

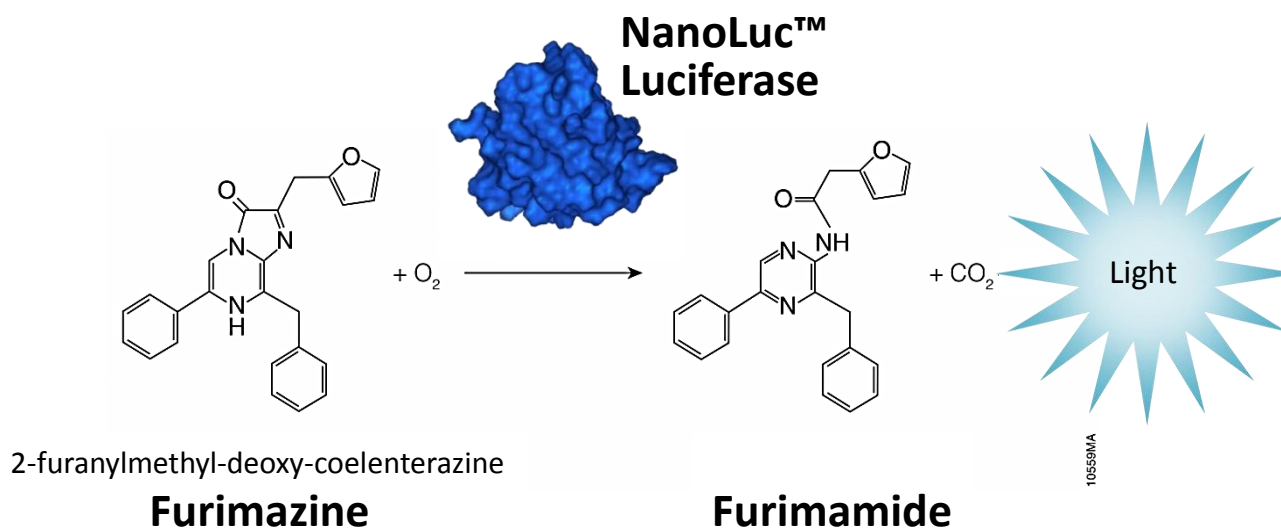
***NanoLuc<sup>®</sup>: A Smaller, Brighter, and  
More Versatile Luciferase Reporter***



**Terry L. Riss, Ph.D.**  
***Senior Product Specialist, Cell Health***  
***Promega Corporation***

## What is NanoLuc™ Luciferase?

*NanoLuc™ (Nluc) is a 19.1 kDa, ATP-independent luciferase that utilizes a novel coelenterazine analog (furimazine) to produce high intensity, glow-type luminescence.*

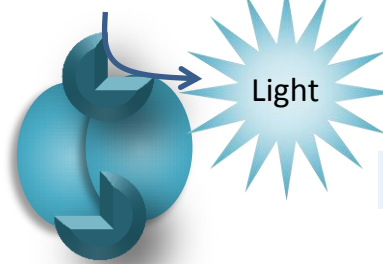


# Evolution of NanoLuc from ocean to lab bench



*Oplophorus gracilirostris* first cataloged in 1881

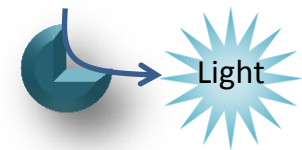
Coelenterazine



130kDa

*Oplophorus* luciferase  
7X brighter than native  
*Renilla* Luciferase  
Shimomura, O., et al. (1978)

Coelenterazine

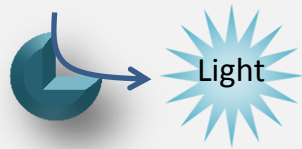


19kOLuc

19kDa subunit is catalytic.  
Light output & stability  
compromised.  
Inouye, S., et al. (2000)

## Promega Advanced Technologies Group

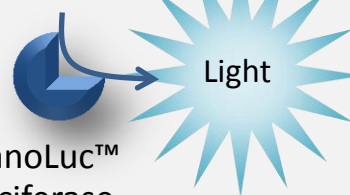
Coelenterazine



19kOLuc

enzyme evolution

Coelenterazine



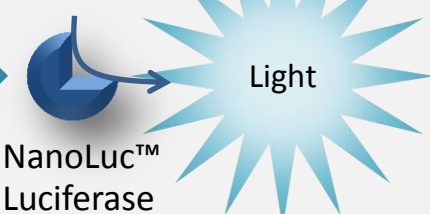
NanoLuc™  
Luciferase

81,000X

Hall, M.P., et al. (2012) ACS Chem Biol 7:1848-1857.

substrate evolution

Furimazine

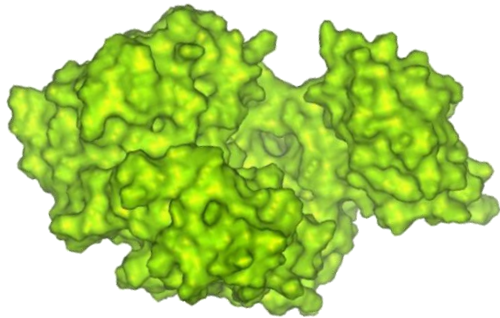


NanoLuc™  
Luciferase

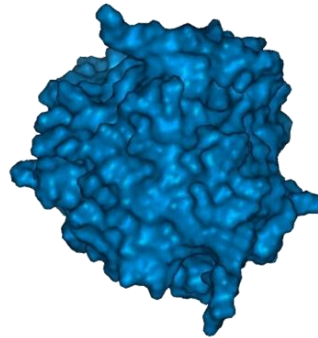
2,500,000X

# NanoLuc™ is very small

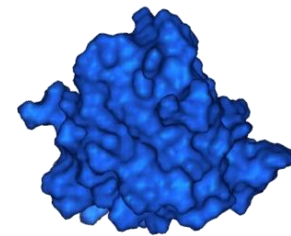
*Firefly (Fluc)*



*Renilla (Rluc)*

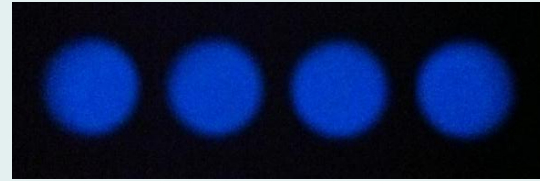


*NanoLuc™ (Nluc)*

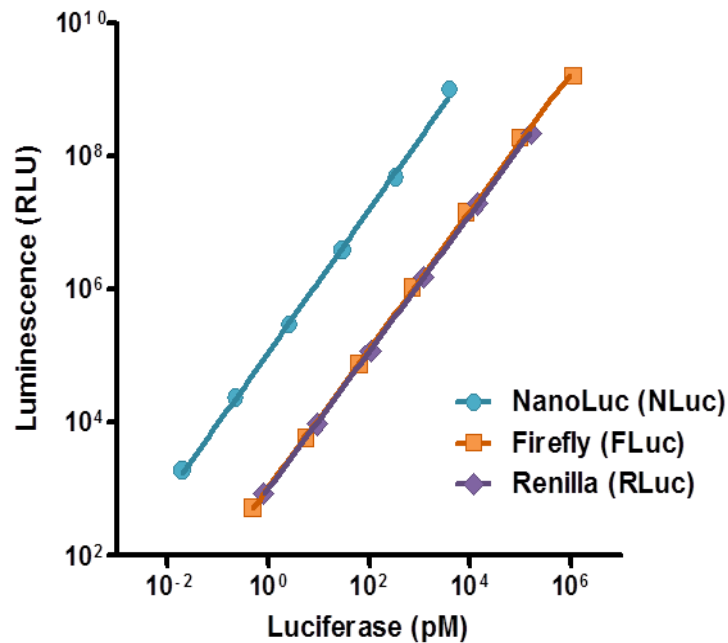


	Amino acids	M.W.	Mol. Vol. Å <sup>3</sup>
<b>Nluc</b>	171	19.1	14
<b>Rluc</b>	312	36.0	32
<b>Fluc</b>	550	60.6	44

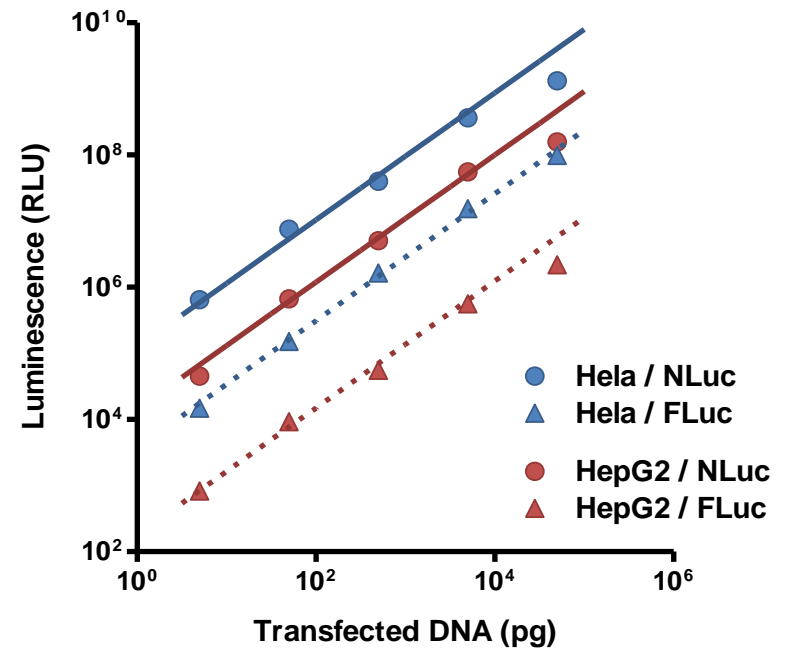
# NanoLuc is very bright



Living HEK293 cells in 96-well plate (50,000 cells per well). Imaged by a hand-held iPhone



Purified enzyme



Expression in mammalian cells

# NanoLuc™ has excellent physical properties



## Thermal stable enzyme

- Retains activity following 30 min incubation at 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

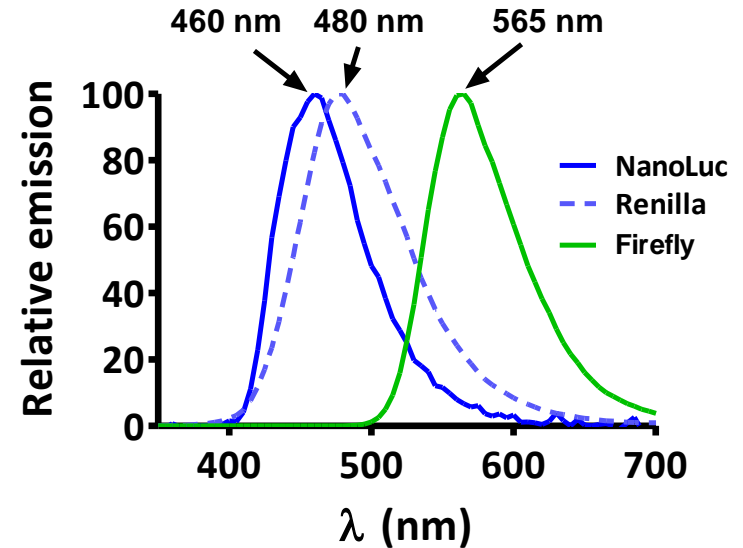
## No post-translational modifications detected in mammalian cells

## No disulfide bonds

- Supports high levels of activity inside living cells

## Uniform distribution in cells

- No apparent compartmental bias in the absence of targeting sequences



unfused NLuc  
Immunofluorescence

# Nano-Glo™ Luciferase Assay Reagent



## Nano-Glo™ Luciferase Assay Reagent:

Furimazine

- Provides maximal brightness

Glow kinetics (no flash reaction)

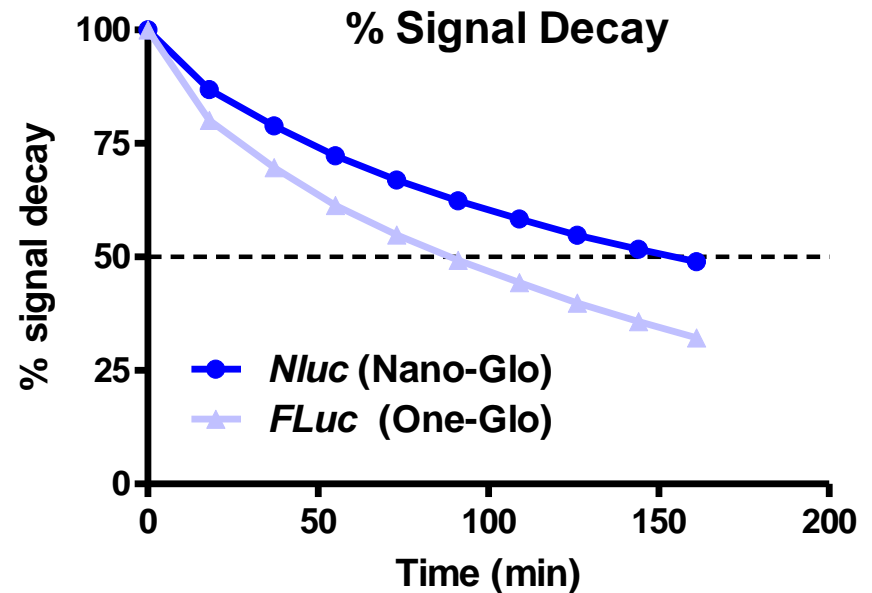
- Half-life routinely >2 hour at room temperature

Low autoluminescence background

- Enhances assay sensitivity

Stable reconstituted reagent:

- ~10% decrease in activity over 8 hrs at RT



## Add-Mix-Measure format like

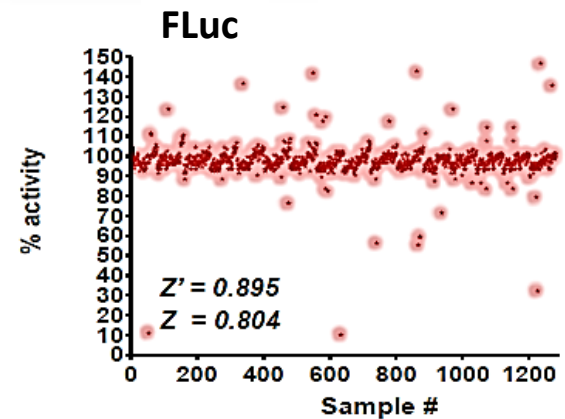
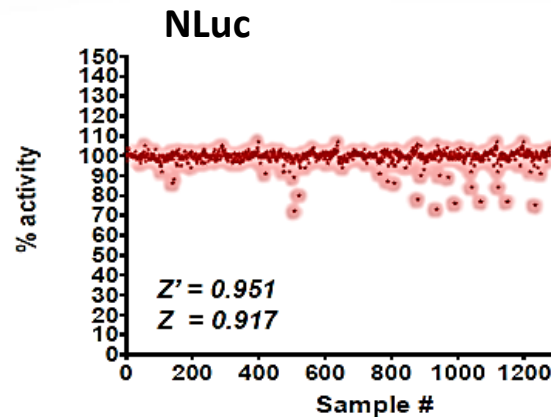
- ONE-Glo™ Luciferase Assay System
- Bright-Glo™ Luciferase Assay System
- Steady-Glo® Luciferase Assay System
- Renilla-Glo™ Luciferase Assay System

# Reduced false hits with NanoLuc<sup>®</sup> Luciferase in HTS



## LOPAC library (Sigma)

- Library of Pharmaceutically Active Compounds
- 1280 compounds
- Small organic ligands w/ well documented pharmacological activities
- Used to screen for non-specific luciferase activity modulators



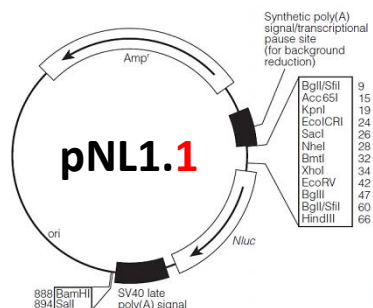
		Level of inhibition			
		≥ 10%	≥ 20%	≥ 30%	≥ 50%
% of library compounds	<b>NanoLuc</b>	<b>1.2%</b>	<b>0.5%</b>	-	-
	<b>Firefly</b>	<b>1.9%</b>	<b>0.7%</b>	<b>0.5%</b>	<b>0.3%</b>

Experimental details: LOPAC library members at 10 μM final concentration; incubation with purified NanoLuc or firefly luciferase for 2 min.; Fluc detection using ONE-Glo™ Luciferase Assay..



# 3 Varieties of NanoLuc<sup>®</sup> Luciferase for you

## Intracellular Formats

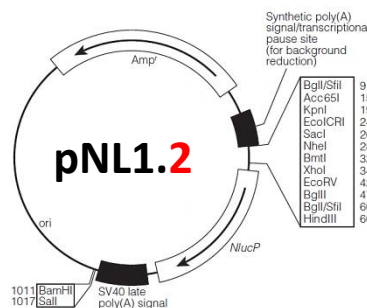


**Nluc (513 bp)**

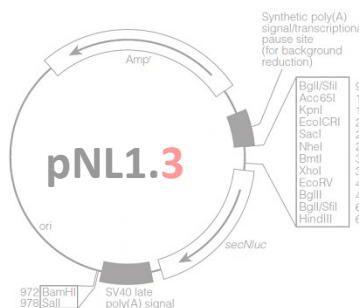
Protein destabilization domain



**NlucP (636 bp)**



## Secretion Format



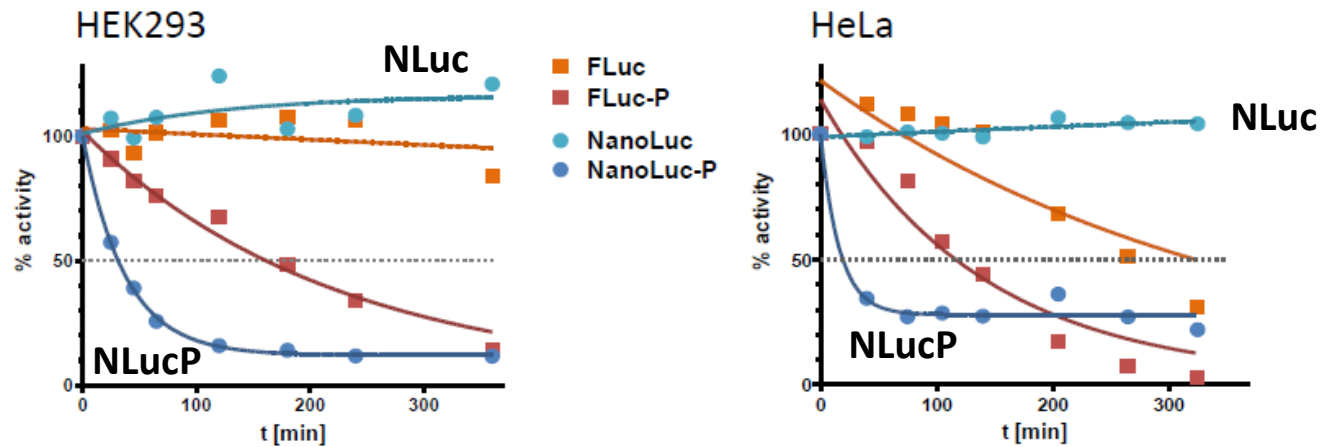
**secNluc (597 bp)**

Secretion signal

# Intracellular stability of NanoLuc™ & Firefly



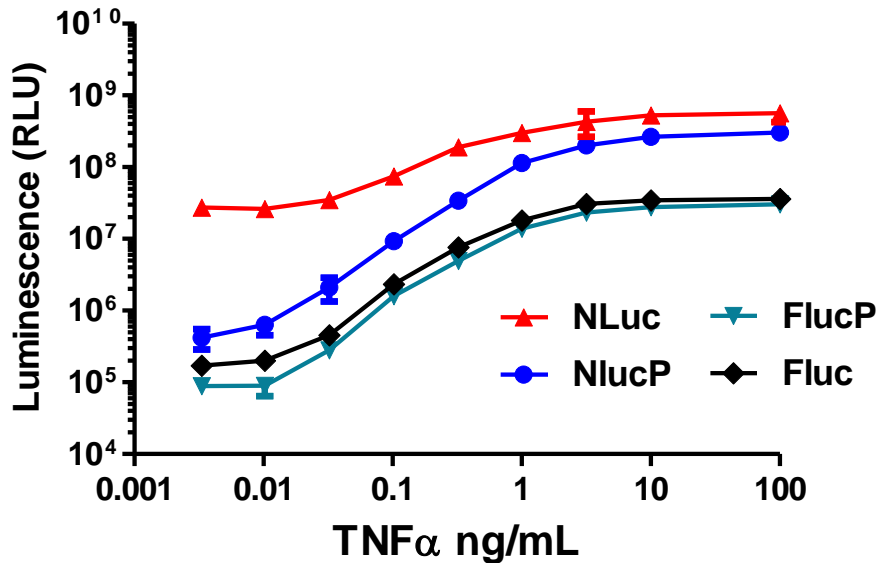
*New protein synthesis blocked by addition of cycloheximide*



Cell line	FLuc	FLucP	NLuc	NLucP
HEK-293	>6 h	2.0 ± 0.4 h	>6 h	18 ± 11 min
HeLa	3.8 ± 1.3 h	1.4 ± 0.2 h	>6 h	20 ± 6 min
U2OS (n=1)	>6 h	2.8 h	>6 h	36 min

**Relative protein stability in cells: NlucP < FlucP < Fluc < Nluc**

# NlucP gives the greatest dynamic response



**Experimental details:** transient transfection of HEK293 cells with NF- $\kappa$ B inducible constructs. rhTNF $\alpha$  treatment for 5 hours.

## Brightness

**NLuc** > **NlucP** > **Fluc** > **FlucP**

(18 experiments)

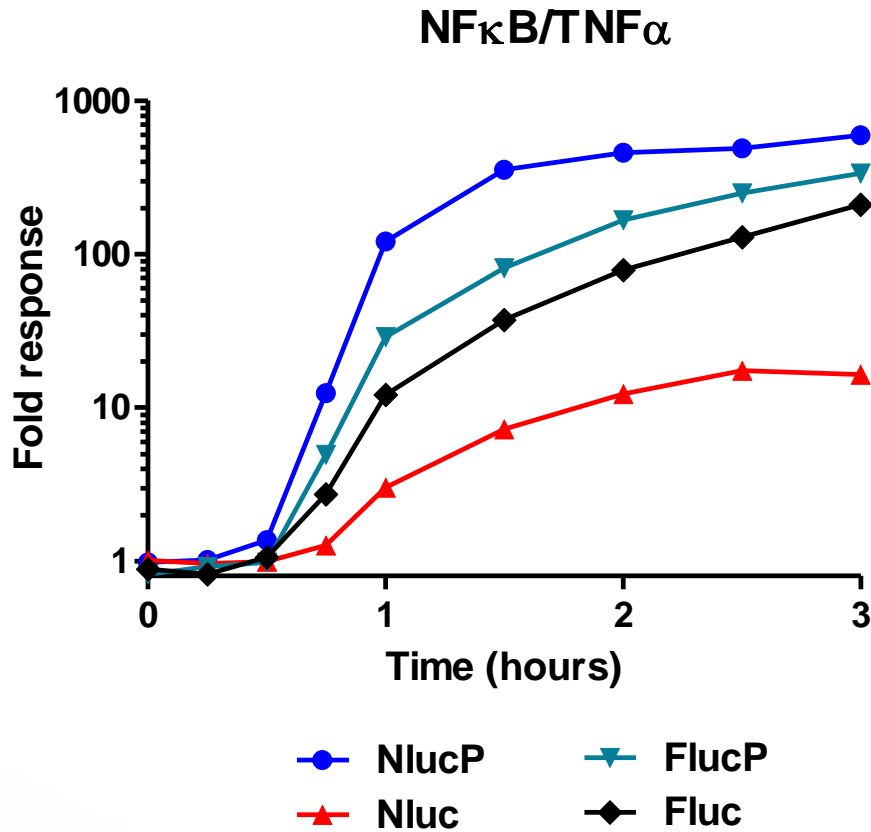
NLuc 13-236 fold brighter than Fluc (79 fold avg.)

NlucP 2-27 fold brighter than FLucP (10 fold avg.)

NLuc 10-78 fold brighter than NlucP (34 fold avg.)

→ Very similar pharmacology/EC50s

# NlucP responds earliest to stimuli



## Relative Response

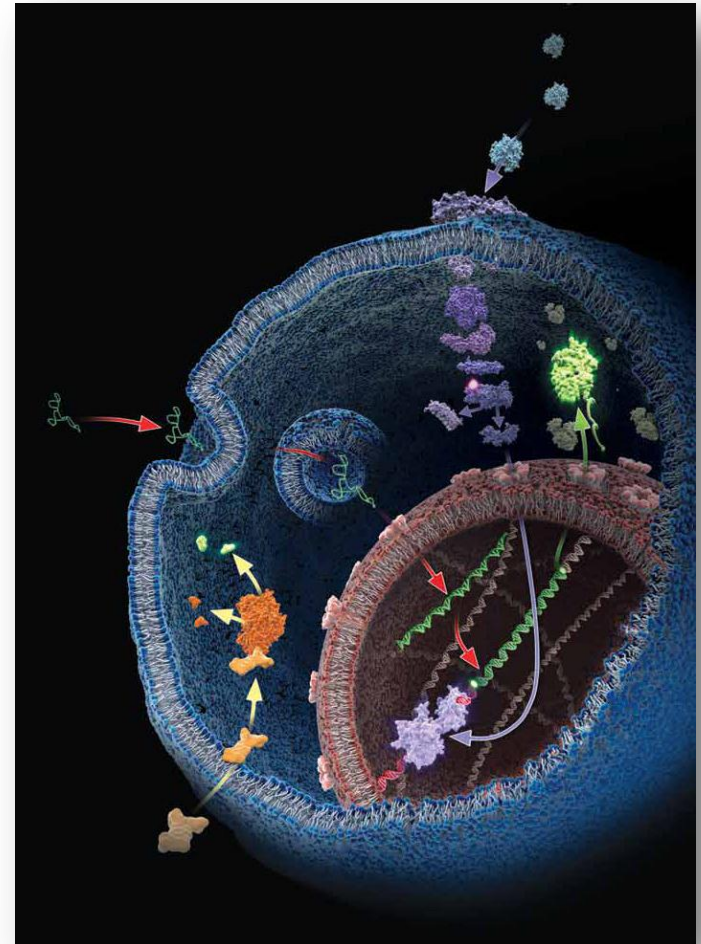
**NlucP** > **FlucP** > **Fluc** > **Nluc**

**Experimental details:** transient transfection of HEK293 cells with NF $\kappa$ B inducible constructs; addition of 100 ng/ml rhTNF $\alpha$  at time zero.

# Summary: *NanoLuc Luciferase as an intracellular reporter*

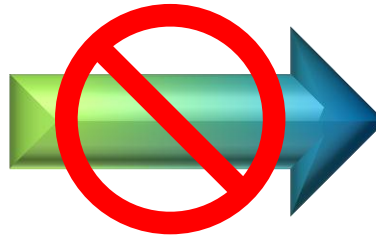
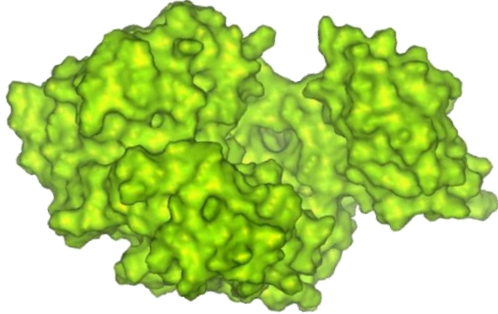


- ✓ NlucP for a faster response
- ✓ NlucP for greatest dynamic range
- ✓ NlucP for measuring weak responses
- ✓ Nluc where maximum brightness is needed.



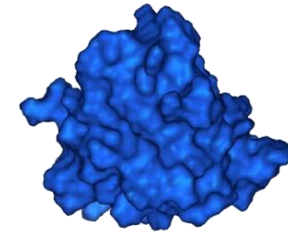
# Should I switch from Firefly to NanoLuc™ Luciferase?

*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**

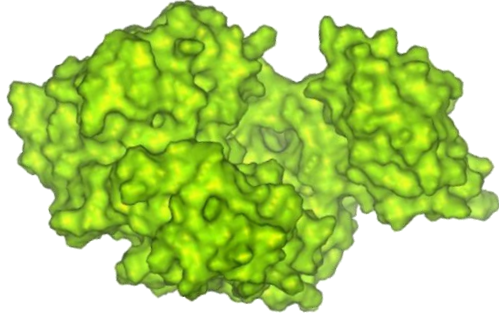
- ✓ Excellent signal:background
- ✓ Excellent dynamic range

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

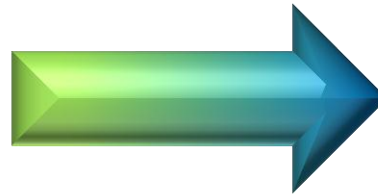
ARE	HSE	ISRE	STAT5	SRE
p53	HRE	SIE	NFAT	SRF
ATF6	XRE	SBE	CRE	
MRE	AP1	TCF-LEF	NF-κB	

# Should I switch from Firefly to NanoLuc™ Luciferase?

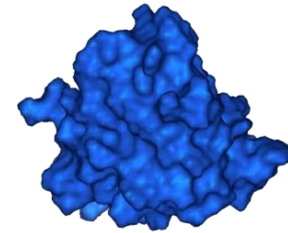
*Firefly (Fluc)*



*Yes, if ...*



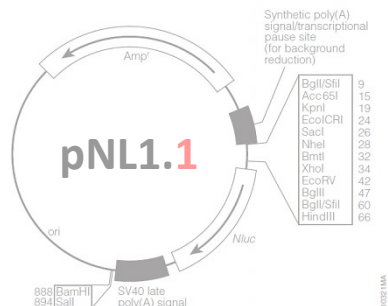
*NanoLuc™ (Nluc)*



- ✓ **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

# 3 Varieties of NanoLuc™ Luciferase for you

## Intracellular Formats

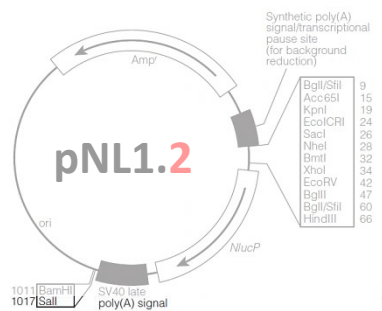


*Nluc* (513 bp)

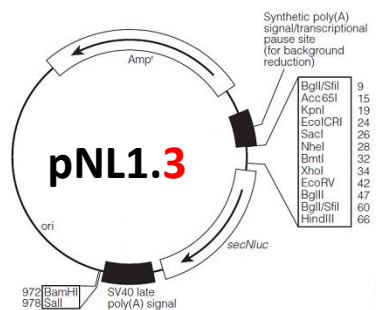
Protein destabilization domain



*NlucP* (636 bp)



## Secretion Format

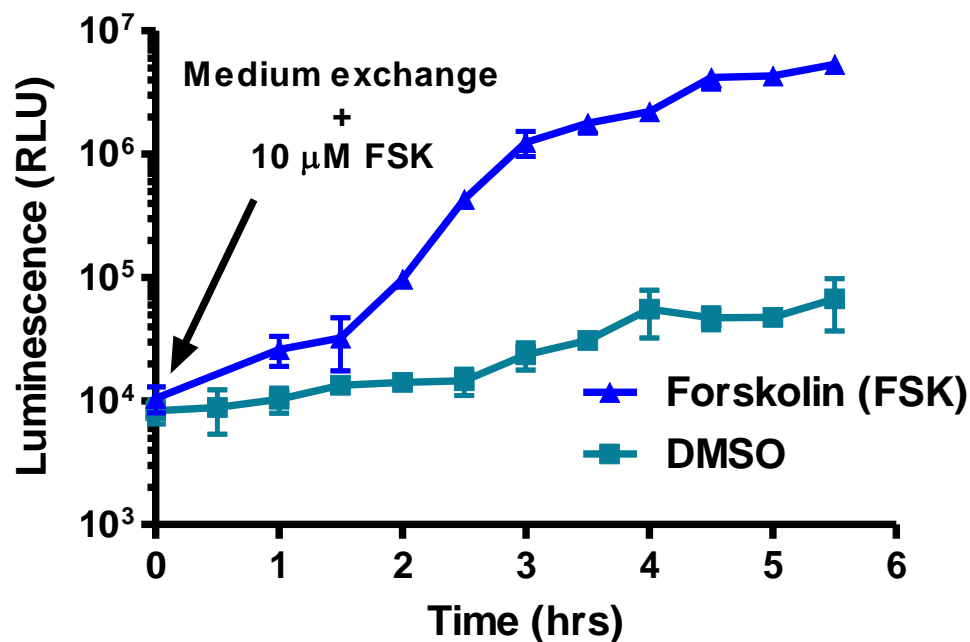


*secNluc* (597 bp)

Secretion signal

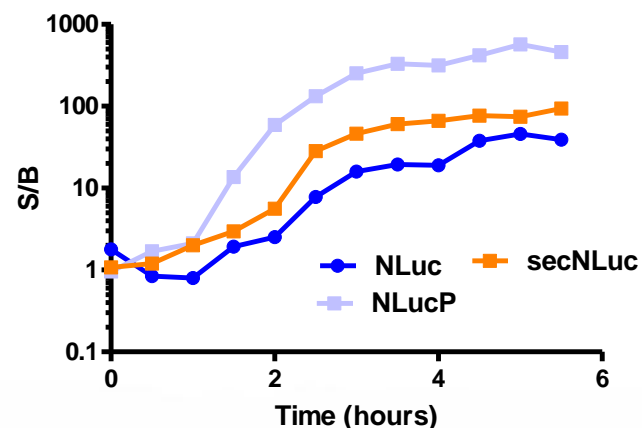


## Secretion based format using secNluc

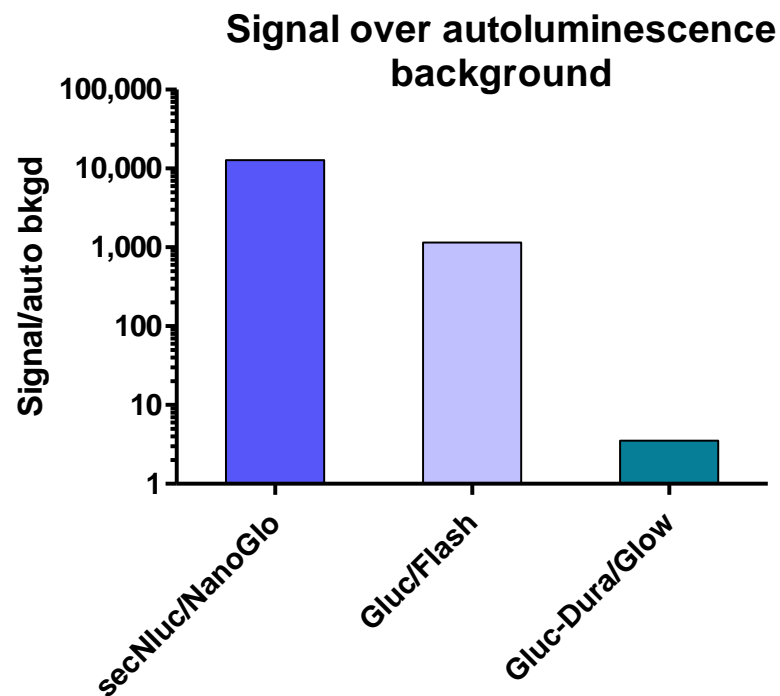
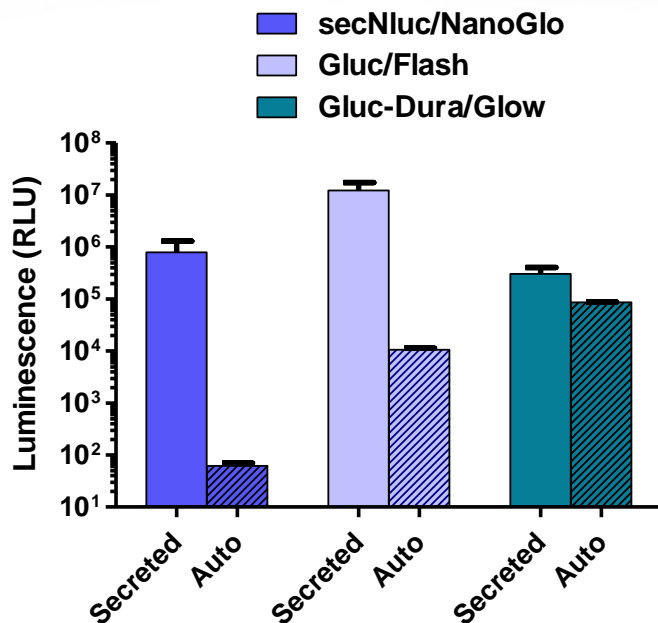


**Experimental details:** transient transfection of HEK293 cells with CREB inducible construct; addition of 10  $\mu$ M forskolin at time zero.

- Sample medium at multiple time points without cell lysis
  - Kinetic studies from the same set of wells
- Half-life of secNluc protein > 4 days at 37°C in medium
- Response dynamics similar to unfused Nluc
- Similar pharmacology vs. Nluc/NlucP



## Gluc kits: bright, but high autoluminescence background



**Gluc kits: high background limits sensitivity & dynamic range**

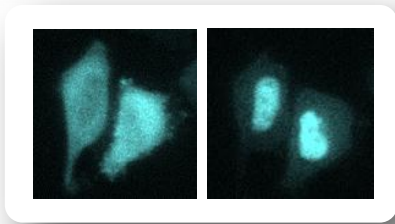
**Experimental details:** HepG2 cells (DMEM +10% FBS) transiently transfected w/CMV promoter constructs; removal of aliquots after 22 hrs; n = 12 per treatment.

# ***NanoLuc™ Luciferase as a protein function probe***

Applications of full-length NanoLuc™ Luciferase.

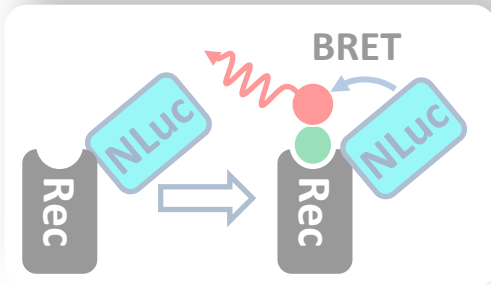
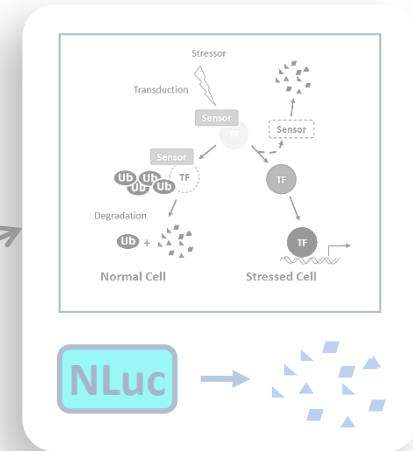


# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments

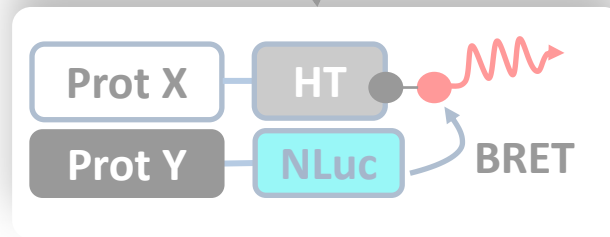


Protein  
Translocation

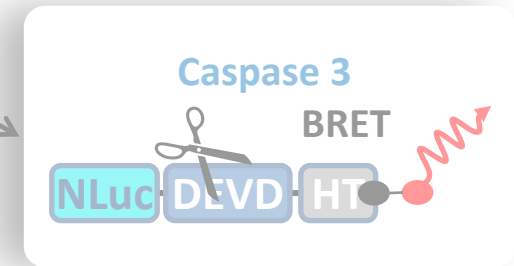
Protein  
Stability



Receptor  
Interactions



Protein-Protein Interactions



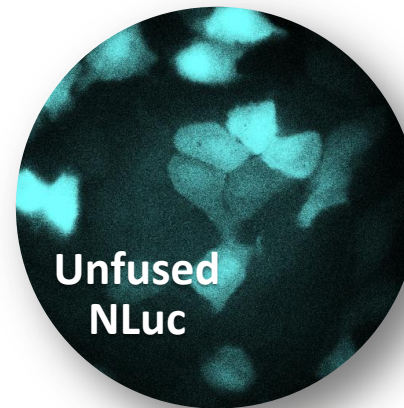
Biosensors

# NanoLuc™ Luciferase excels in bioluminescent imaging applications



Nluc brightness leads to short exposure times:

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



*NanoLuc & LV200 featured @ASCB 2012 Olympus Product Showcase*

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

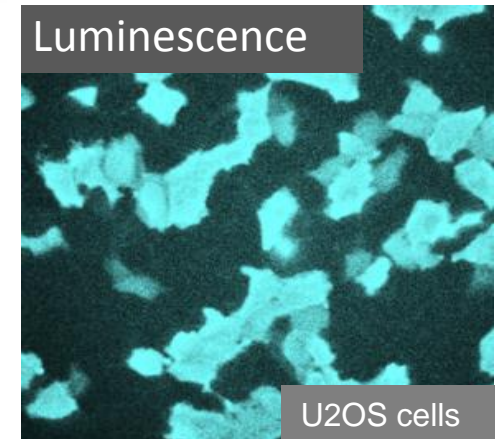
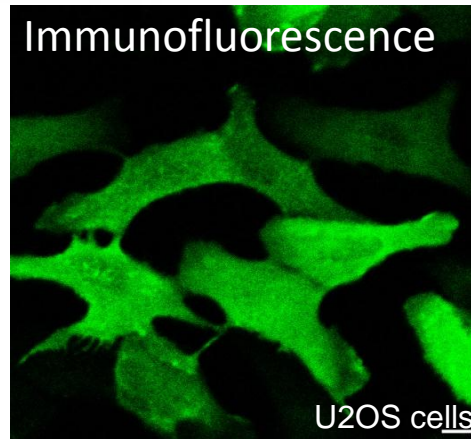


Olympus LV200  
Bioluminescence Imager

# NanoLuc is well behaved in mammalian cells

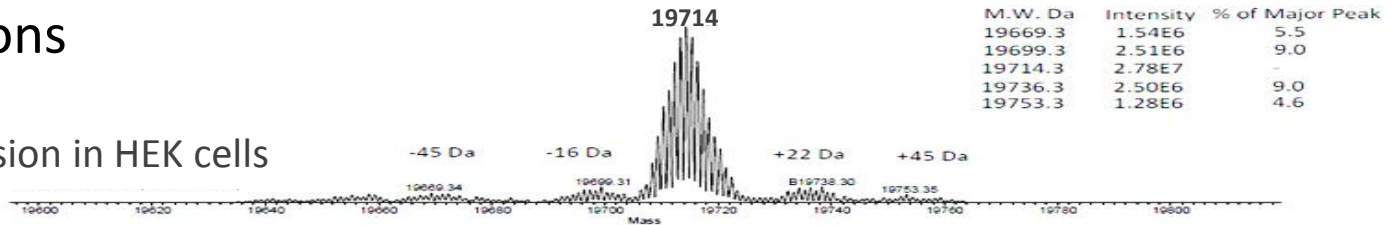


- Uniform intracellular distribution

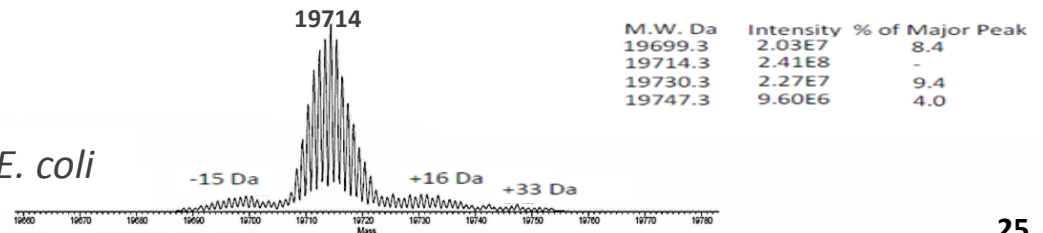


- No post-translational modifications

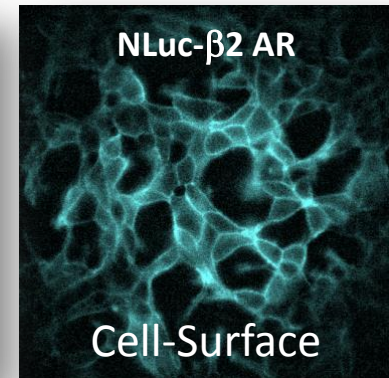
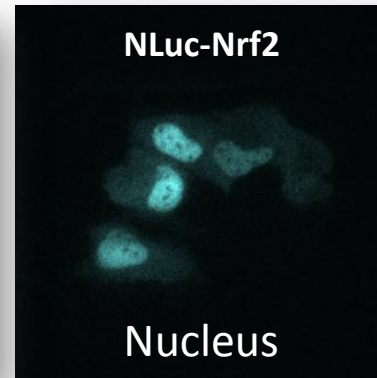
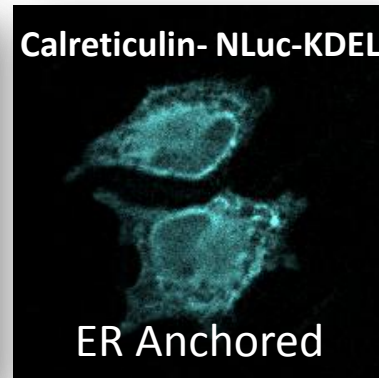
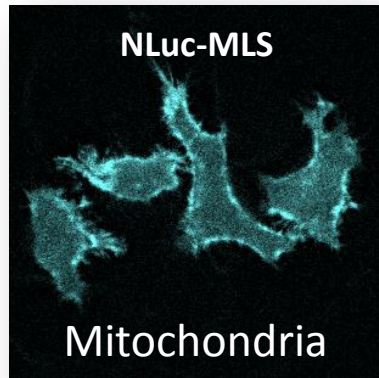
Expression in HEK cells



Expression in *E. coli*



# *NanoLuc™ Fusions can be designed to go anywhere*

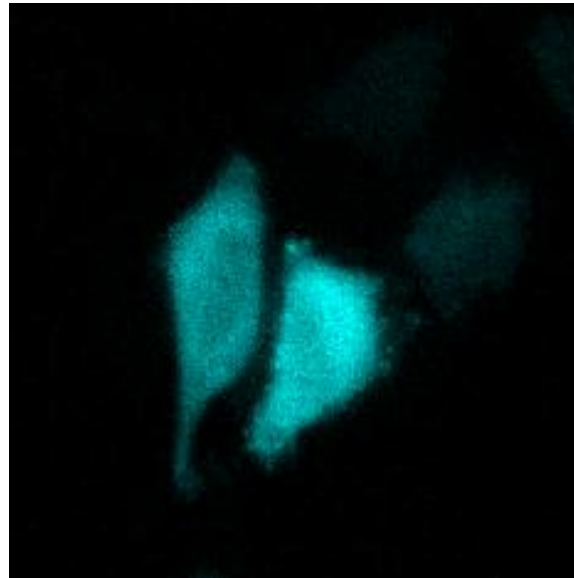


NanoLuc™ Luciferase fusions could be a useful tool to investigate cell biology

# ***Bioluminescence imaging of protein translocation***



cytoplasm → nucleus



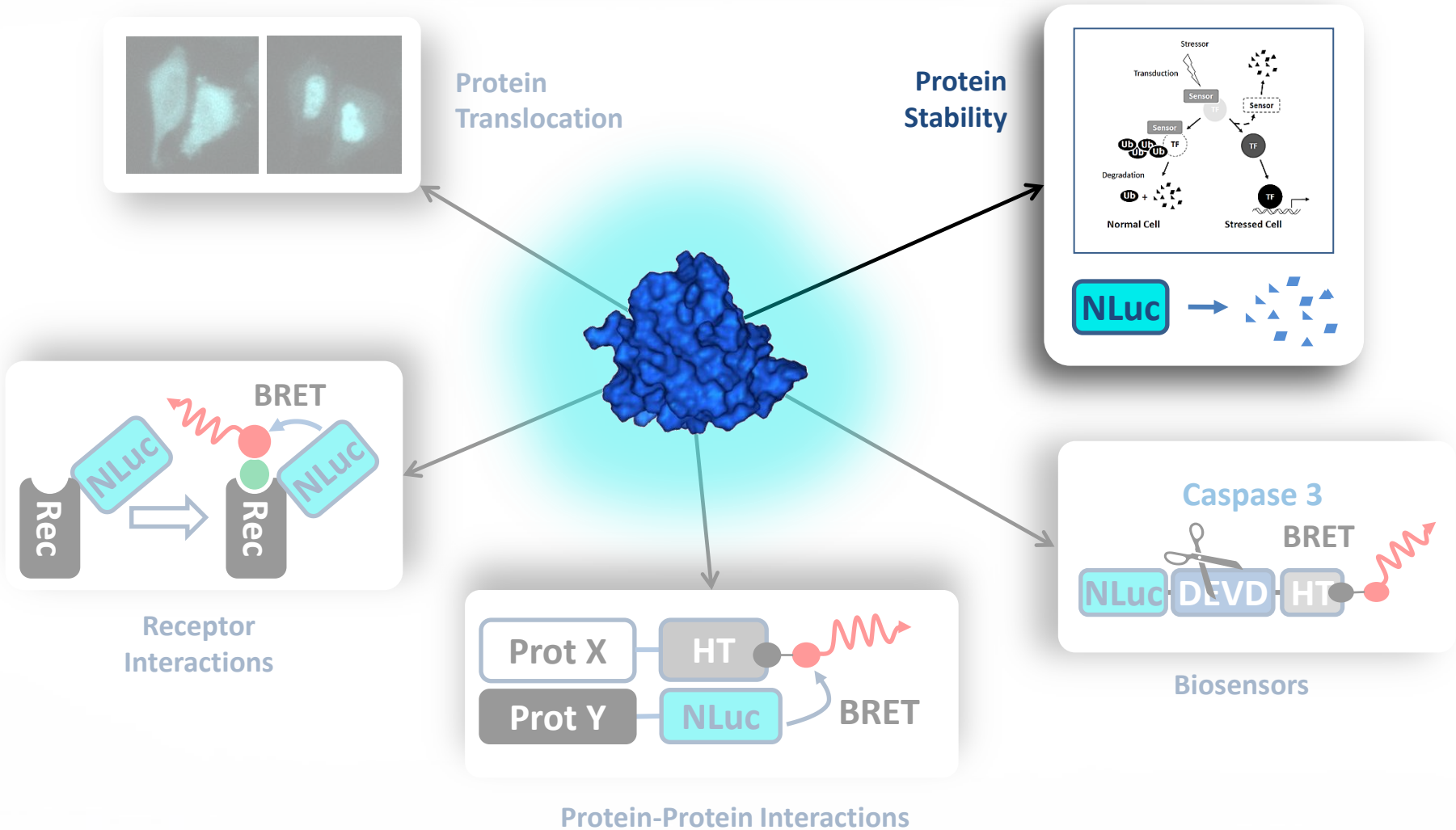
*Time lapse: 13 minutes*

**NanoLuc fusion to  
Glucocorticoid Receptor**

HeLa cells; 500nM dexamethasone treatment

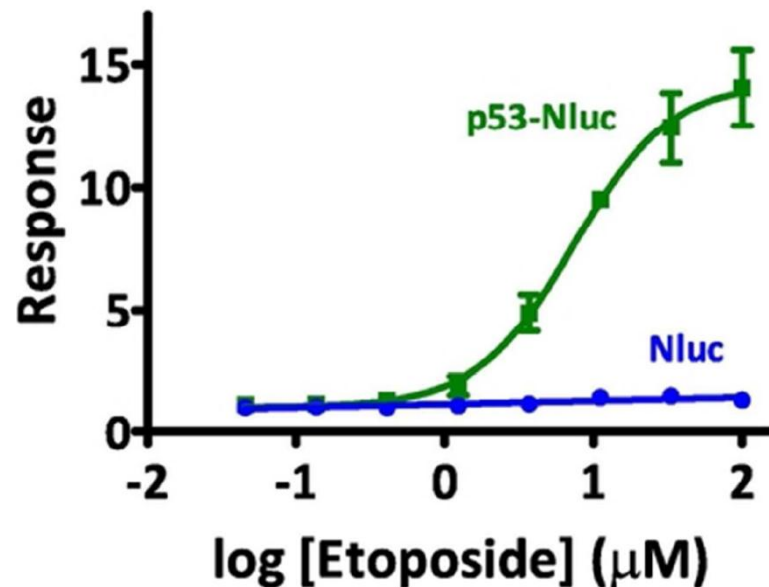
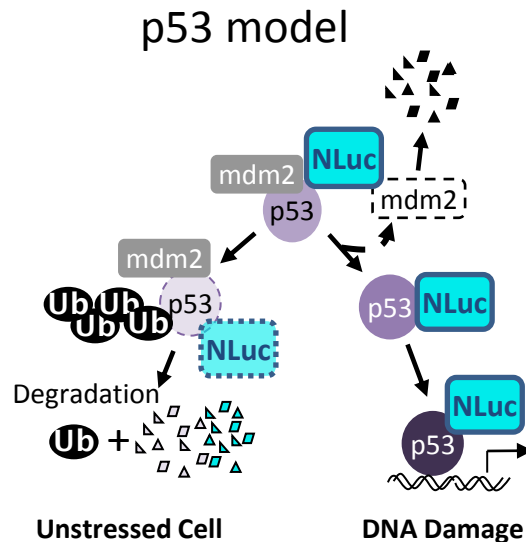


# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments



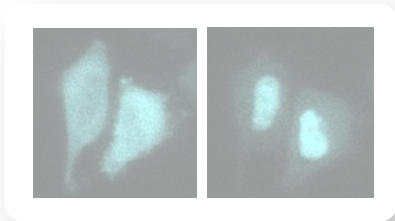
# Monitoring Protein Stability with NanoLuc™ Luciferase

*NanoLuc™ Luciferase can be added to a protein as a probe for protein stability.*



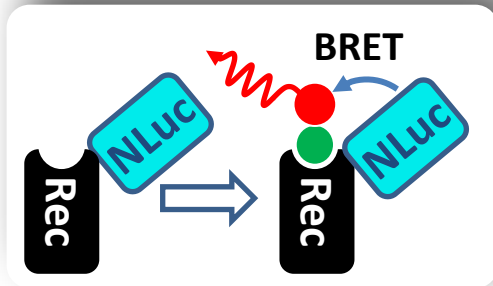
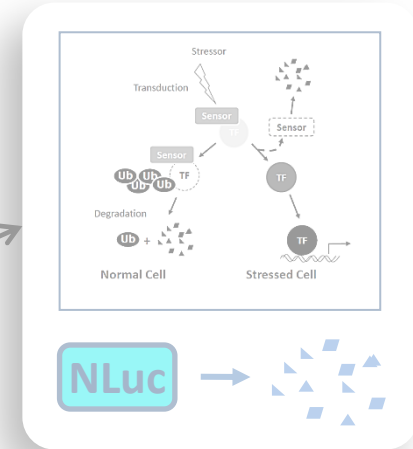
The fusion can be used as a probe of stability

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

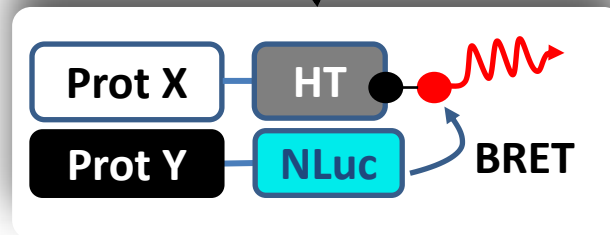


Protein  
Translocation

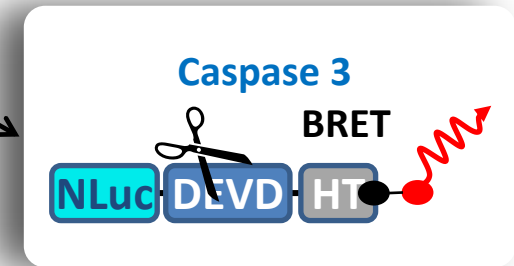
Protein  
Stability



Receptor  
Interactions



Protein-Protein Interactions

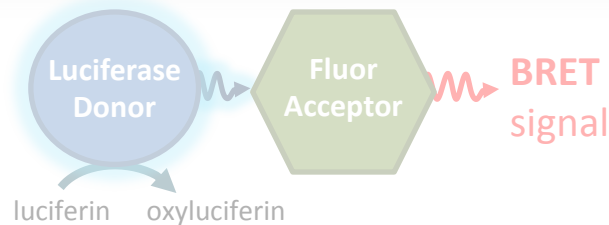
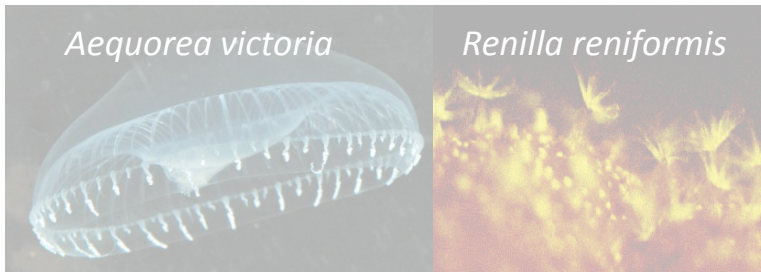


Biosensors

# Bioluminescence Resonance Energy Transfer (BRET)

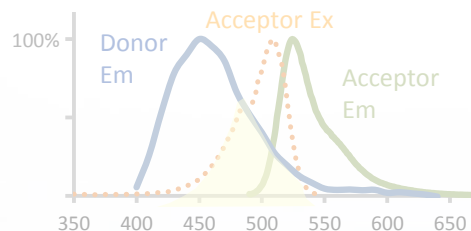


## ...in Nature

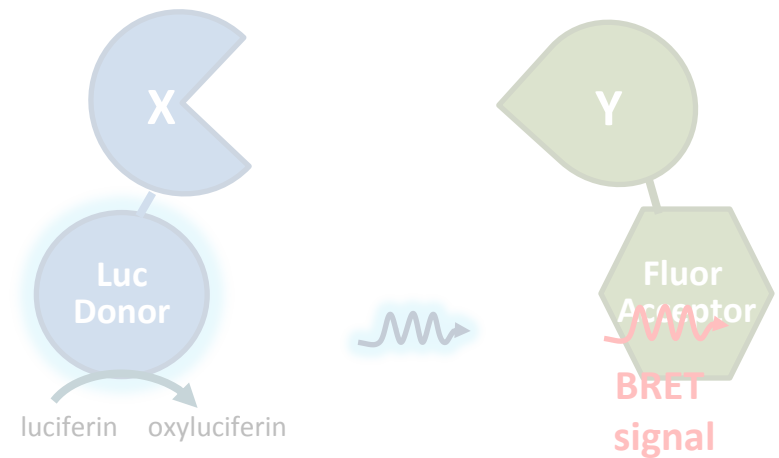
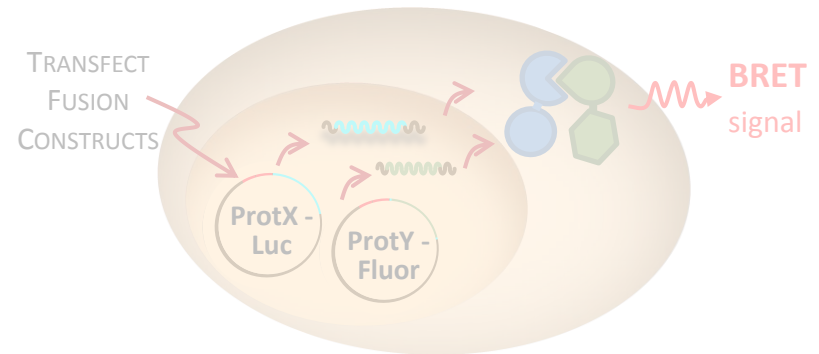


### Requirements for BRET:

- Donor & Acceptor must be close (10-100 Å)
- Donor emission spectrum must overlap with Acceptor excitation spectrum



## ...in the Lab



first described by C.H. Johnson & colleagues in 1999 **32**

# Could NanoLuc™ work better as a BRET donor?



## ***BRET-beneficial properties of NanoLuc Luciferase:***

### **100x brighter than Rluc...**

- ✓ less spectral overlap of Donor & Acceptor needed
- ✓ Better S:B and dynamic range

### **Blue-shifted emission...**

- ✓ Lots of space on the spectrum for the Fluor Acceptor

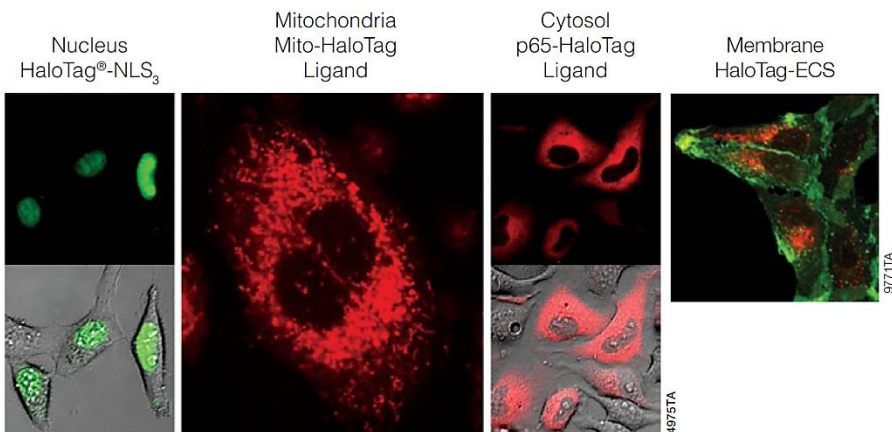
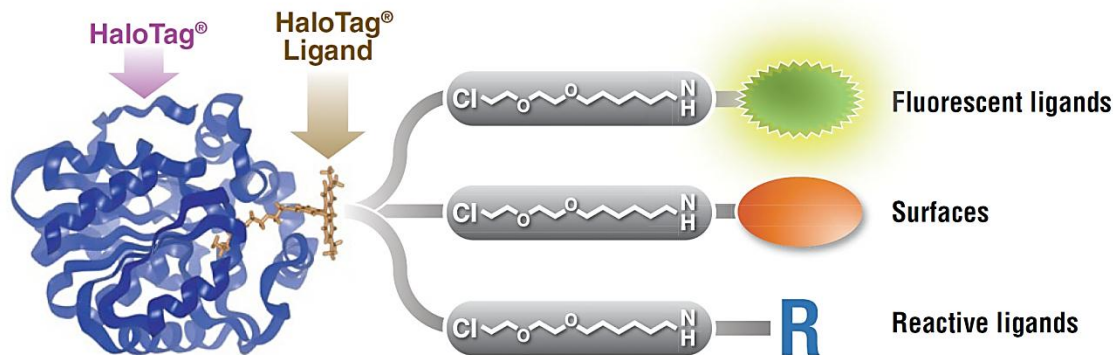
*Nluc 460nm; Rluc 480nm; Fluc 560nm*

# We have a potential acceptor fusion protein:

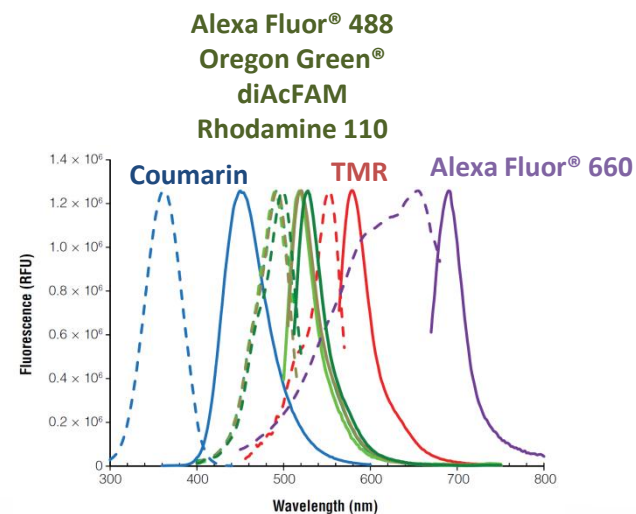
## HaloTag®

34.1kDa protein engineered from halophilic bacterial hydrolase

- Forms **covalent attachment** to functional ligand
- Add ligand to cells expressing HaloTag fusion



Goes anywhere in the cell

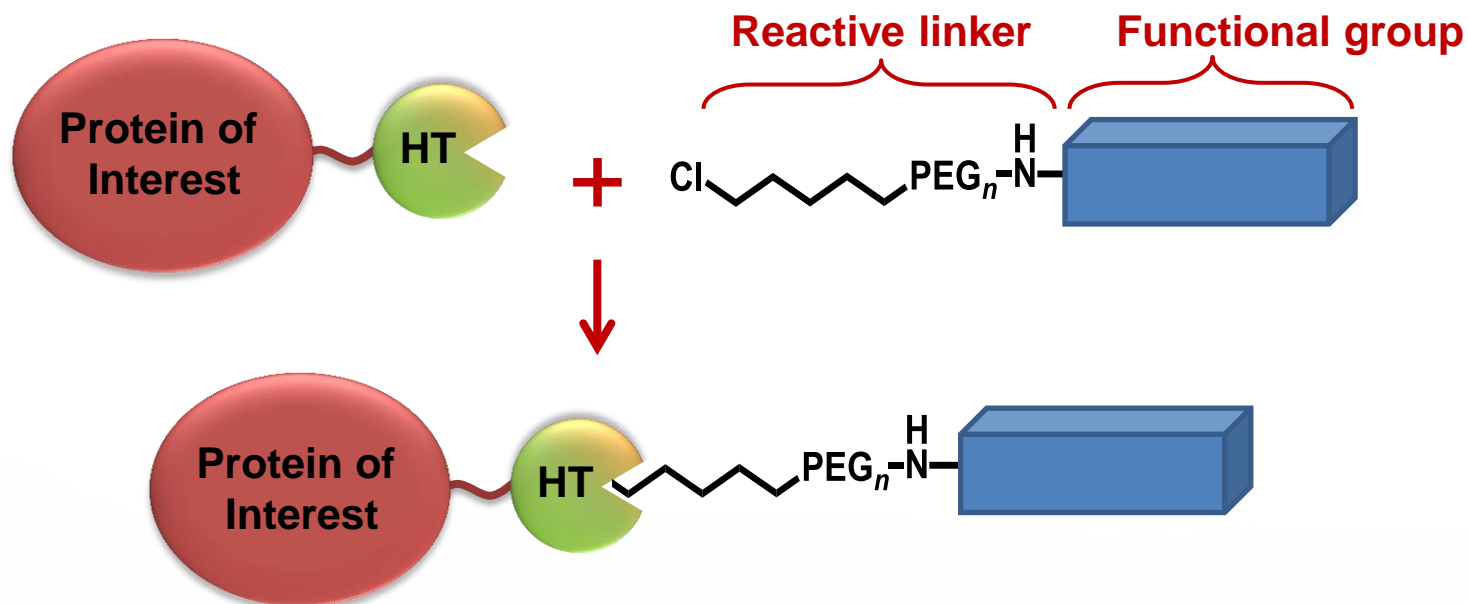


Variety of fluors ready-to-use

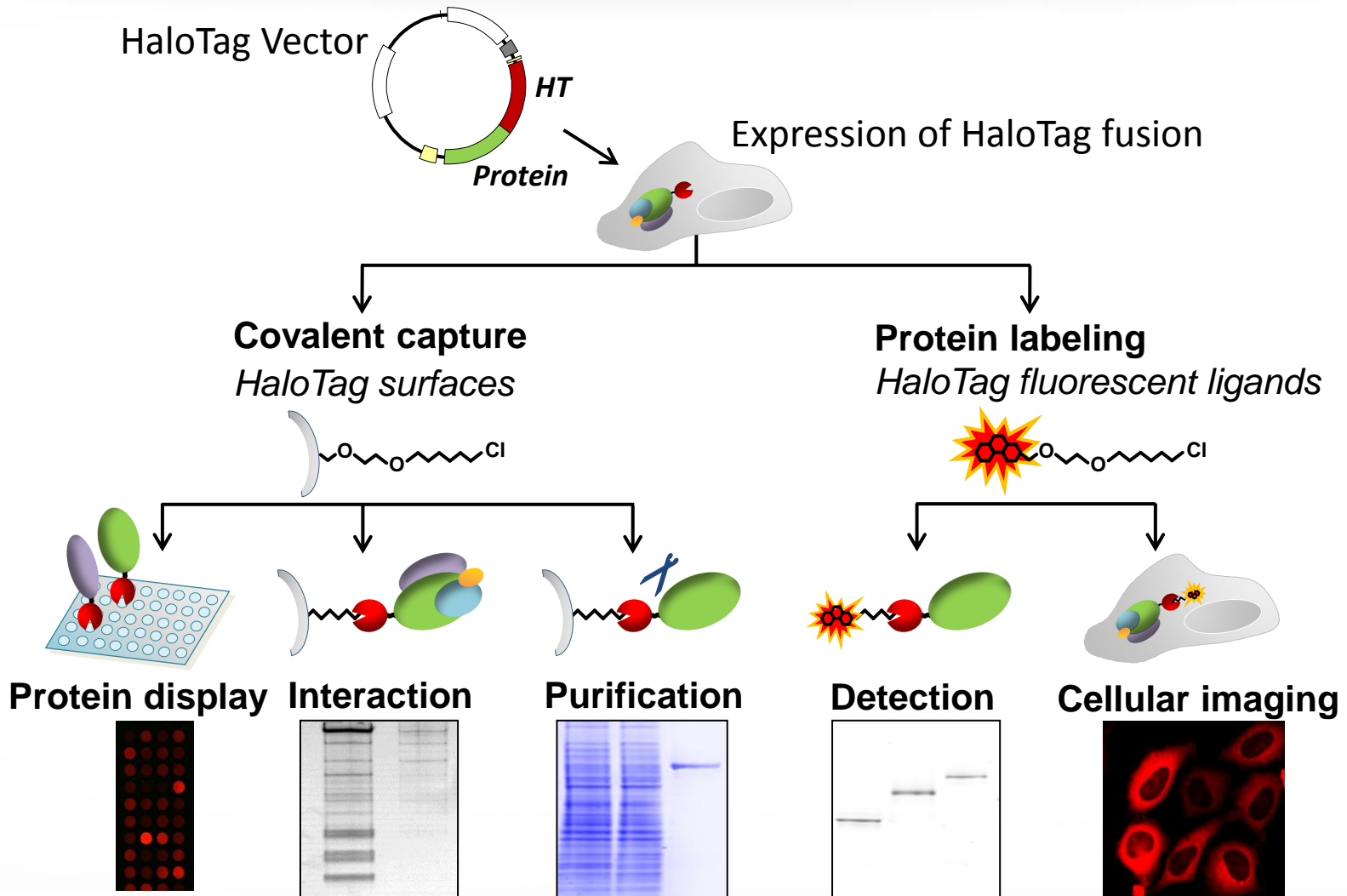
# HaloTag technology – affinity tag alternative

Protein-mediated covalent attachment:

- *Strong binding – covalent bond is essentially irreversible*
- *High specificity – modified catalytic mechanism*
- *High binding rate – optimized protein structure*

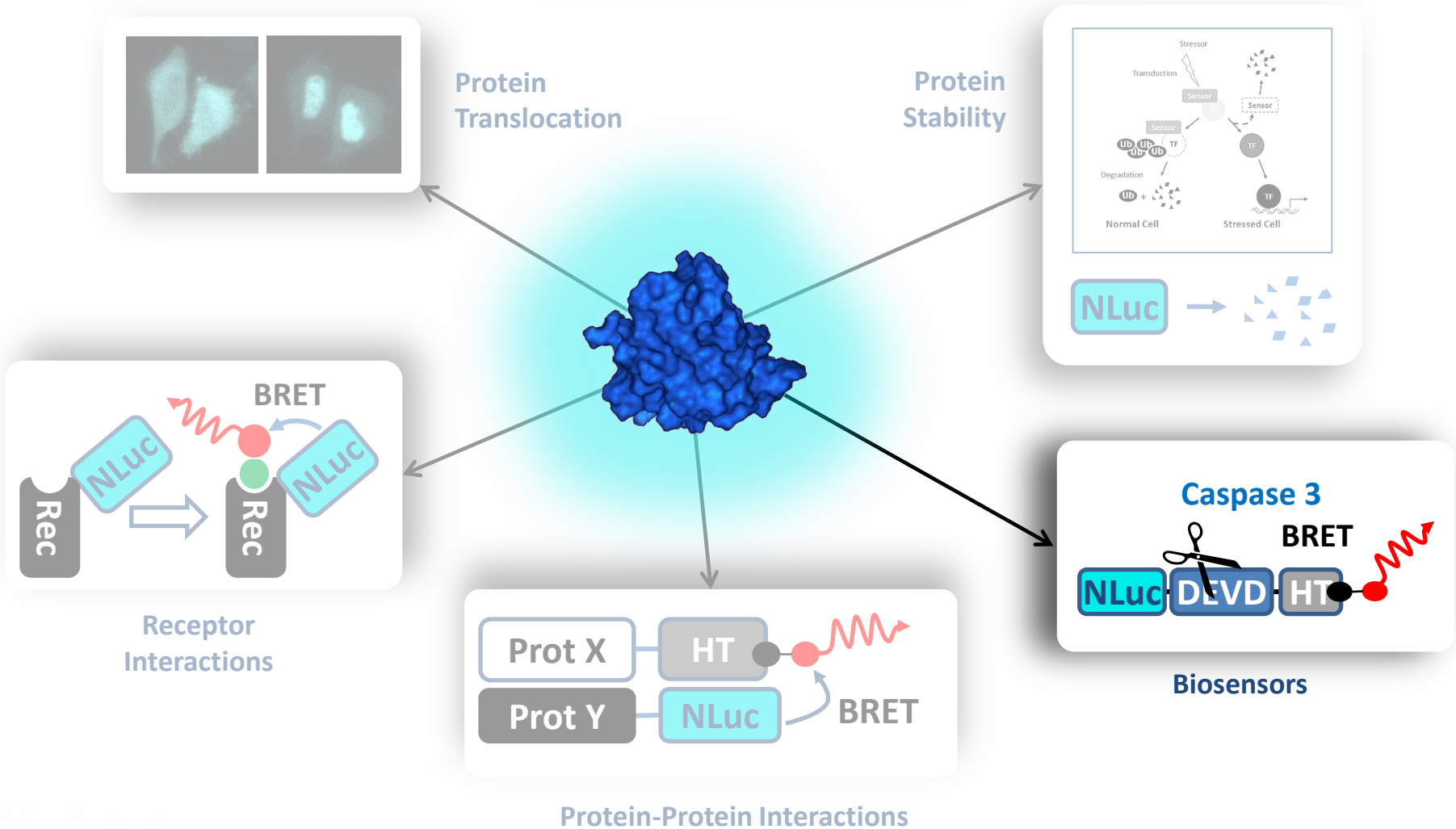


# Multiple uses with one genetic construct

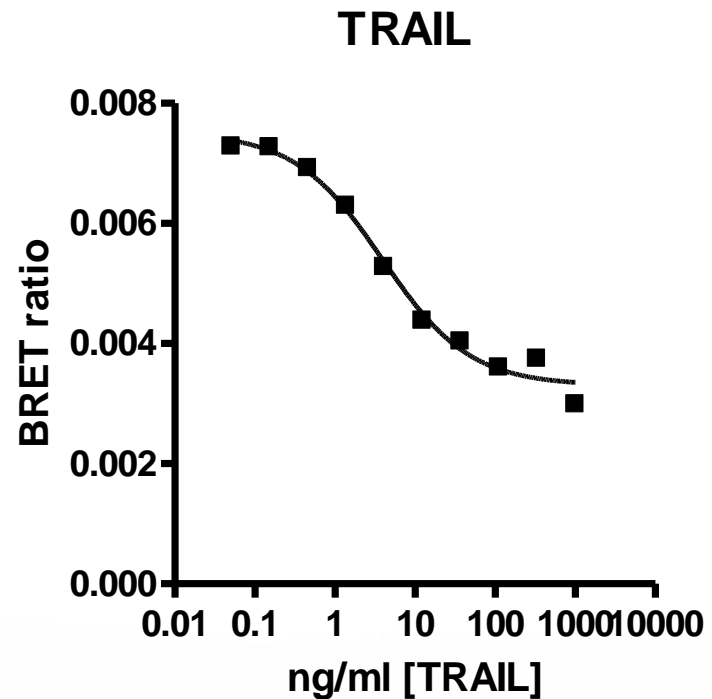
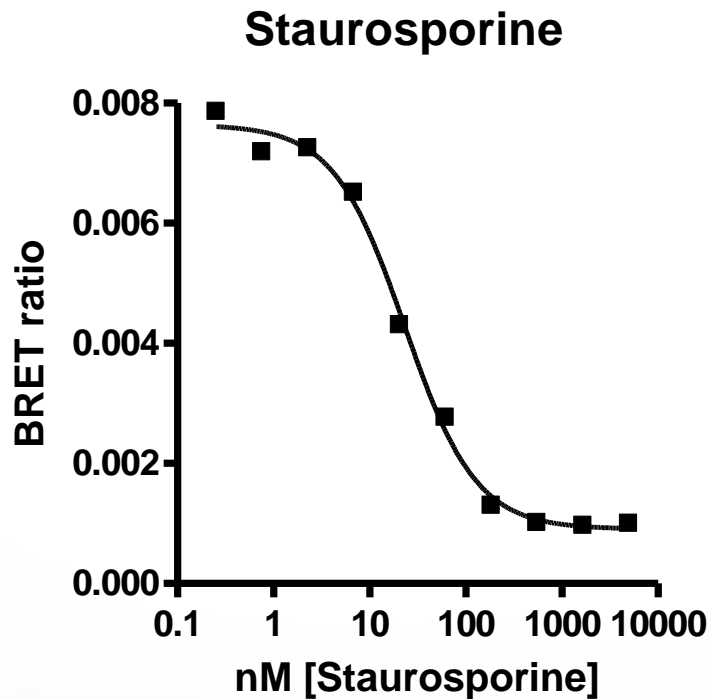
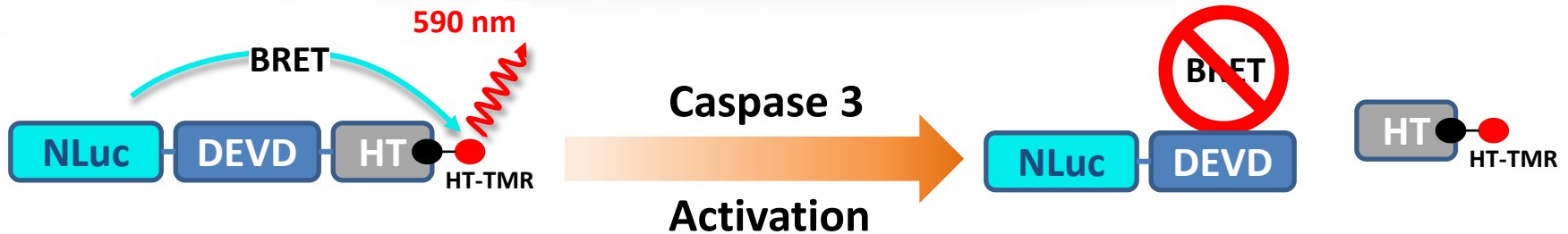




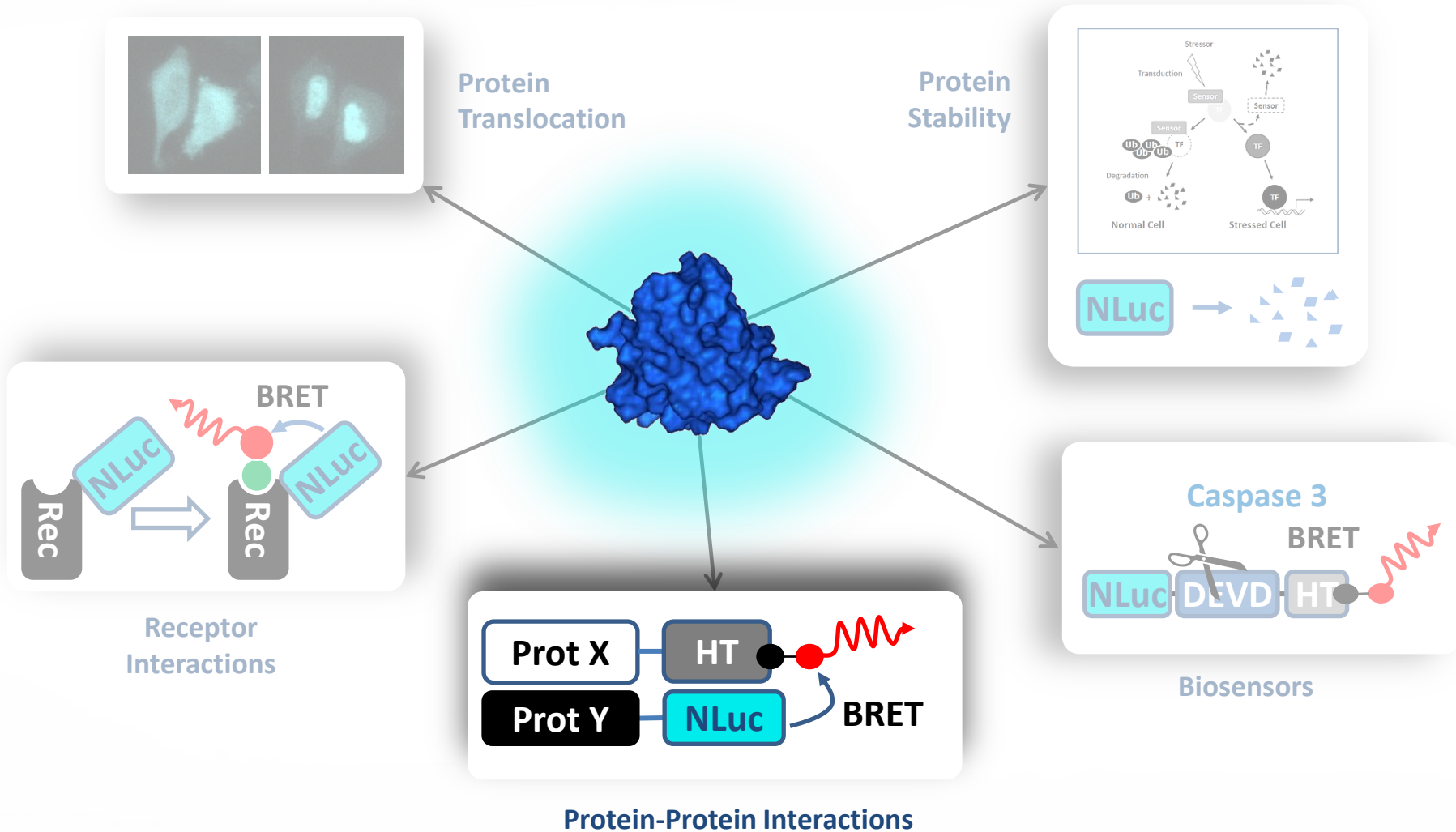
# NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments



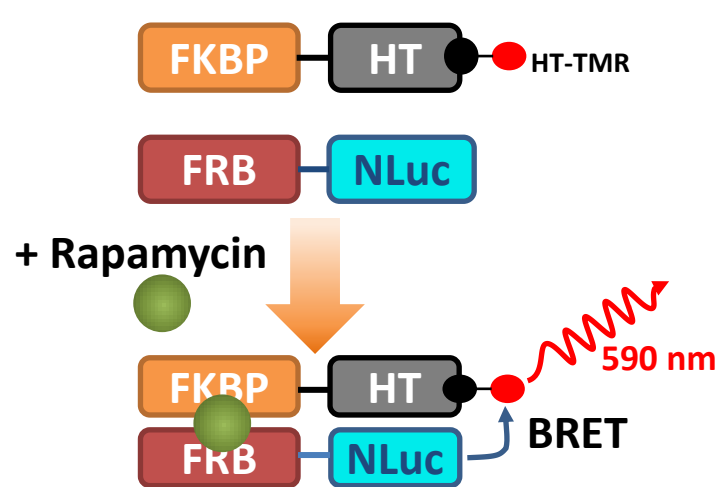
# Biosensor: Disrupting BRET with a cleavable bridging peptide sequence



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



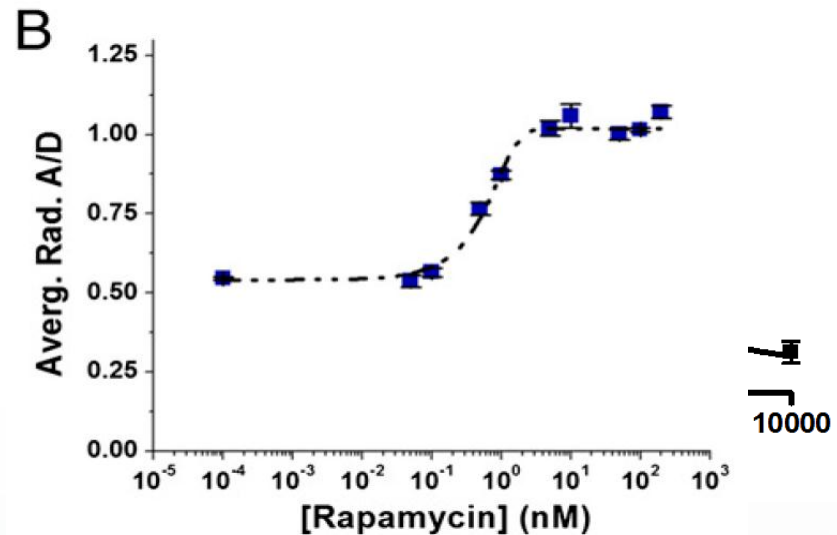
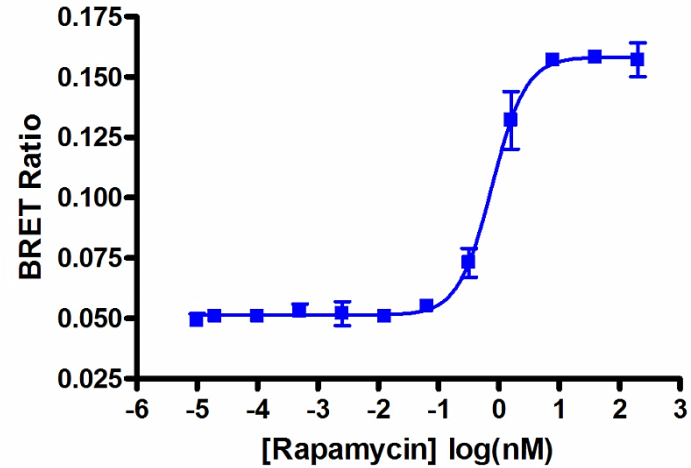
# Can NLuc:HT Pair be used for Protein-Protein BRET?



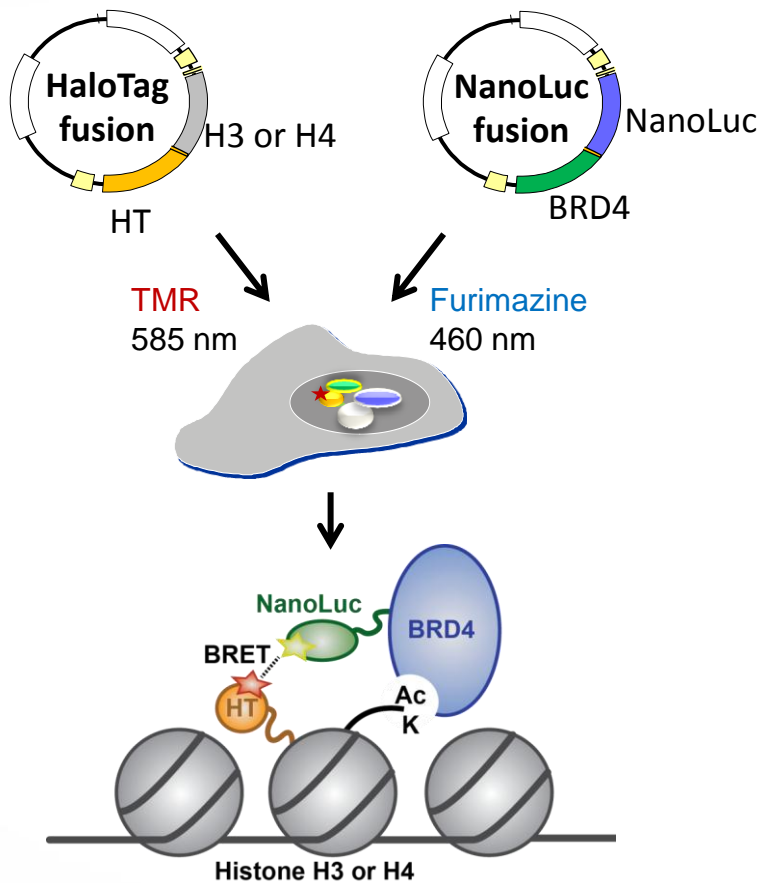
Same model system used with BRET 6 System

RLuc8.6 → TurboFP

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

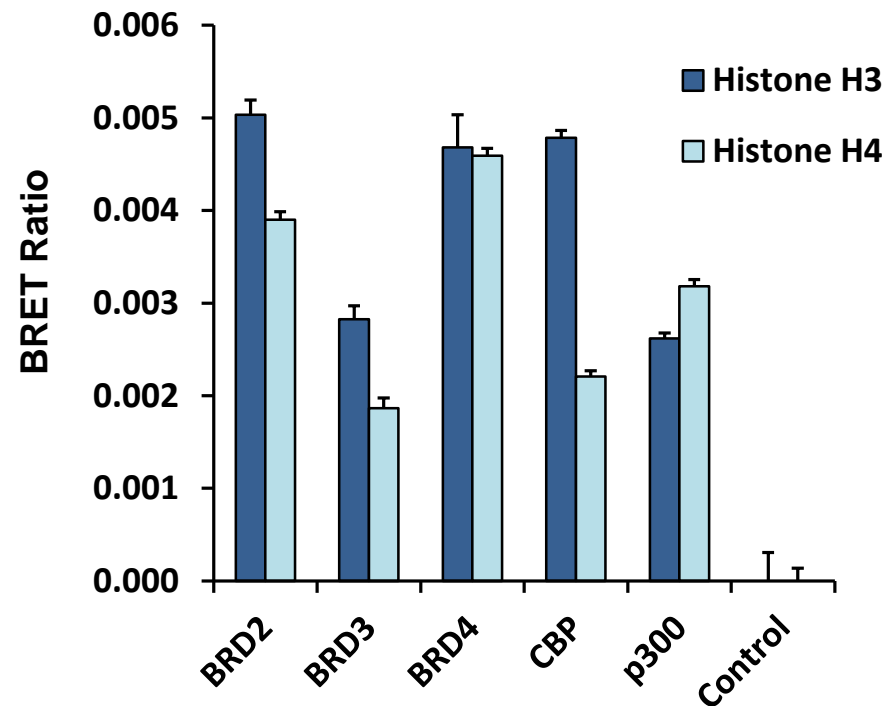


# BRET assay for detecting bromodomain interactions with histones



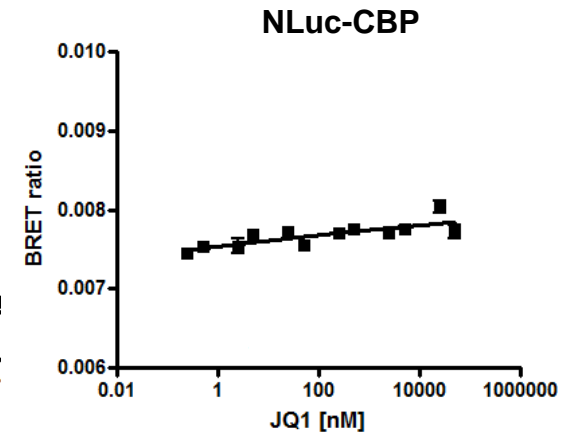
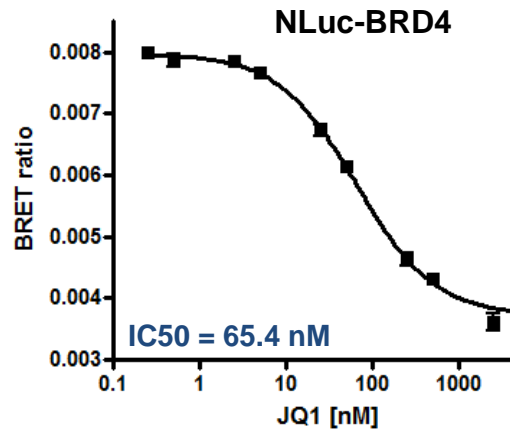
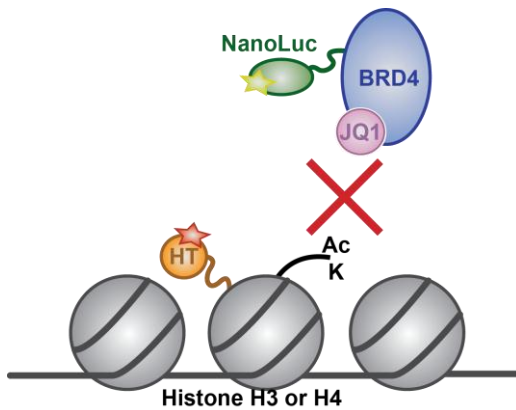
BRET (600/480nm) Measurement

## BRET Measurement in HCT116 Cells

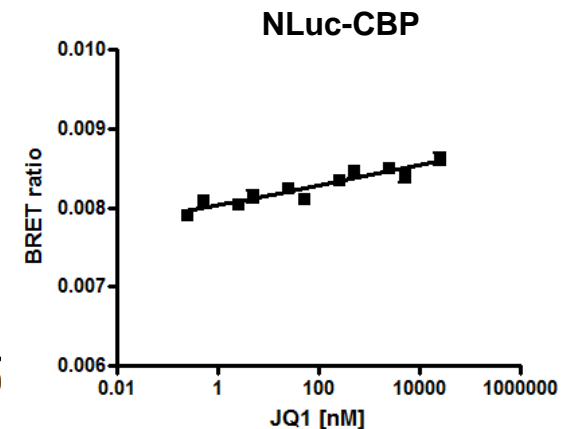
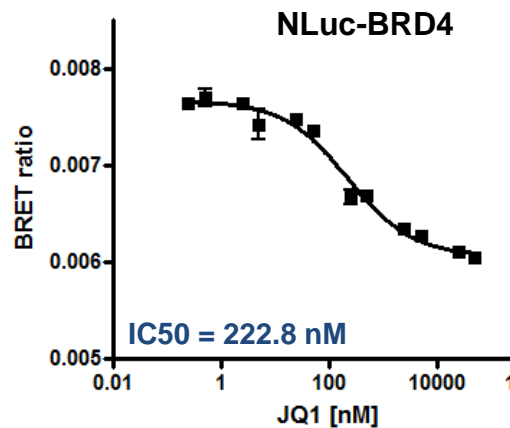


# Measuring effect of BET inhibitor on histone-bromodomain interactions

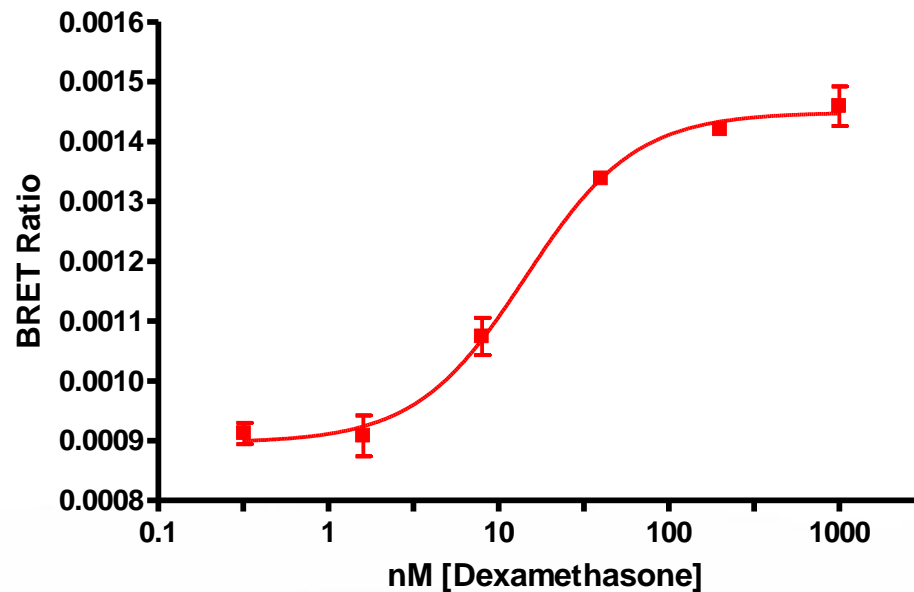
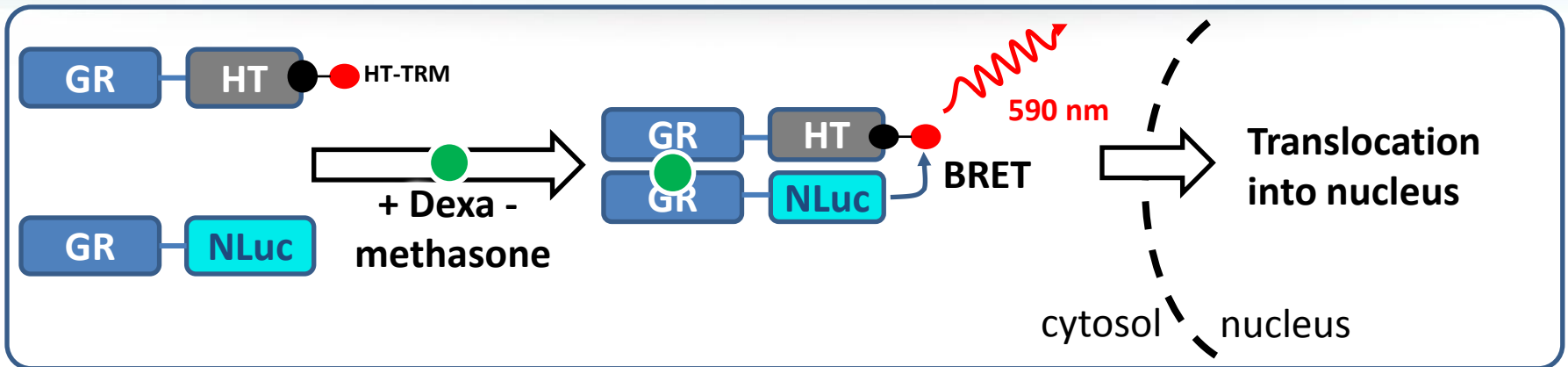
## Histone H3.3-HaloTag



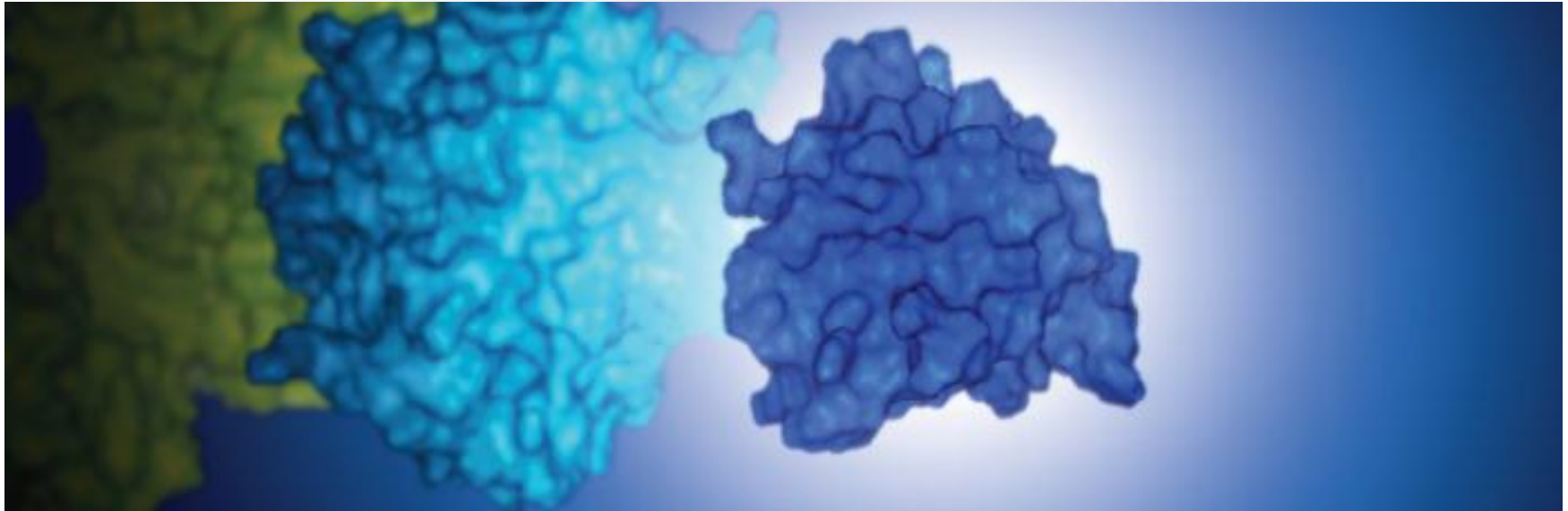
## Histone H4-HaloTag



# Glucocorticoid Receptor Dimerization



# *Bright Future for NanoLuc™ Luciferase fusions*



Brightness improves  
bioluminescent imaging

Versatility to go anywhere in cell

Versatility to allow stability  
measurements

Brightness allows BRET with  
HaloTag® Fusions

- Biosensors
- Protein:Protein Interactions

Brightness allows BRET with  
fluorescent ligands

- Ligand binding assays



# Nano-Glo™ Luciferase Assay System



Just make enough 1X Nano-Glo Assay Reagent from the 50X substrate and buffer to meet your needs. No need to make up all the reagent at once.

## **Cat # N1110: 10ml (100 assays)**

- 200µl Nano-Glo™ Luciferase Assay Substrate
- 10ml Nano-Glo™ Luciferase Assay Buffer

## **Cat # N1130 10x10ml (1,000 assays)**

- 10 x 200µl Nano-Glo™ Luciferase Assay Substrate
- 10 x 10ml Nano-Glo™ Luciferase Assay Buffer

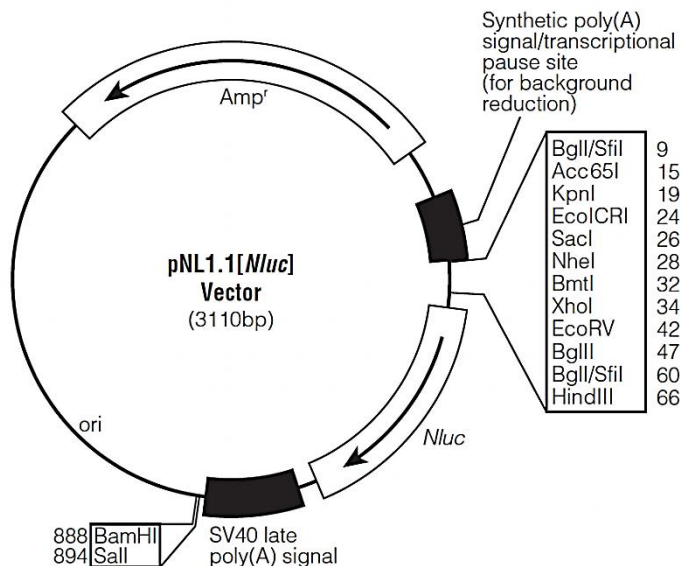
## **Cat.# N1120 100ml (1,000 assays)**

- 2 x 1ml Nano-Glo™ Luciferase Assay Substrate
- 100ml Nano-Glo™ Luciferase Assay Buffer

## **Cat.# N1150 10x100ml (10,000 assays)**

- 5 x 4ml Nano-Glo™ Luciferase Assay Substrate
- 10 x 100ml Nano-Glo™ Luciferase Assay Buffer

# Currently Available Vectors



The pNL vectors are built on the pGL4 Vector backbone allowing transfer of inserts from pGL4 series vectors to the pNL vectors.

Plasmid	Reporter	Marker	Promoter	MCS	Catalog #
pNL1.1	Nluc	-	-	Y	N1001
pNL1.2	NlucP	-	-	Y	N1011
pNL1.3	secNluc	-	-	Y	N1021
pNL2.1	Nluc	Hygro	-	Y	N1061
pNL2.2	NlucP	Hygro	-	Y	N1071
pNL2.3	secNluc	Hygro	-	Y	N1081
pNL3.1	Nluc	Hygro	minP	Y	N1031
pNL3.2	NlucP	-	minP	Y	N1041
pNL3.3	secNluc	-	minP	Y	N1051
pNL1.1.CMV	Nluc	-	CMV	N	N1091
pNL1.3.CMV	secNluc	-	CMV	N	N1101
pNL3.2.NF-κB-RE	NlucP	Hygro	NF-κB-RE/minP	N	N1111

10321MA

## ***Want to know what vectors are coming?***

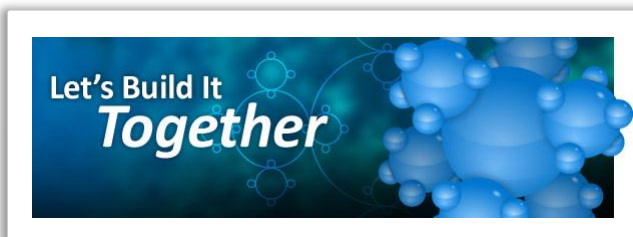
New vectors using Promega exclusive technologies are typically offered as a custom product before they become a catalog product.



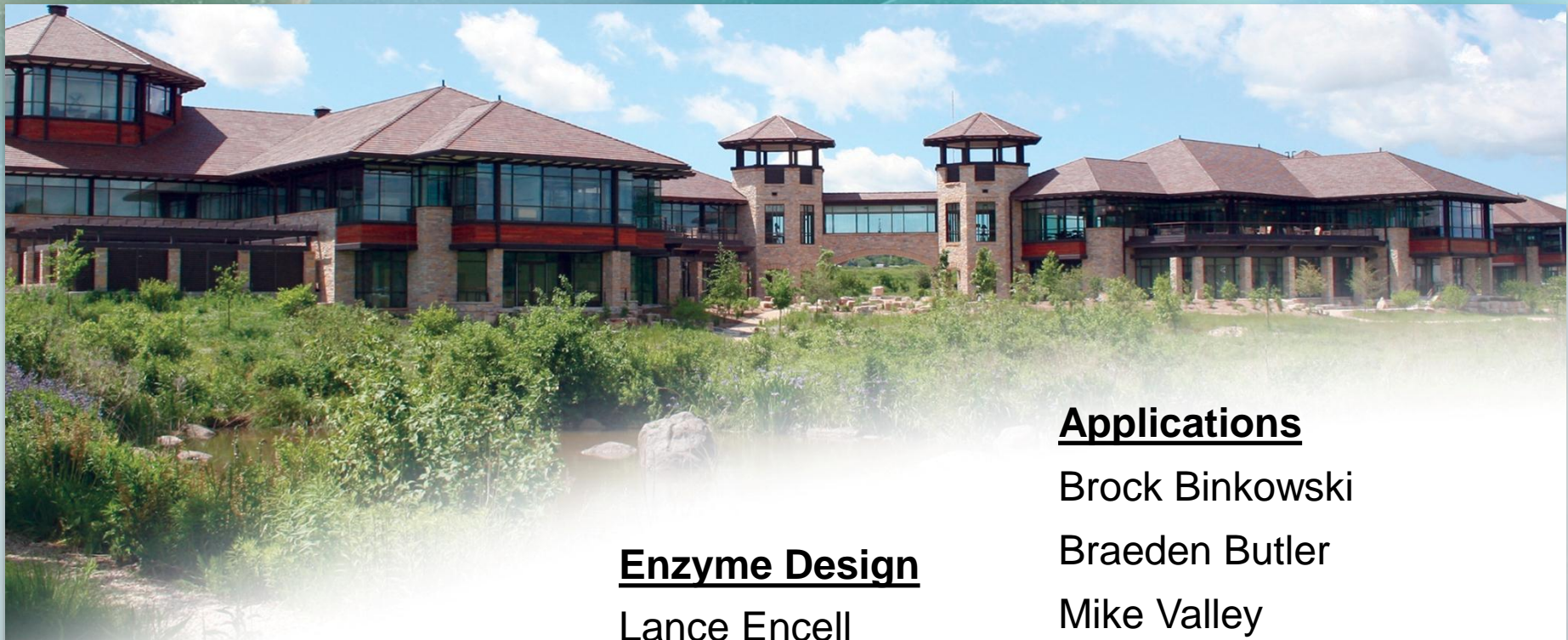
Visit the Current Research Materials page at:

**[www.promega.com/cam](http://www.promega.com/cam)**

Stay on the cutting edge and get research tools before they are available to everyone!



You can also request construction of a special vector to meet your exact needs with ***Custom Assay Services***



**Keith Wood**

**Substrate Design**

Poncho Meisenheimer

James Unch

Ruslan Arbit

Hui Wang

Dieter Klaubert

**Enzyme Design**

Lance Encell

Monika Wood

Mary Hall

Kris Zimmerman

Paul Otto

Hélène Benink

Gedi Vidugiris

Mike Slater

**Applications**

Brock Binkowski

Braeden Butler

Mike Valley

Matt Robers

Thomas Machleidt

Chris Eggers

Frank Fan

Danette Daniels

Marie Schwinn

Jacqui Mendez

***Questions Welcome***