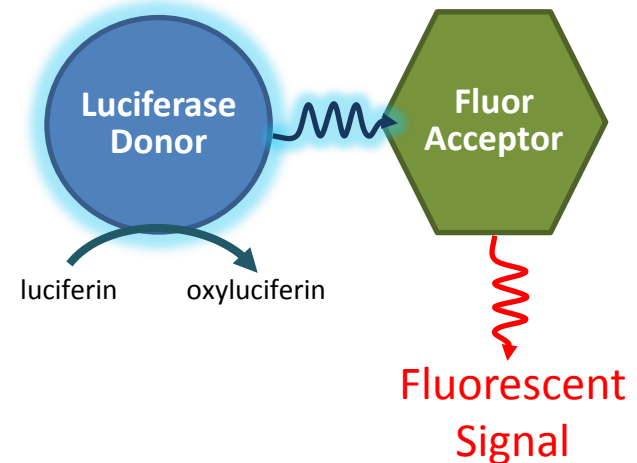
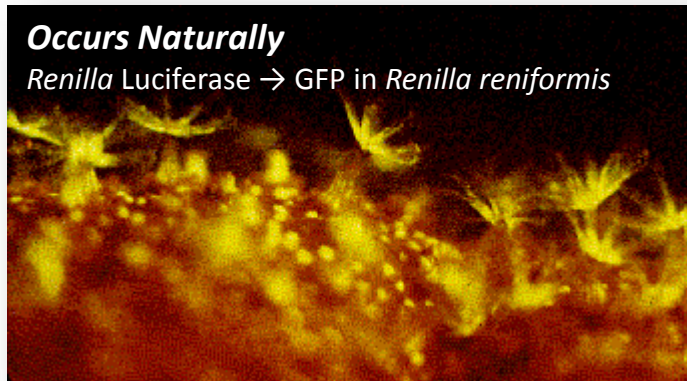


# Monitoring Protein Stability with NanoLuc<sup>TM</sup> Luciferase



The fusion can be used as a probe of stability

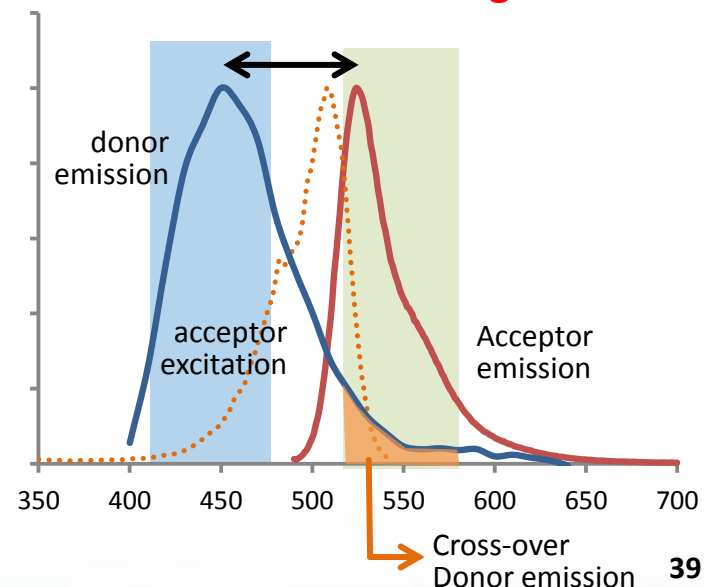
# Bioluminescence Resonance Energy Transfer (BRET)



Important characteristics for BRET applications in research:

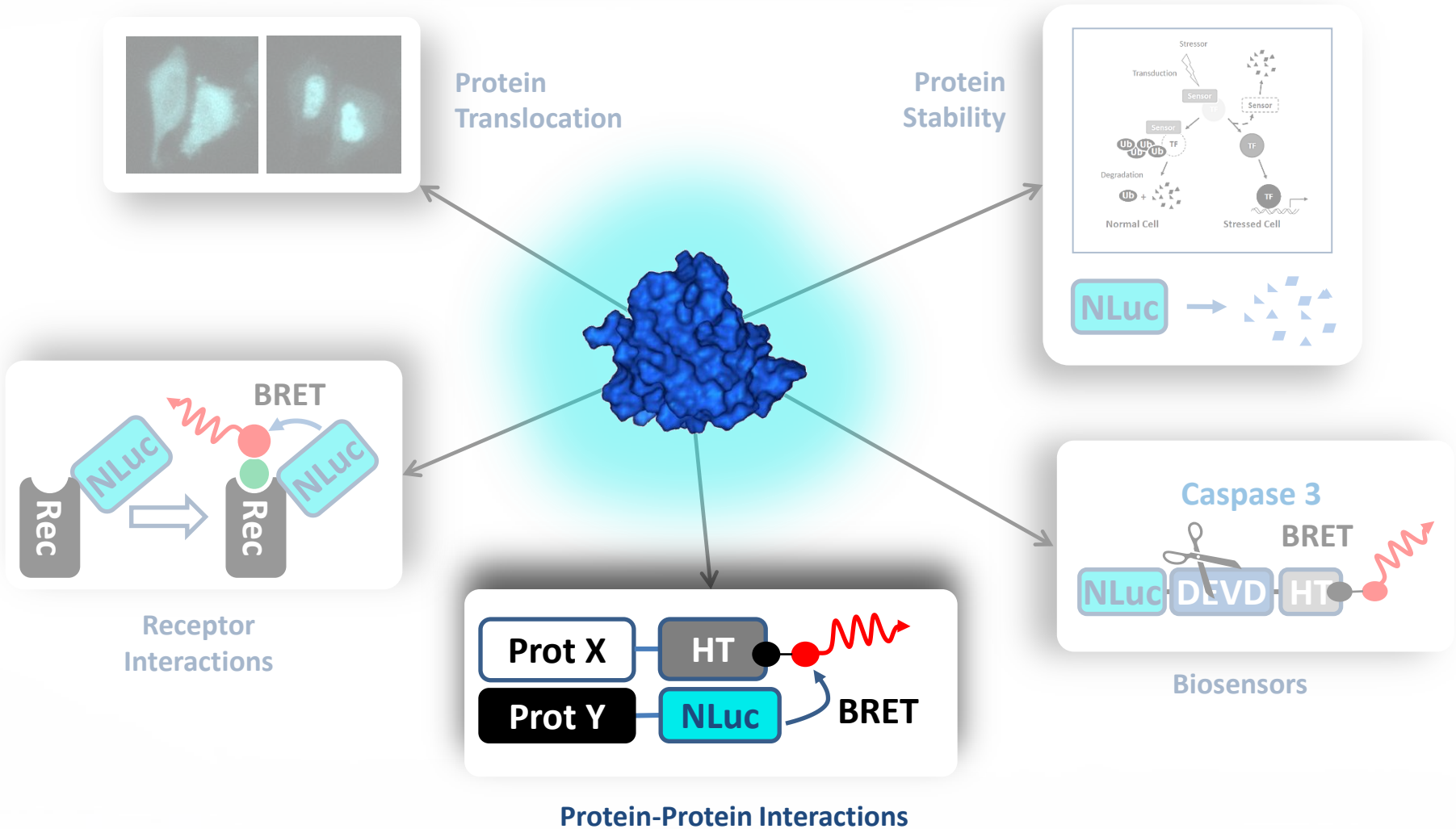
- Donor emission must overlap with acceptor excitation spectra
- Donor & Acceptor must be close (<10nm)
- Acceptor emission must be discernable from Donor emission
- Output intensity is dependent upon donor intensity

*Based on publications from S.S. Gambhir at Stanford*



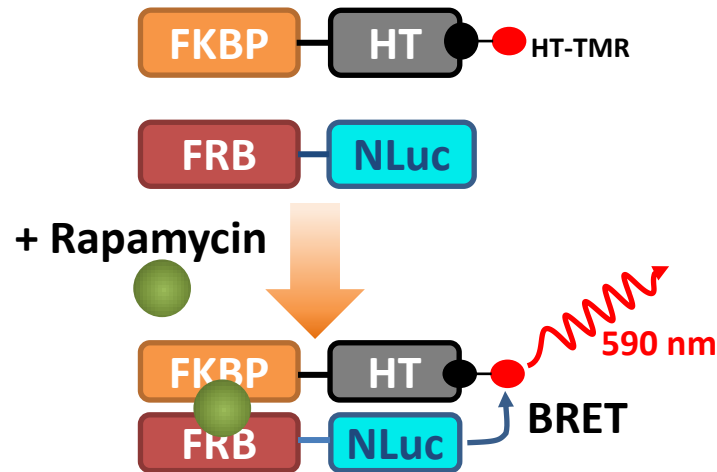


# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments





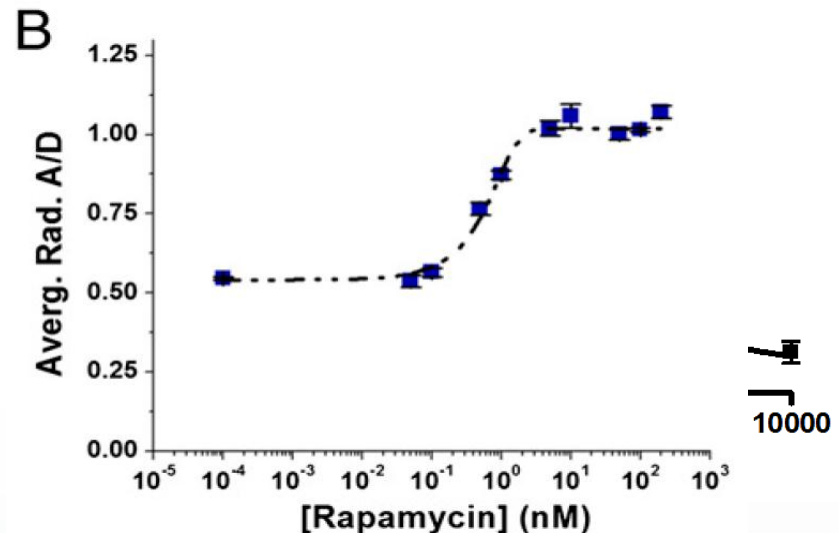
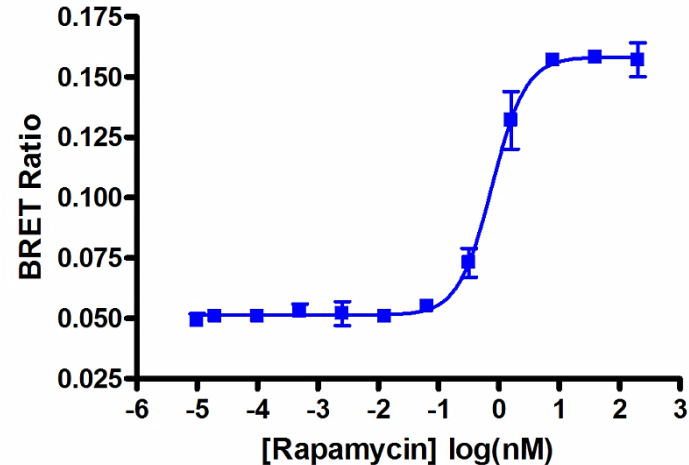
# Can NLuc:HT Pair be used for Protein-Protein BRET?



Same model system used with BRET 6 System

**RLuc8.6** → **TurboFP**

Dragulescu-Andrasi, A., et al  
(2011) *PNAS* **108**, 12060-5.



# *Summary*

- Bioluminescence
- NanoLuc luciferase
- Genetic reporter
  - Dual luciferase assays
  - Animal imaging
- Protein reporter
  - Location
  - Stability
  - Interactions (NanoBRET, NanoBiT)

תודה  
 Dankie Gracias  
 Спасибо شكراً  
 Merci Takk  
 Köszönjük Terima kasih  
 Grazie Dziękujemy Děkojame  
 Ďakujeme Vielen Dank Paldies  
 Kiitos Tänname teid 谢谢  
**Thank You** Tak  
 感謝您 Obrigado Teşekkür Ederiz  
 Σας ευχαριστούμε 감사합니다  
 Bedankt ඔබට  
 Děkuje vám  
 ありがとうございます  
 Tack