

Applications of a smaller, brighter, more versatile luciferase:

NanoLuc™ Luciferase Technology

Kyle Hooper, Ph.D.

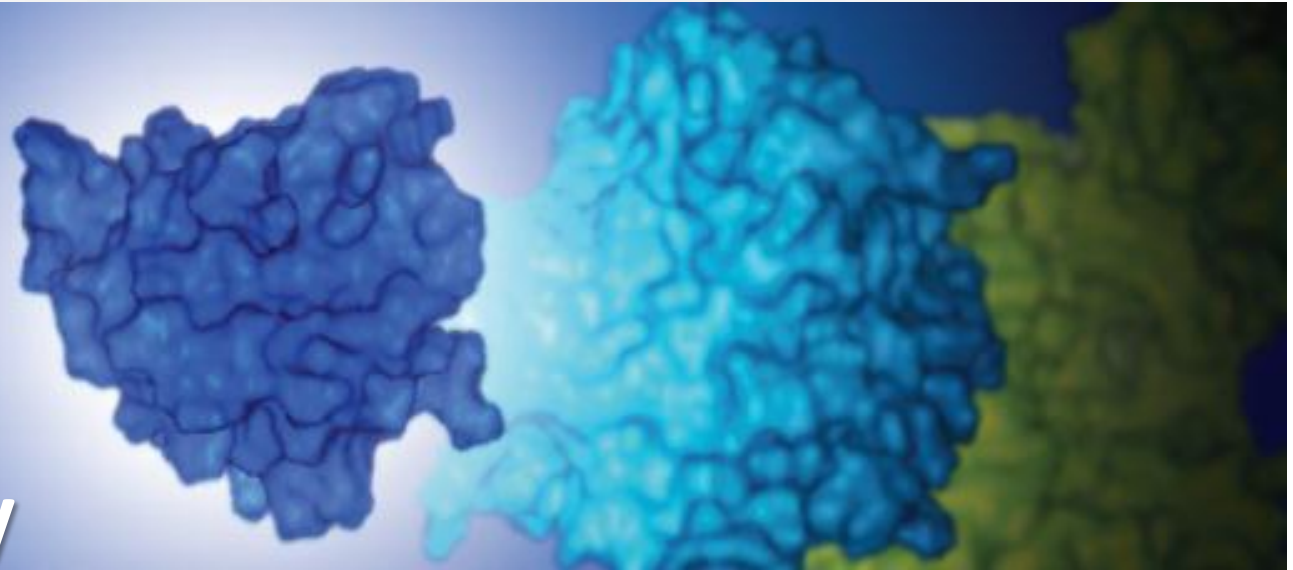
Fall 2012



Today, I will speak about...



NanoLuc™ Luciferase Technology



Origins of NanoLuc Luciferase

NanoLuc Luciferase as a reporter

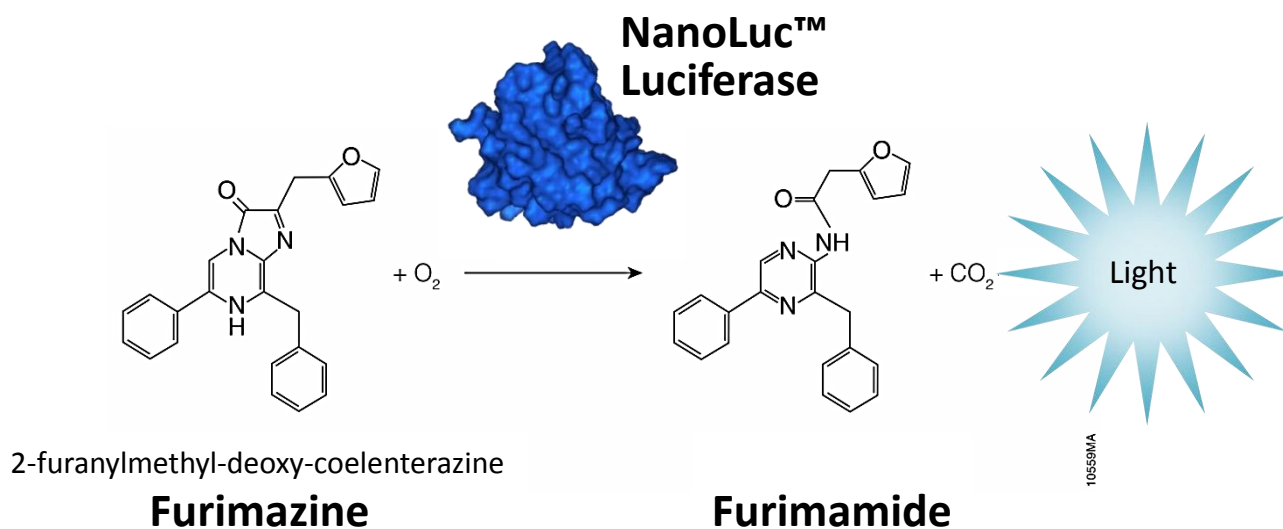
- Full-Length **Nluc**
- Destabilized, full-length **NlucP**
- Secreted, full-length **secNluc**

NanoLuc Luciferase as a fusion partner

- Protein Translocation
- Protein Stability
- Protein:Protein Interactions
- Receptor Interactions
- Biosensors

What is NanoLuc™ Luciferase?

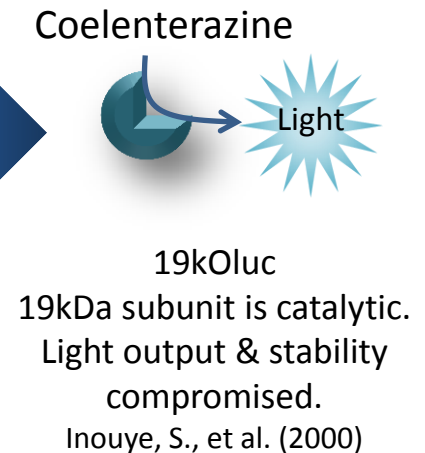
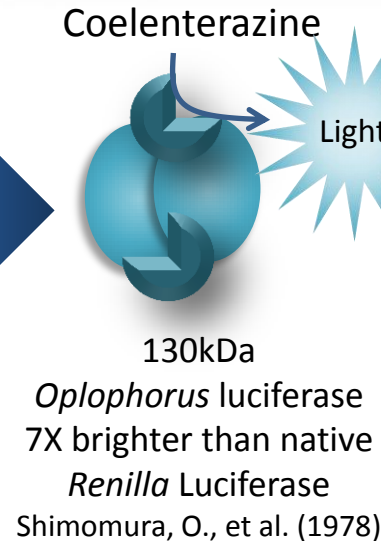
NanoLuc™ (Nluc) is a 19.1 kDa, ATP-independent luciferase that utilizes a novel coelenterazine analog (furimazine) to produce high intensity, glow-type luminescence.



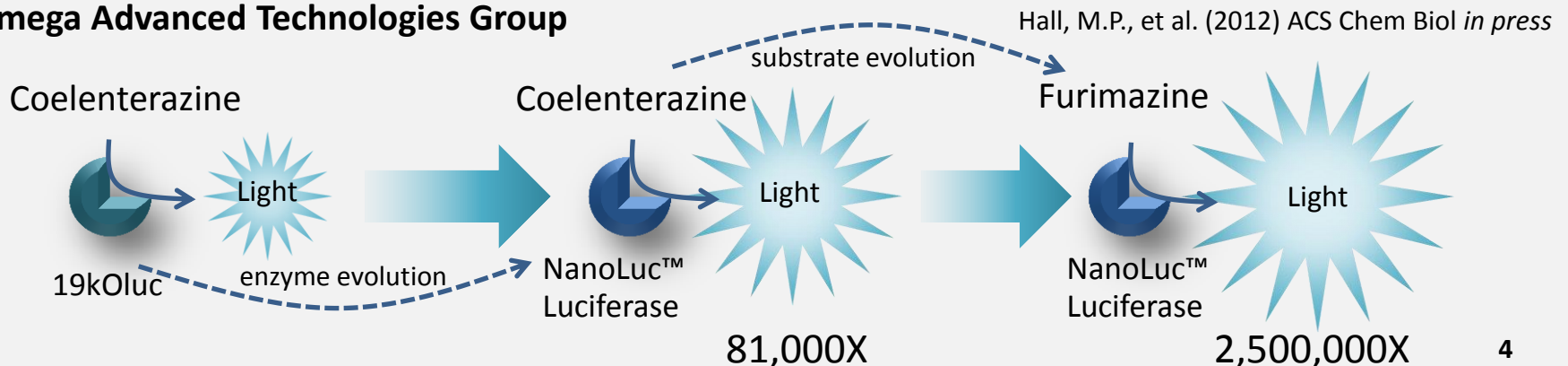
Evolution of NanoLuc from ocean to lab bench



Oplophorus gracilirostris first cataloged in 1881



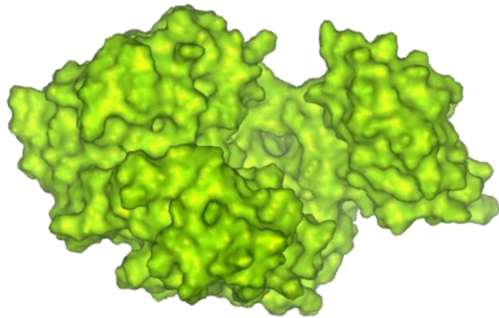
Promega Advanced Technologies Group



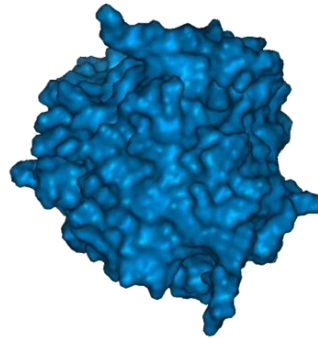
NanoLuc™ is very small



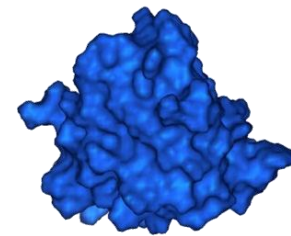
Firefly (Fluc)



Renilla (Rluc)

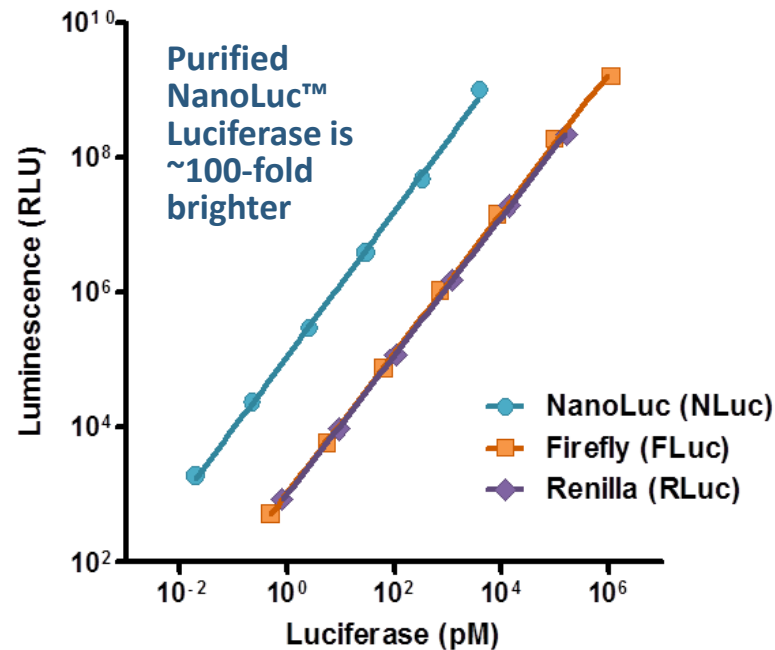


NanoLuc™ (Nluc)

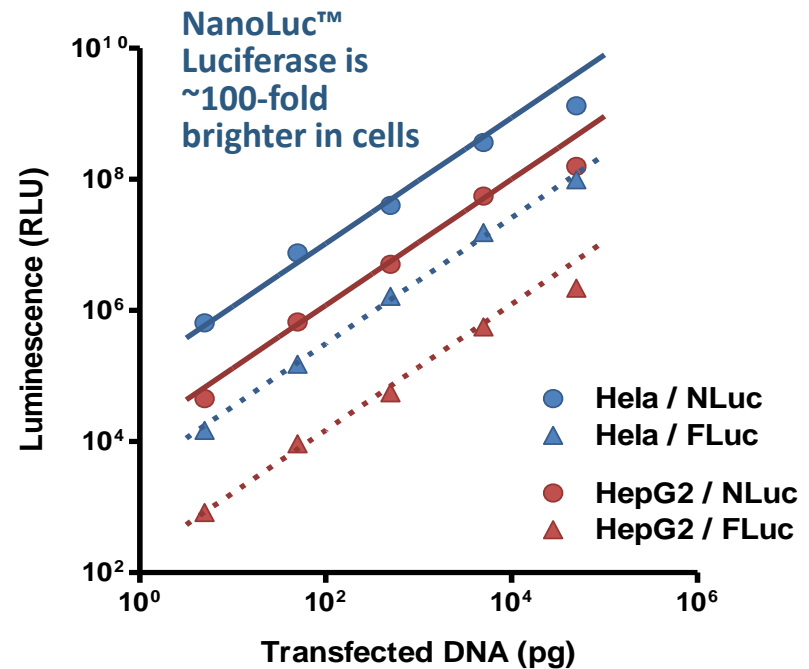


	Amino acids	M.W.	Mol. Vol. Å ³
Nluc	171	19.1	14
Rluc	312	36.0	32
Fluc	550	60.6	44

NanoLuc is bright



Recombinant NLuc/Nano-GloTM Assay
Recombinant FLuc/ONE-GloTM Assay
Recombinant RLuc/Renilla-GloTM Assay



CMV-driven NLuc/Nano-GloTM Assay
CMV-driven FLuc/ONE-GloTM Assay

NanoLucTM 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

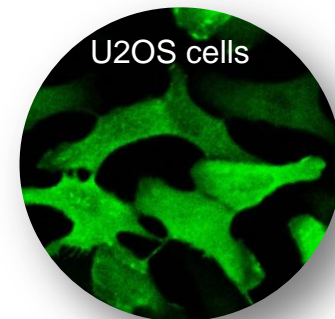
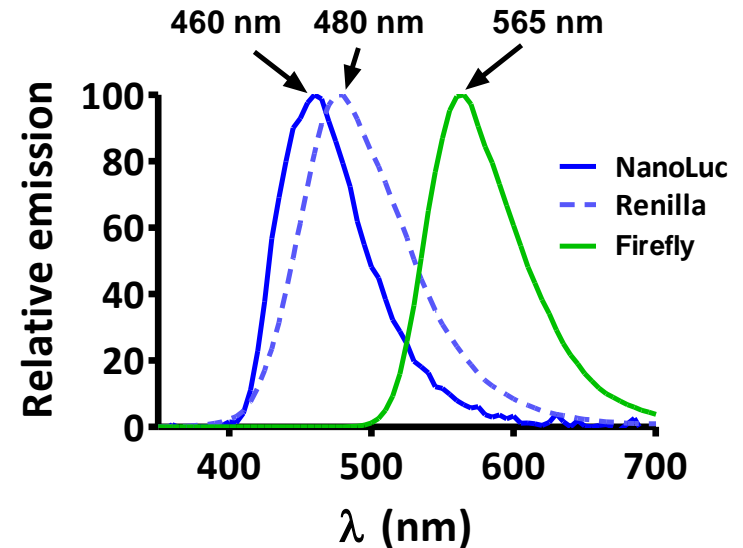
No post-translational modifications detected in mammalian cells

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)

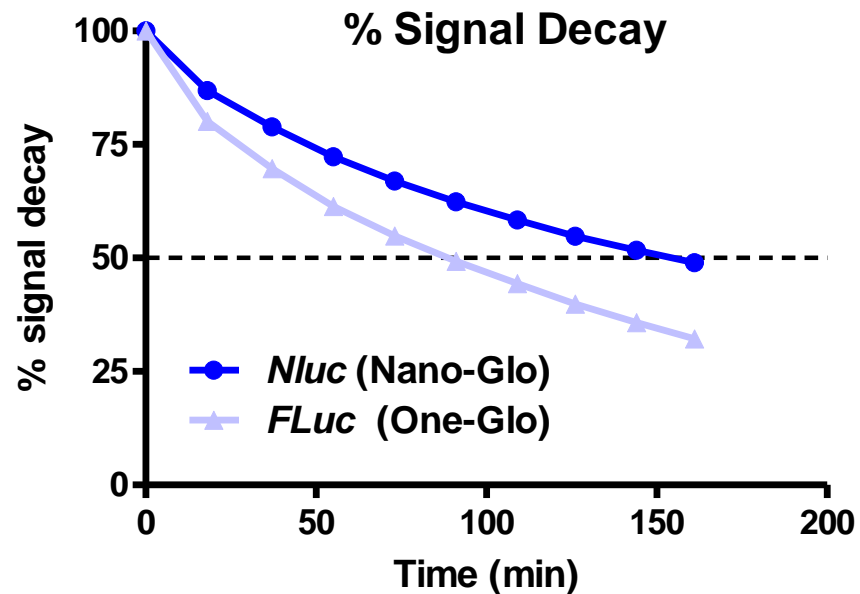
- Half-life routinely >2 hour at room temperature

Low autoluminescence background

- Enhances assay sensitivity

Stable reconstituted reagent:

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



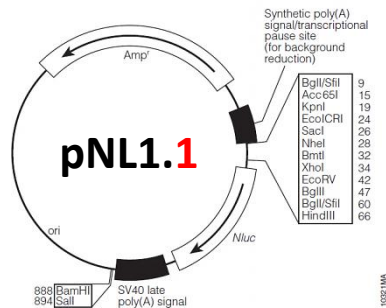
Add-Mix-Measure format like

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

3 Varieties of NanoLuc™ Luciferase for you



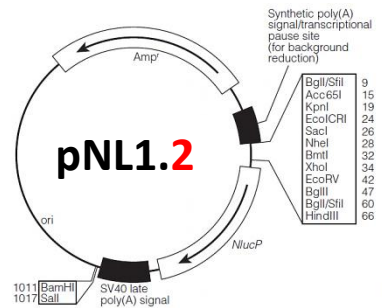
Intracellular Formats



NanoLuc™ Luciferase

Nluc (513 bp)

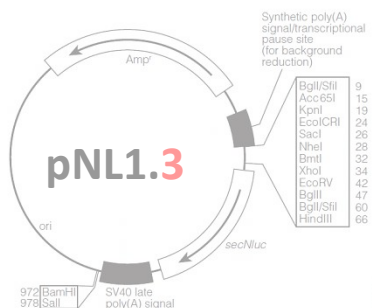
Protein destabilization domain



NanoLuc™ Luciferase **PEST**

NlucP (636 bp)

Secretion Format



IL6 **NanoLuc™ Luciferase**

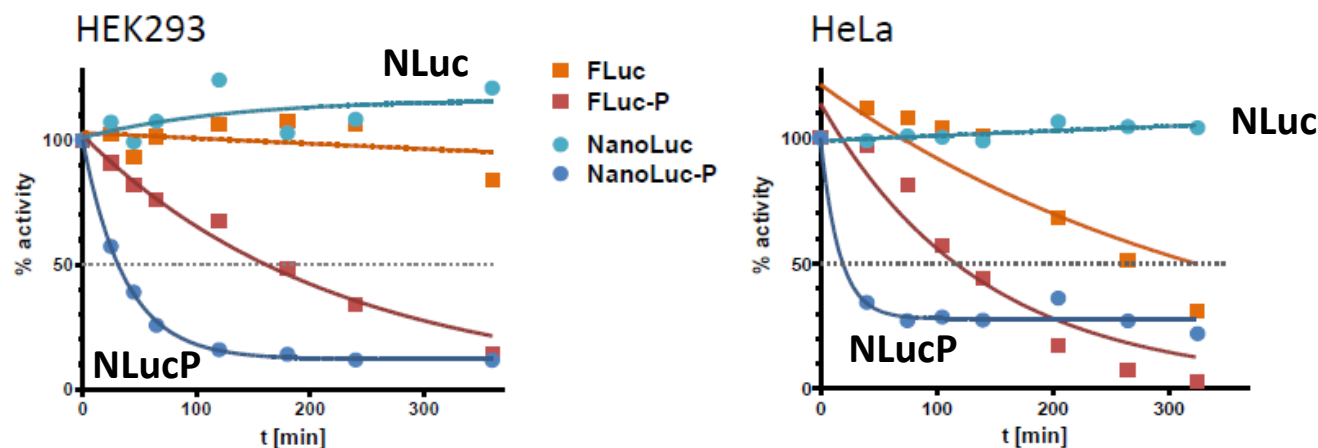
secNluc (597 bp)

Secretion signal

Intracellular stability of NanoLuc™ & Firefly



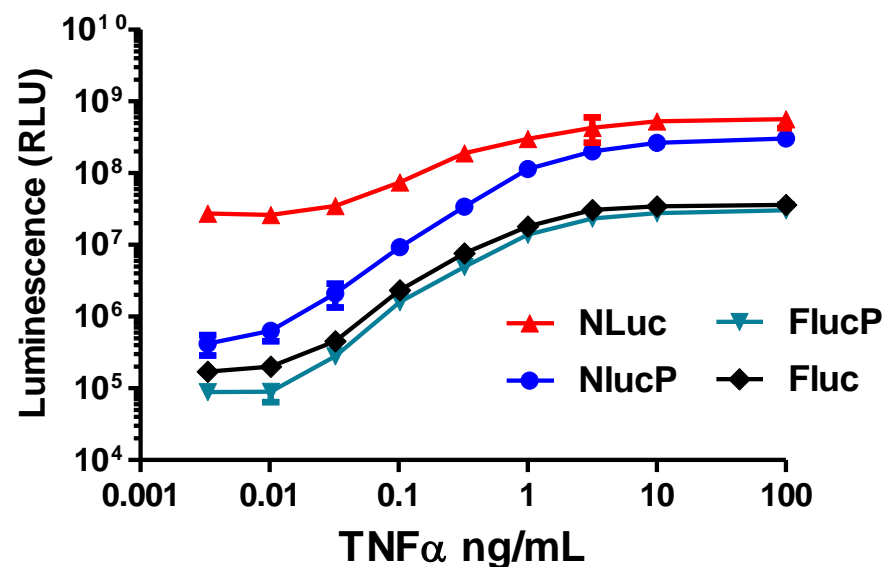
New protein synthesis blocked by addition of cycloheximide



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-κB inducible constructs. rhTNFα treatment for 5 hours.

Brightness

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

(18 experiments)

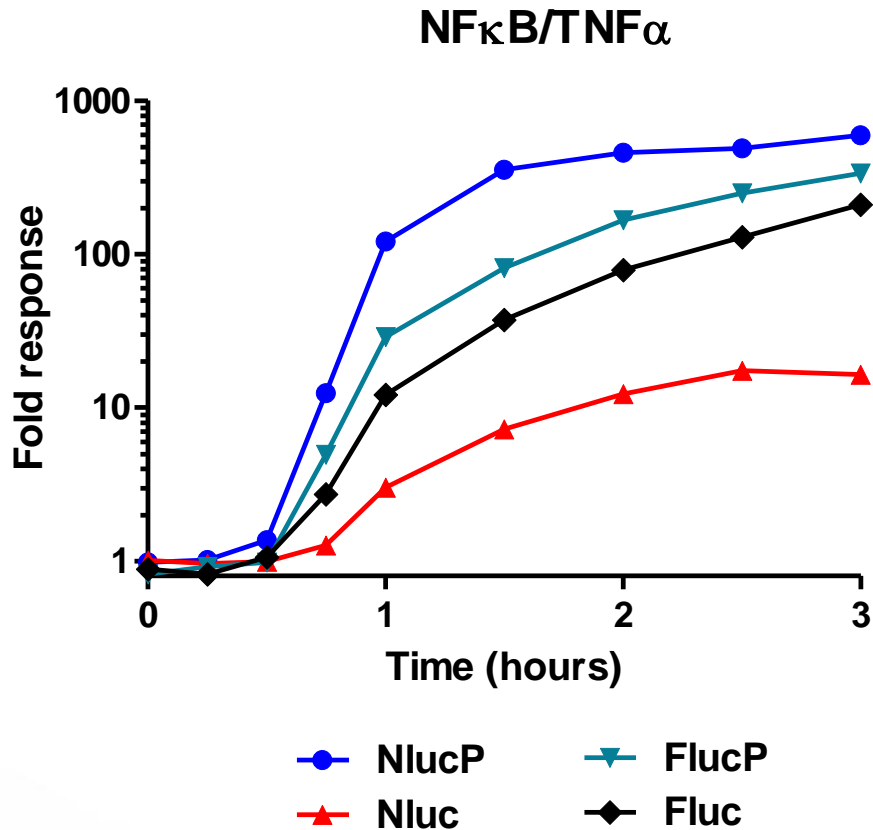
NLuc 13-236 fold brighter than Fluc (79 fold avg.)

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

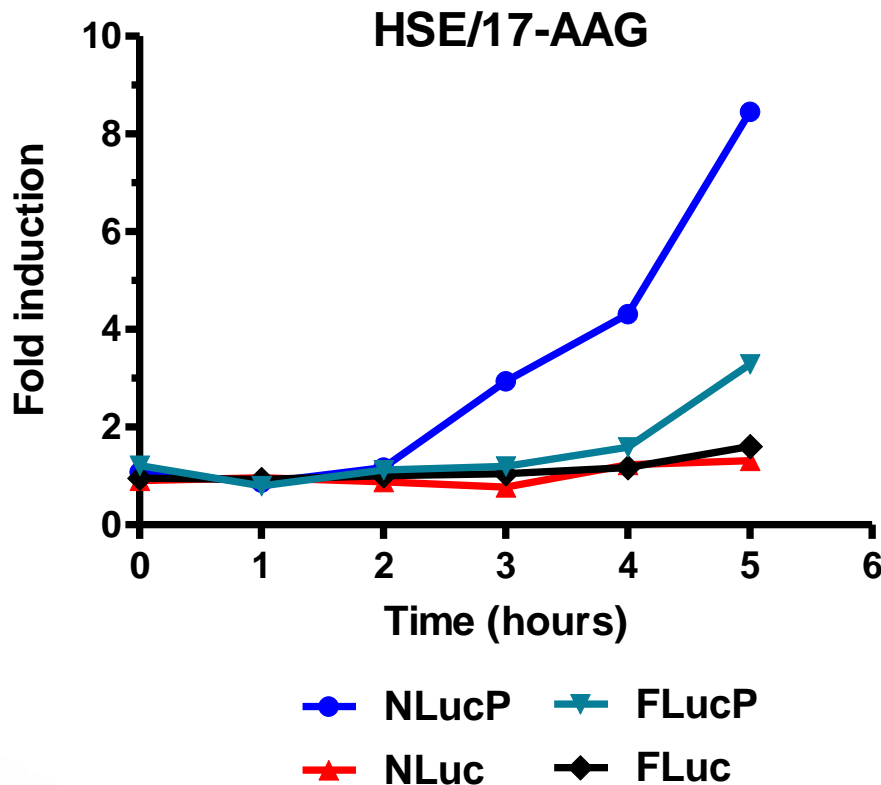


Relative Response

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

Experimental details: transient transfection of HEK293 cells with NF κ B inducible constructs; addition of 100 ng/ml rhTNF α at time zero.

NLucP allows study of weakly induced responses



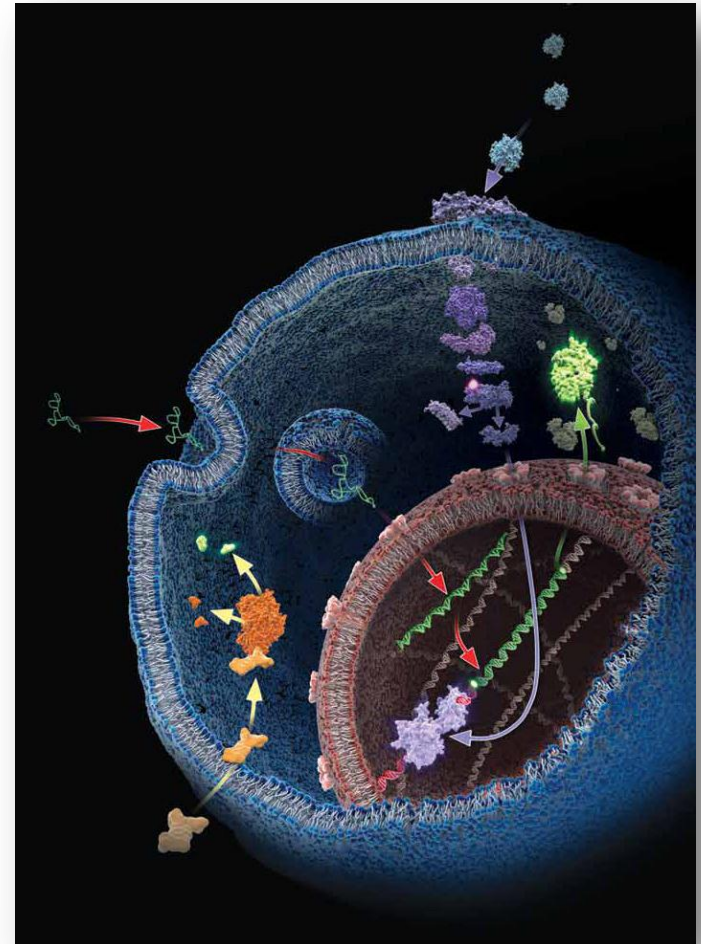
Relative Response

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

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

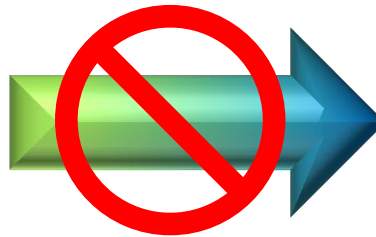
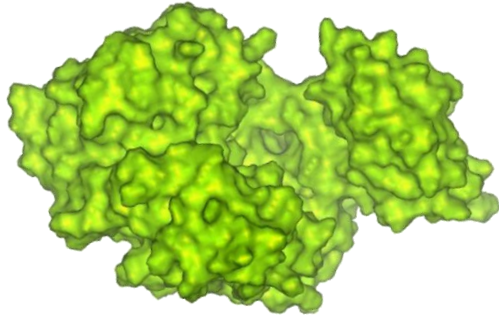
NanoLuc Luciferase as an intracellular reporter

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



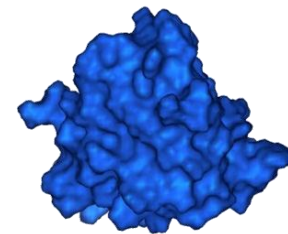
Should I switch from Firefly to NanoLuc™ Luciferase?

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

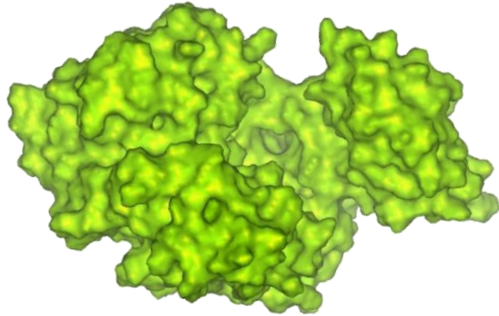
- ✓ Excellent signal:background
- ✓ Excellent dynamic range

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

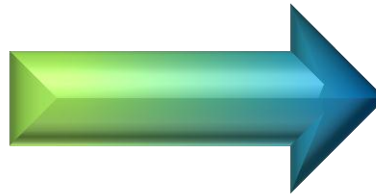
ARE	HSE	ISRE	STAT5	SRE
p53	HRE	SIE	NFAT	SRF
ATF6	XRE	SBE	CRE	
MRE	AP1	TCF-LEF	NF-κB	

Should I switch from Firefly to NanoLuc™ Luciferase?

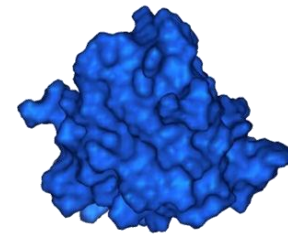
Firefly (Fluc)



Yes, if ...



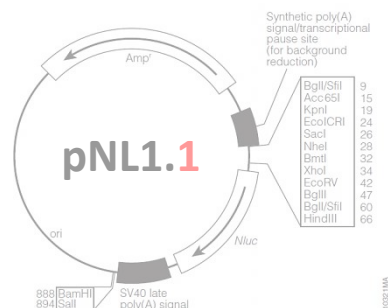
NanoLuc™ (Nluc)



- ✓ **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

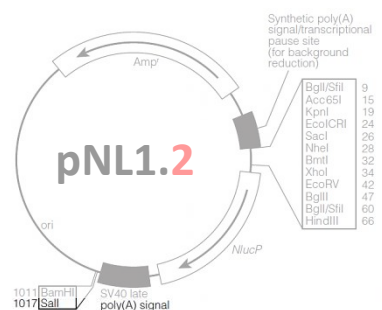
Intracellular Formats



NanoLuc™ Luciferase

Nluc (513 bp)

Protein destabilization domain

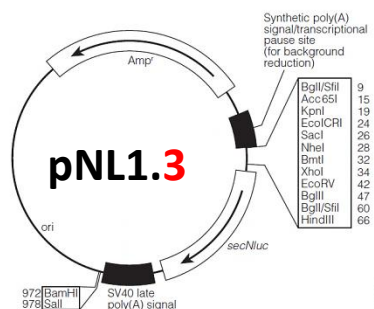


NanoLuc™ Luciferase

NlucP (636 bp)

PEST

Secretion Format



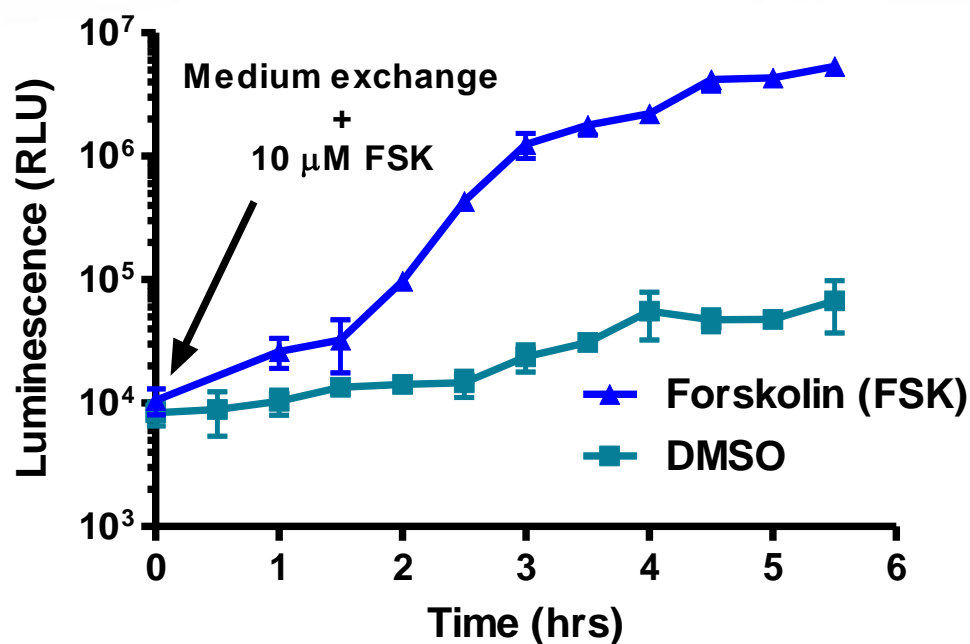
IL6

NanoLuc™ Luciferase

secNluc (597 bp)

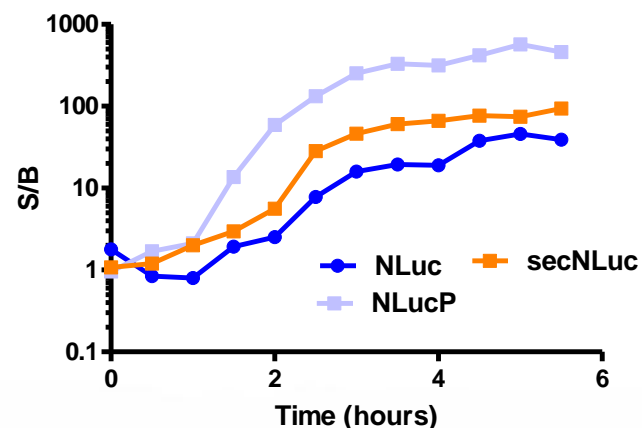
Secretion signal

Secretion based format using secNluc

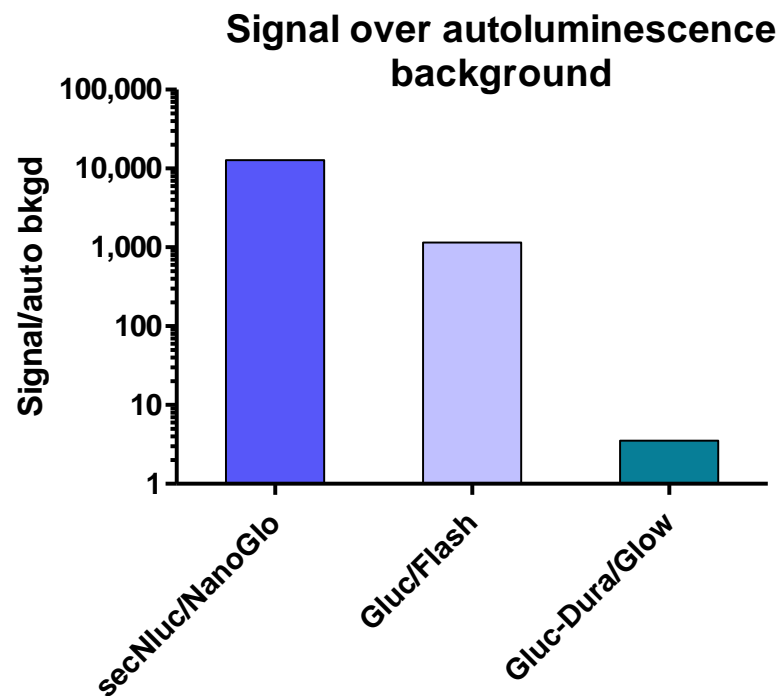
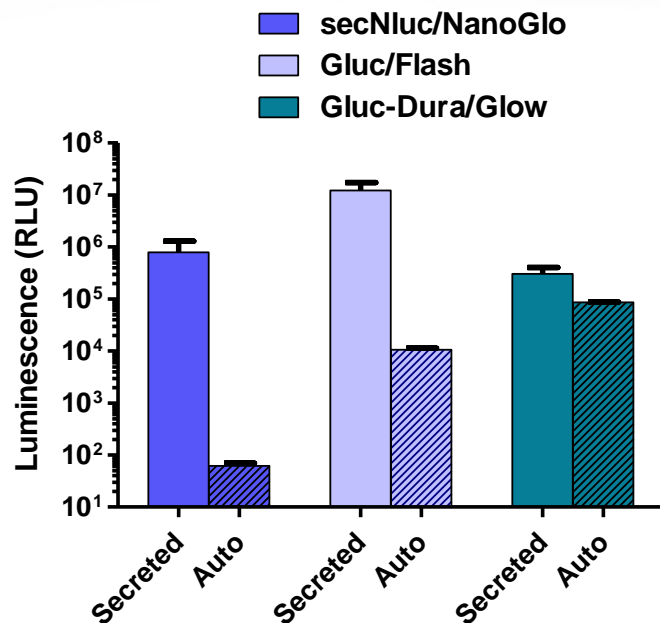


Experimental details: transient transfection of HEK293 cells with CREB inducible construct; addition of 10 μ M forskolin at time zero.

- 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



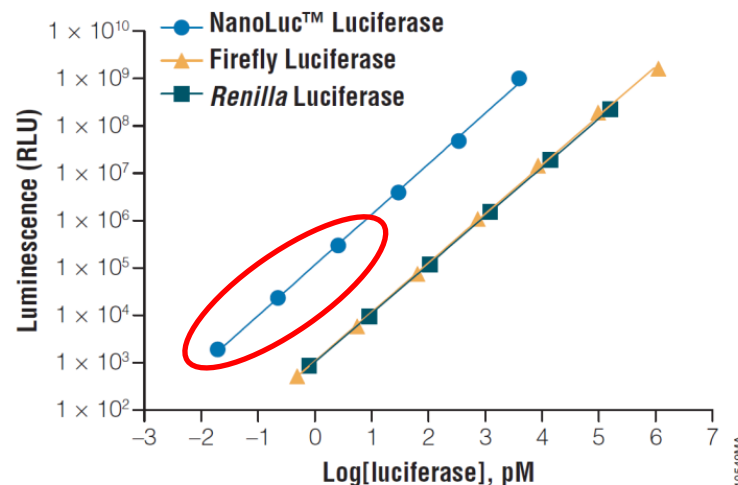
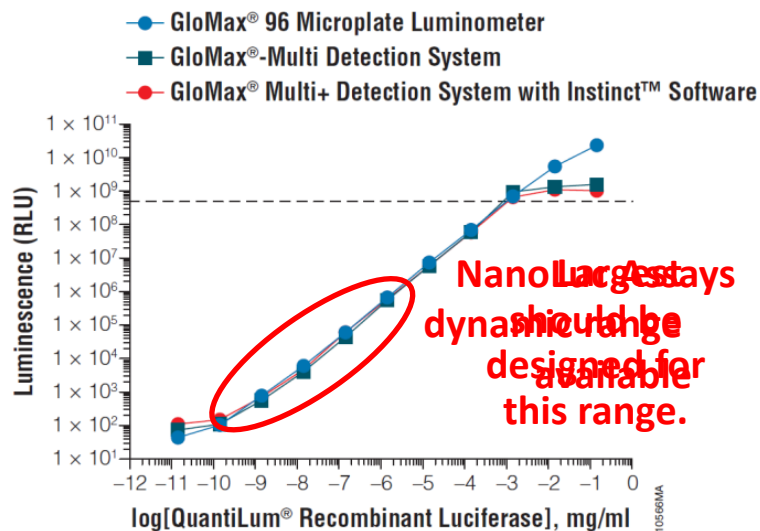
**Gluc kits: high background limits sensitivity
& dynamic range**

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

Handling the bright signal



R&D used GloMax[®] Instruments for NanoLuc[™] Development



GloMax[®]-Multi+
Microplate Multimode Reader



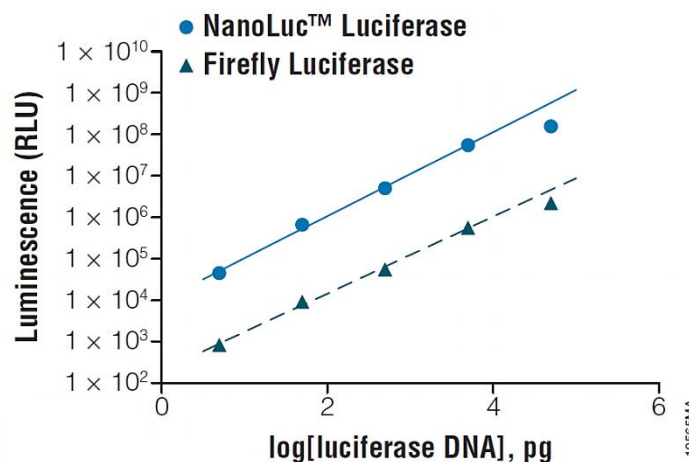
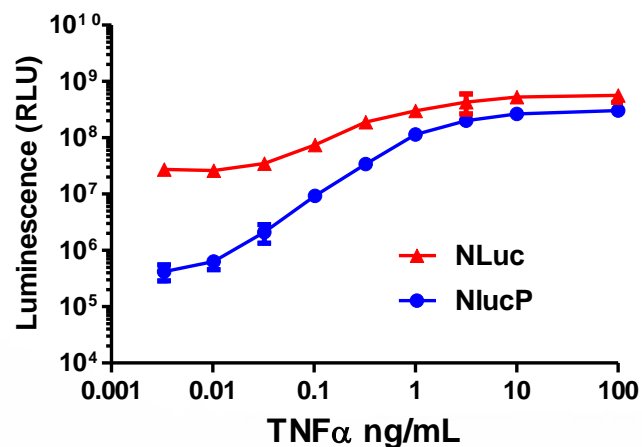
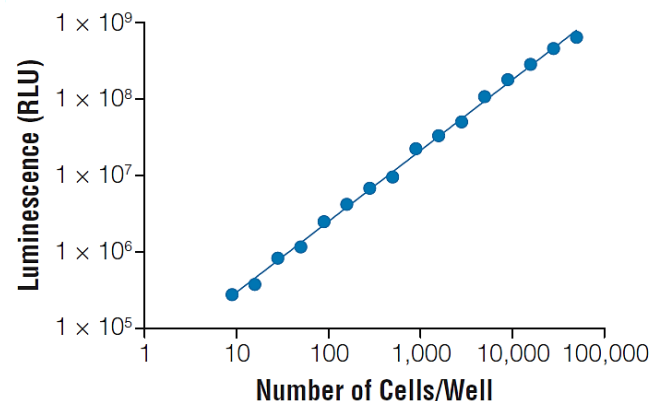
GloMax[®]-Multi
Microplate Multimode Reader



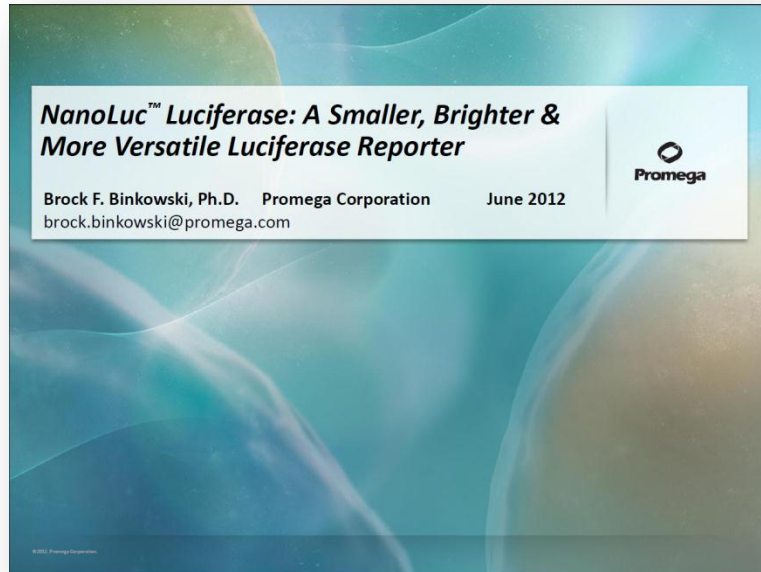
GloMax[®] 96
Microplate Luminometer

Getting NanoLuc™ Luciferase Signals in range

- ✓ Transfect fewer cells
- ✓ Transfect less DNA
- ✓ Switch to NlucP
- ✓ Use a weaker constitutive promoter



More details on NanoLuc™ Development



Webinar recording by lead R&D scientist for development of the NanoLuc™ Vectors and Nano-Glo™ Assay System

**Brock F. Binkowski, Ph.D.
Sr. Research Scientist II**

Broadcast from June 2012
~30 minutes long

www.promega.com/webinars

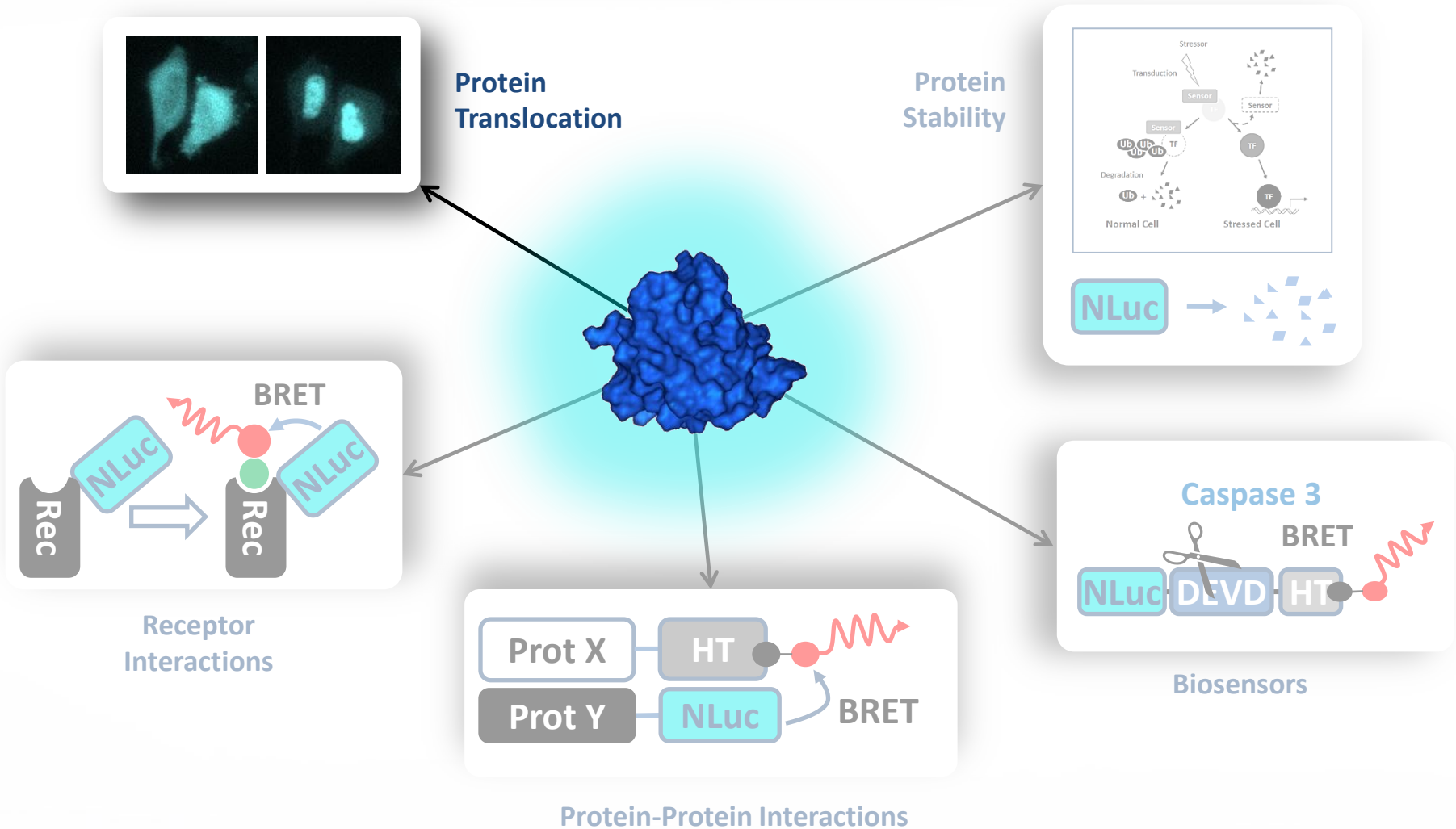
Click *Previous Webinars* in the grey box

NanoLuc™ Luciferase as a protein function probe

Applications of full-length NanoLuc™ Luciferase.



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

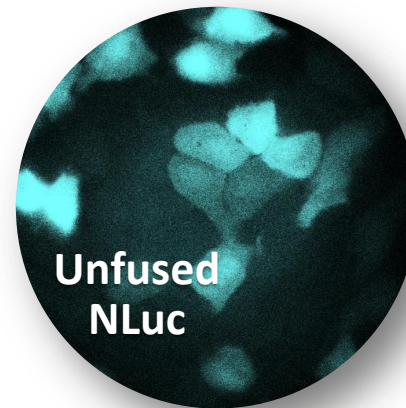


NanoLuc™ Luciferase excels in bioluminescent imaging applications



Nluc brightness leads to short exposure times:

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



*NanoLuc &
LV200 featured*

@ASCB 2012

*Olympus
Product Showcase*

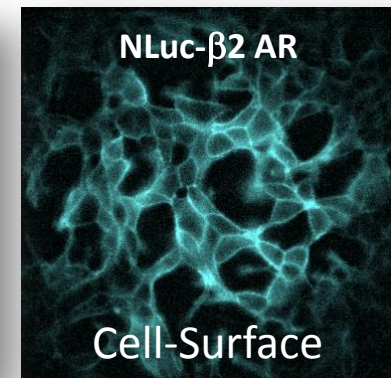
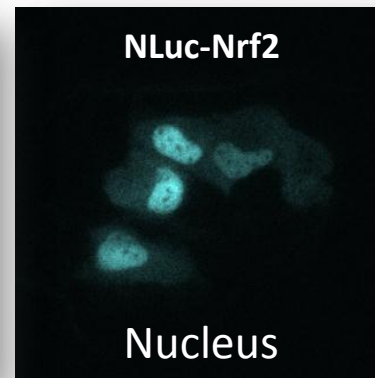
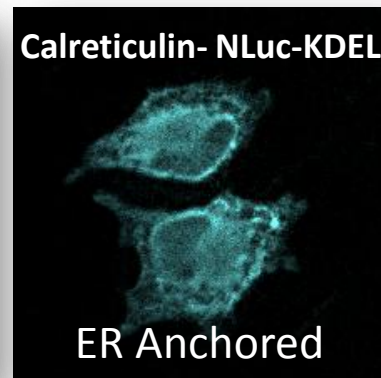
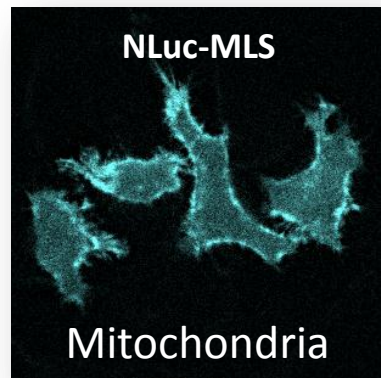
Why bother? Fluorescence works.

- 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™ Fusions can go anywhere...

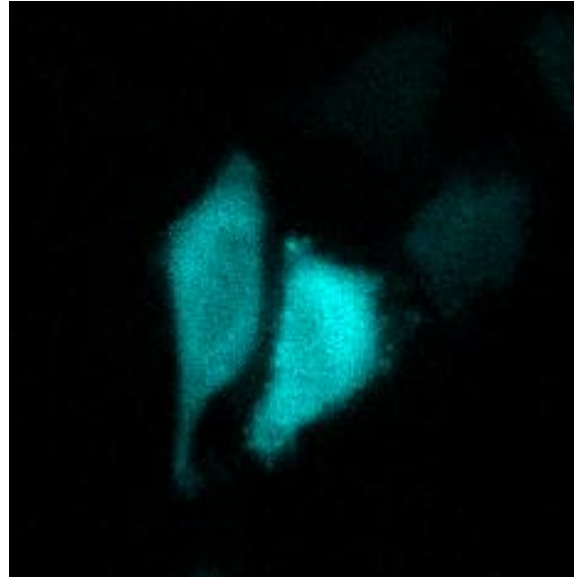


NanoLuc™ Luciferase fusions could be a useful tool to investigate cell biology

Bioluminescence imaging of protein translocation



cytoplasm → nucleus

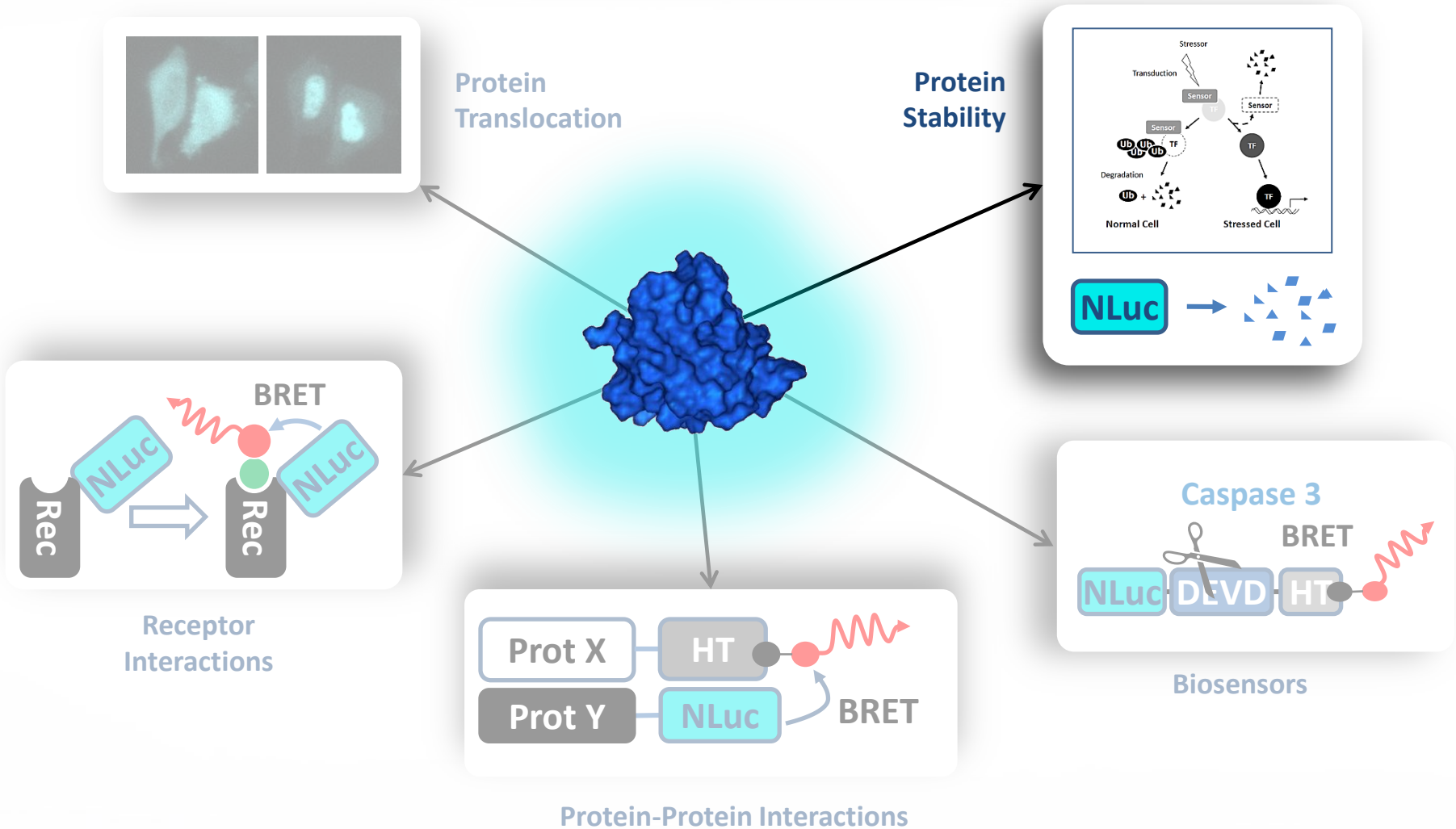


Time lapse: 13 minutes

NanoLuc fusion to
Glucocorticoid Receptor

HeLa cells; 500nM dexamethasone treatment

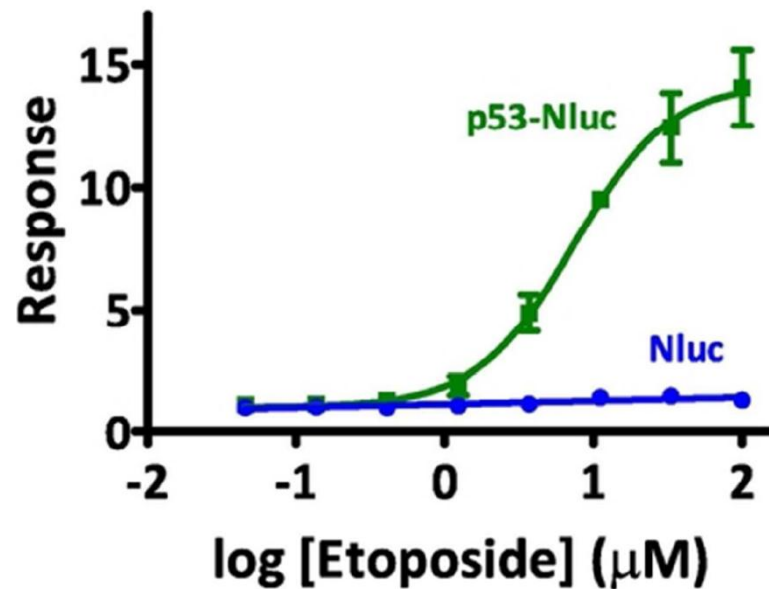
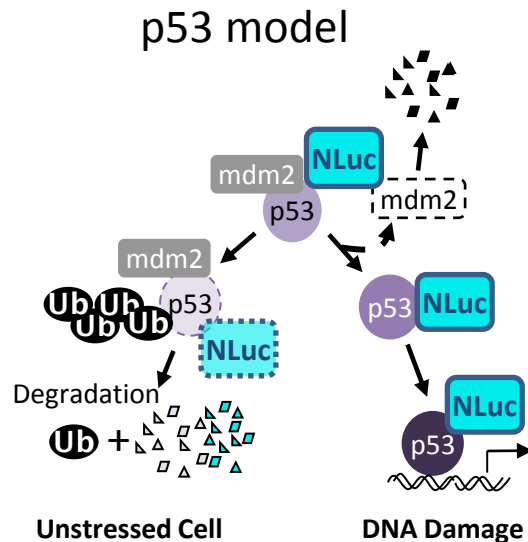
NanoLuc™ Luciferase as a fusion partner: Proof of concept experiments



Monitoring Protein Stability with NanoLucTM Luciferase

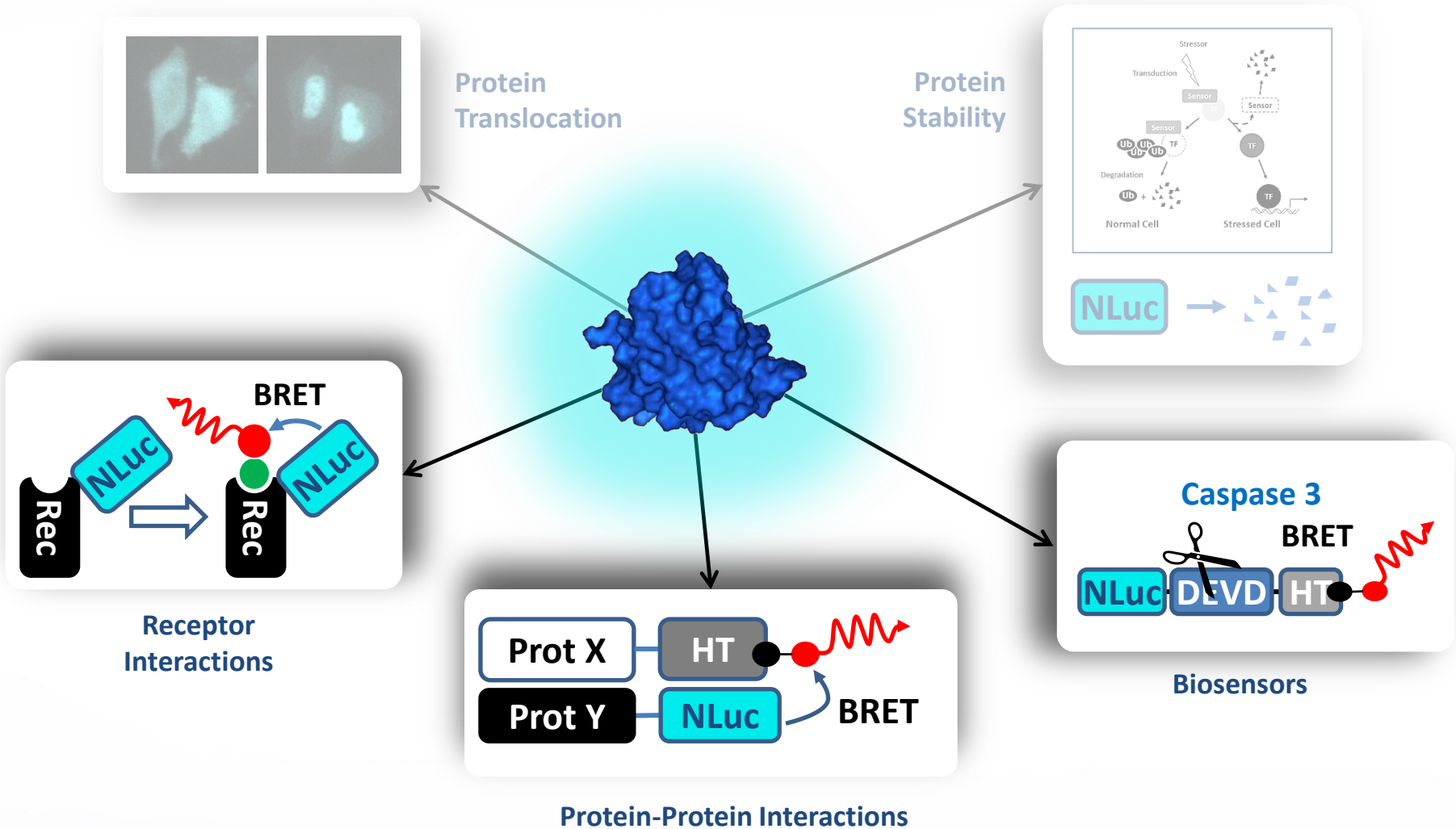


Can NanoLucTM Luciferase be added to a protein as a probe for protein stability?



The fusion can be used as a probe of stability

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

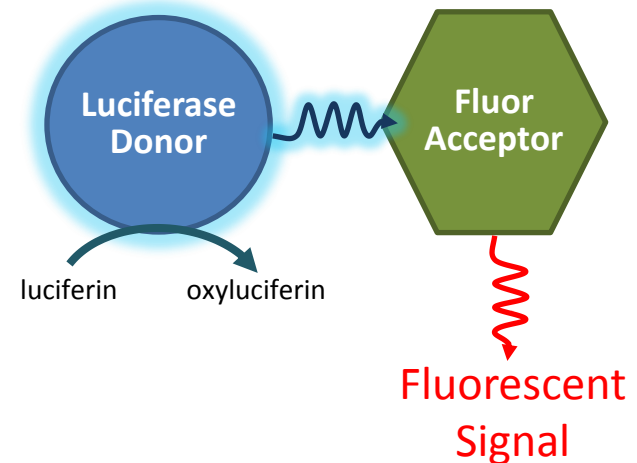
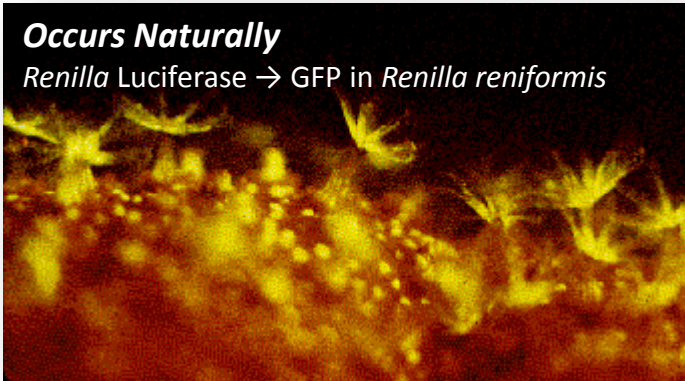


Bioluminescence Resonance Energy Transfer (BRET)



Occurs Naturally

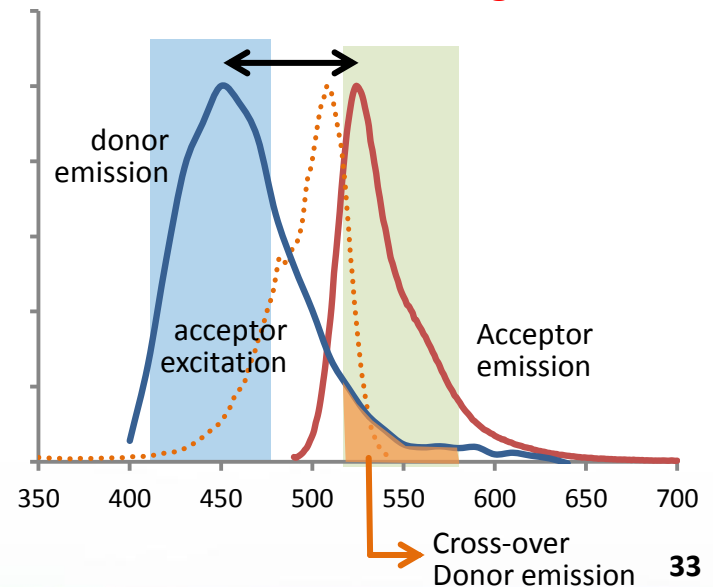
Renilla Luciferase → GFP in *Renilla reniformis*



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

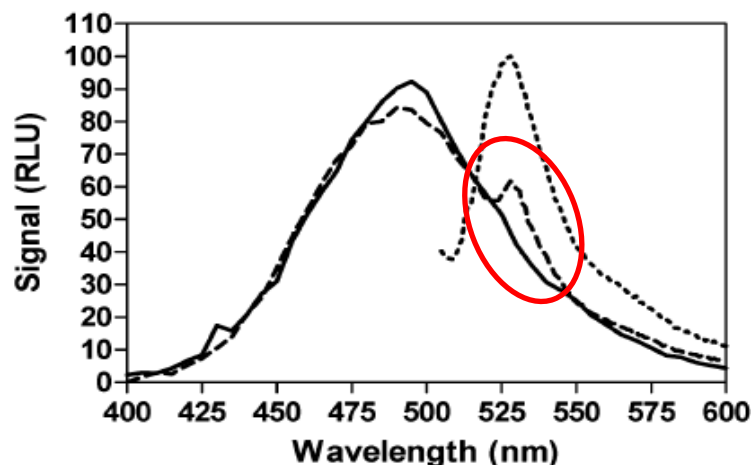


Could NanoLuc™ Luciferase work better as the luminescent donor?



RLuc → GFP

Donor brightness is a key limiter to current BRET technologies.



More spectral overlap needed to get sufficient signal



BRET-beneficial aspects of NanoLuc Luciferase:

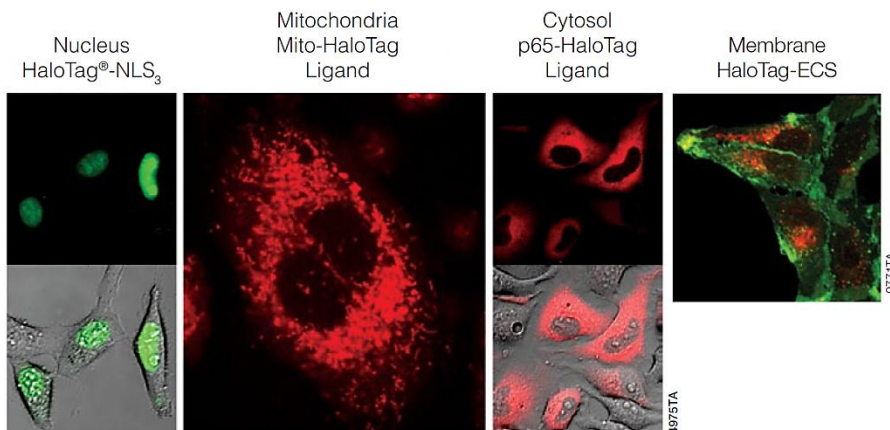
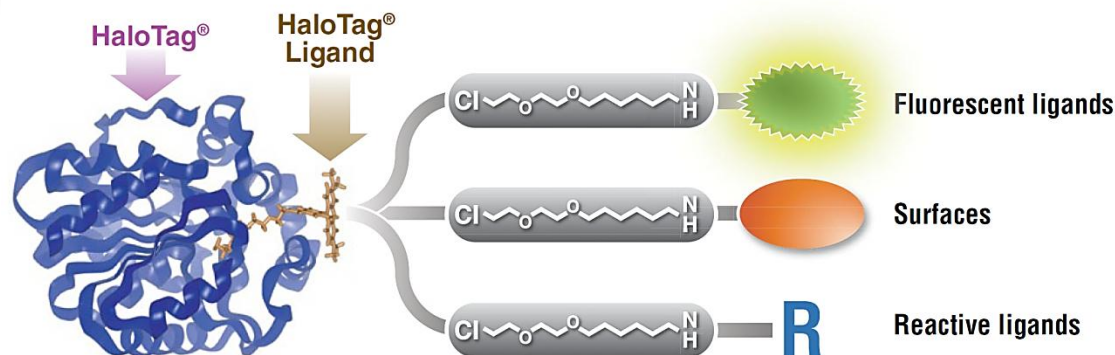
- ~100-fold brighter than RLuc
- ✓ need less spectral overlap with fluor
- ✓ gain greater spectral separation

We have a potential acceptor fusion protein:

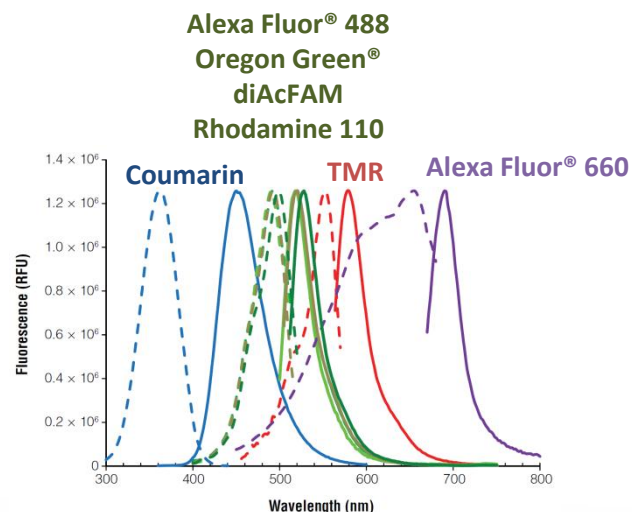
HaloTag® Fusion Protein

34.1kDa protein engineered from halophilic bacterial hydrolase.

- Engineered to lock into enzyme: substrate intermediate for **covalent attachment**.
- No homolog in mammalian cells.

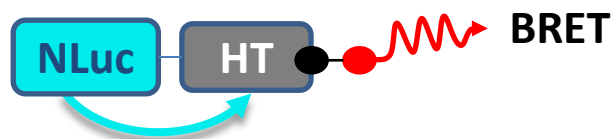


Goes anywhere in the cell

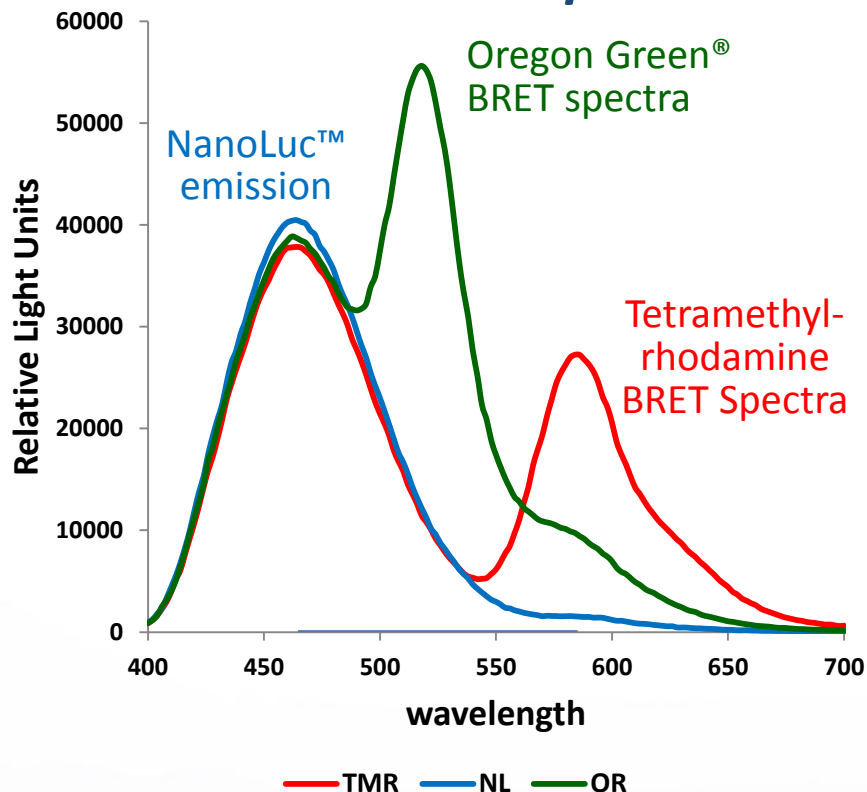


Variety of fluors ready-to-use

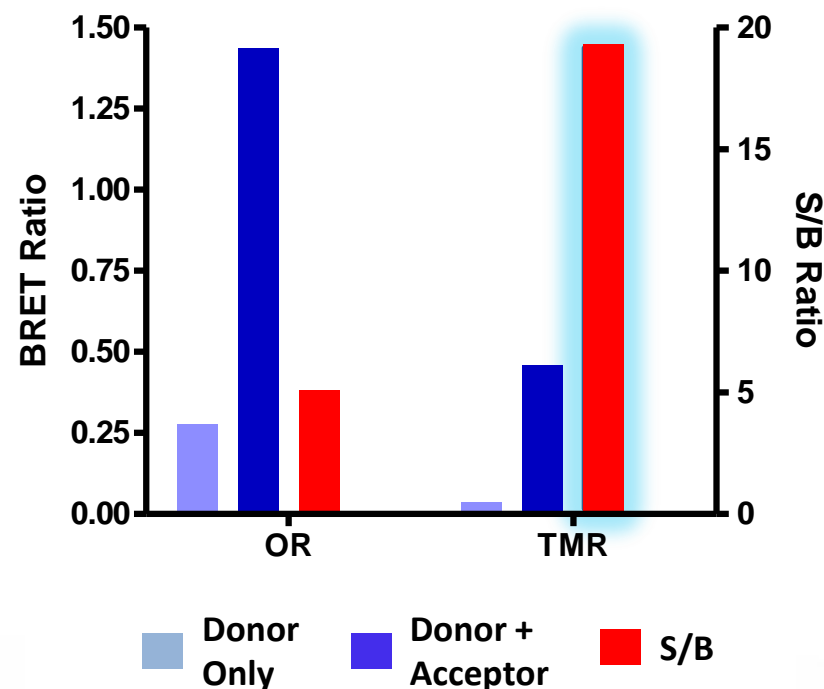
Testing the concept: NanoLuc™ & HaloTag® Fusion can perform BRET



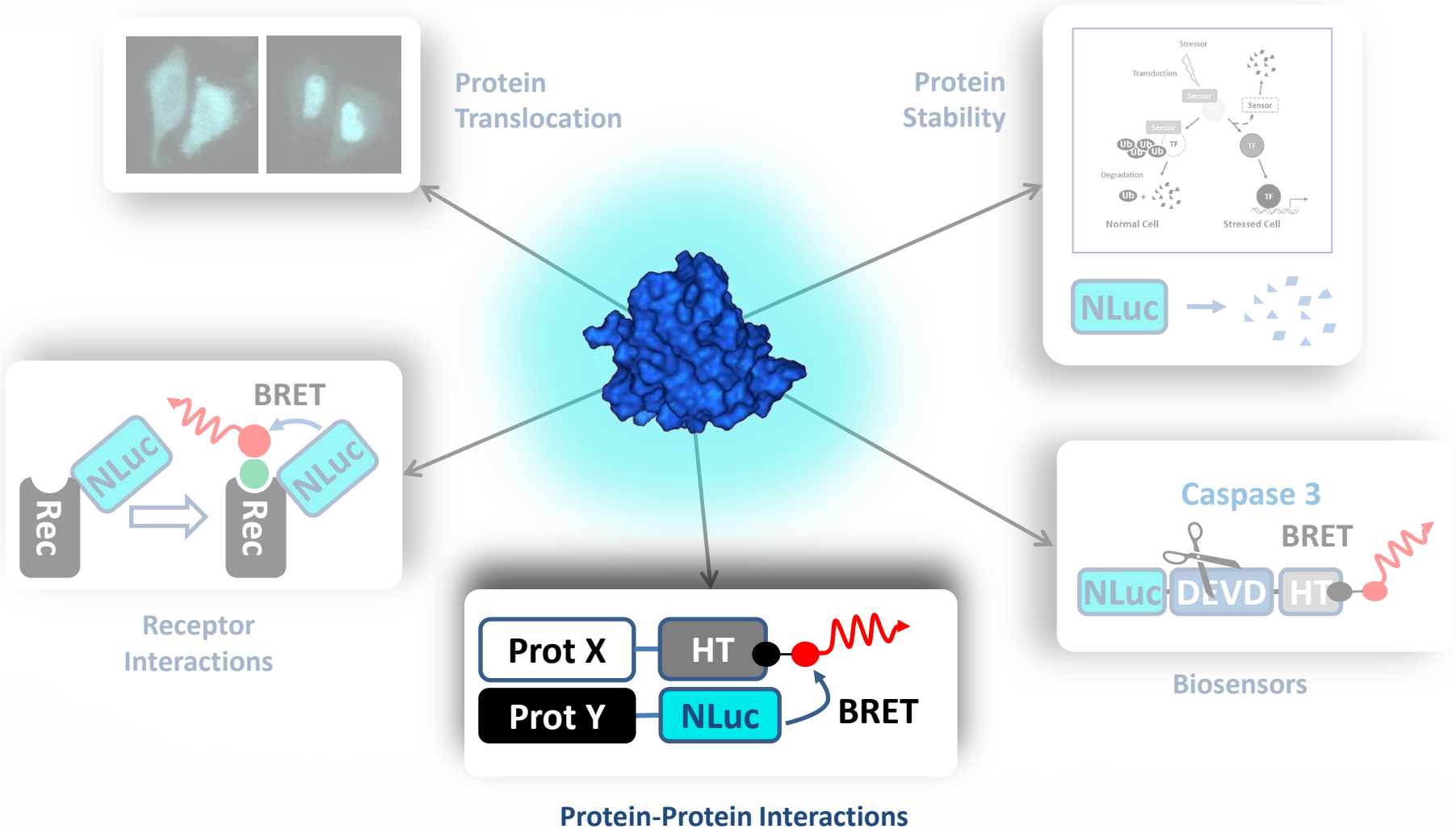
BRET Emission spectrum



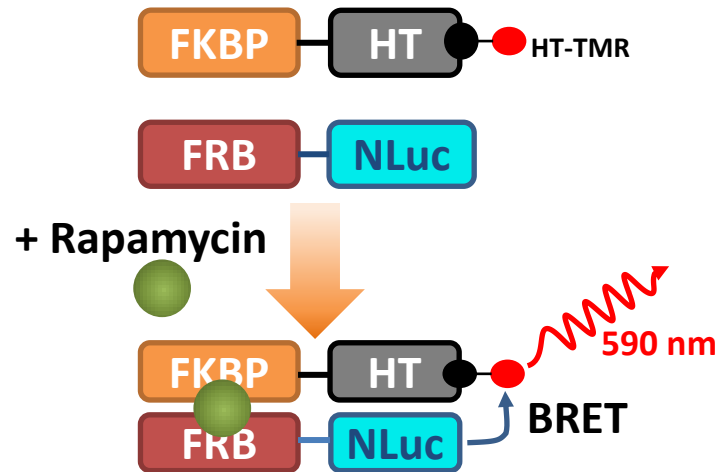
BRET Ratio ($= Em_{acc}/Em_{don}$)



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



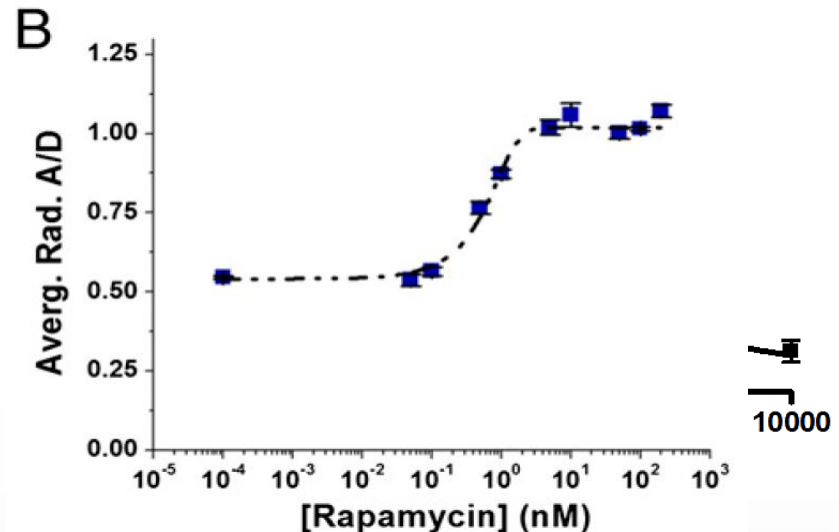
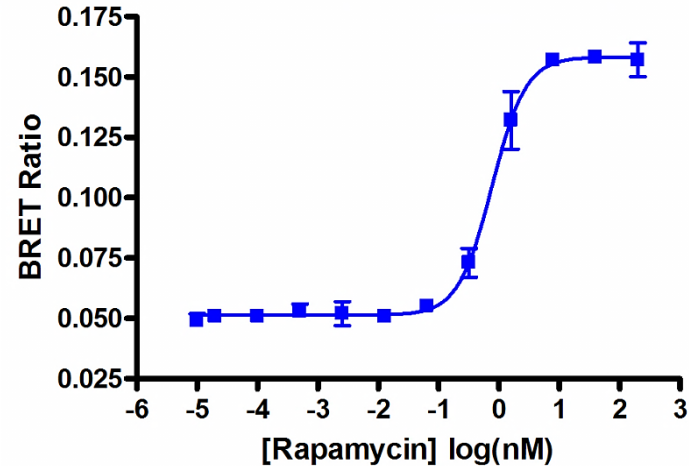
Can NLuc:HT Pair be used for Protein-Protein BRET?



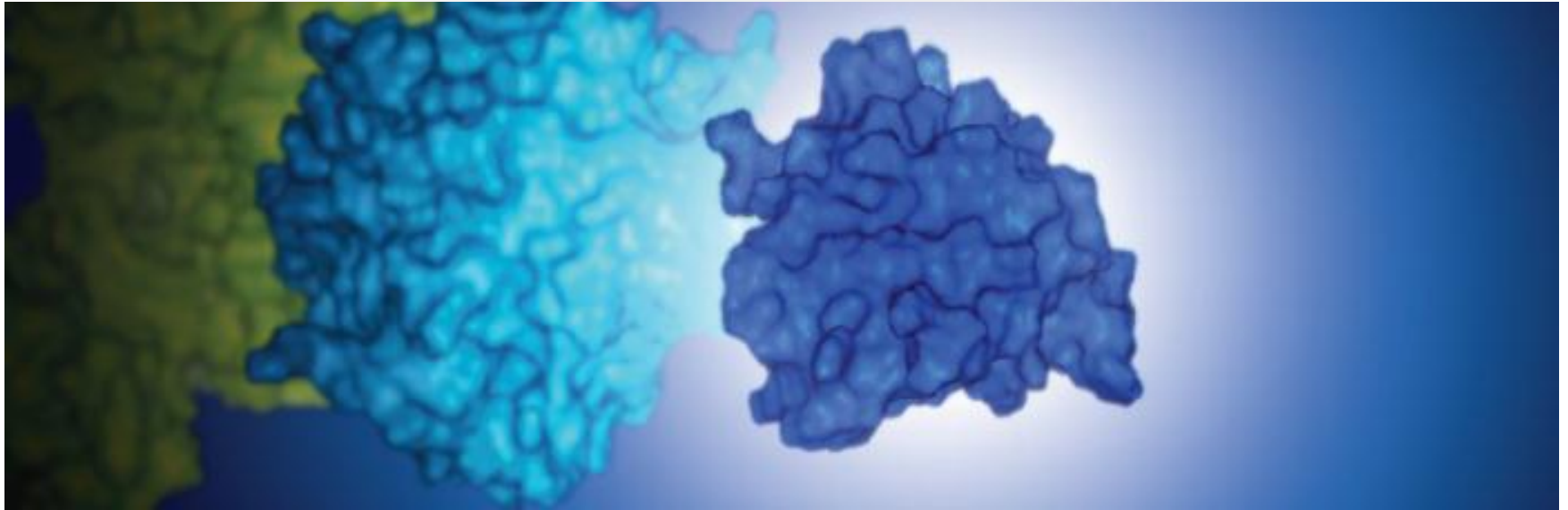
Same model system used with BRET 6 System

RLuc8.6 → TurboFP

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



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

***Get a FREE Sample of the
NanoLuc[™] Vectors.
Ask for the < Source Code >***



pNL 1.1 (*NLuc*)

pNL 1.2 (*NLucP*)

pNL 1.3 (*secNLuc*)

www.promega.com/nanoluc

Plus more information about all the pNL vectors

Technical Services Scientists ready to help



By phone: 800-356-9526

- Available 7am-6pm Central M-F

Online Chat @ promega.com

- Available 7am-6pm Central M-F
- Global Chat with Branch office tech serv scientists, too, after hours (language dependent)

techserv@promega.com

- Guaranteed answers within 24hr
- Most responses within 2hrs



North American Branch Technical Support Team

Questions?

