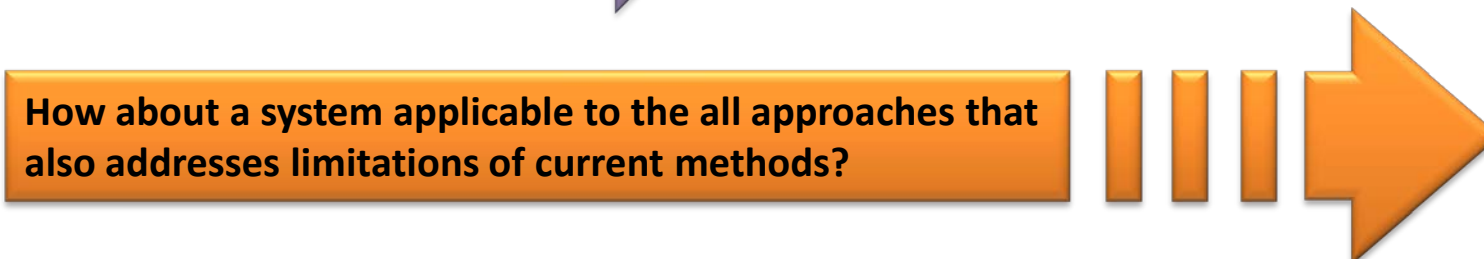
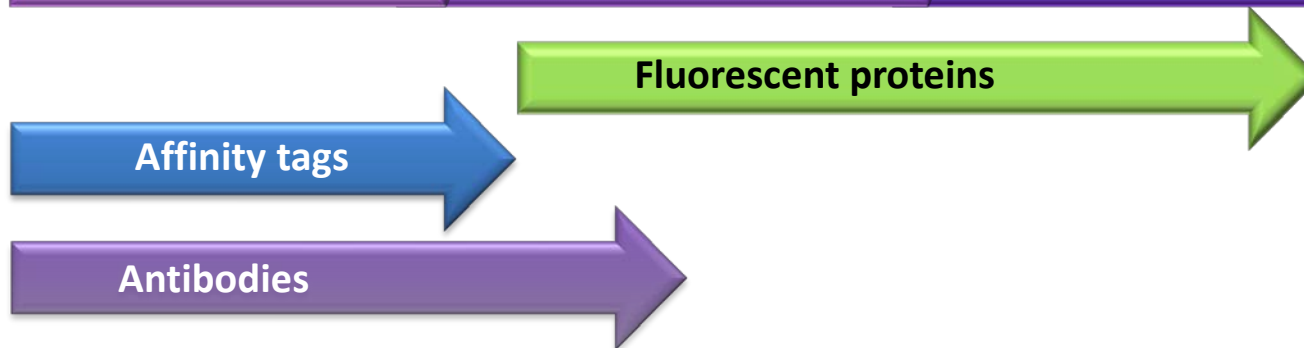


# ***Overcoming Challenges of Protein Analysis in Mammalian Systems***

***Danette L. Daniels, Ph.D.***



# Current Technologies for Protein Analysis



- Minimal interference with protein of interest
- Detection/real-time imaging
- High Signal/background
- Efficient capture/isolation
- Differential labeling

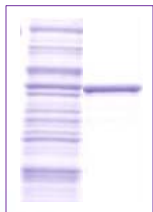
# HaloTag Platform



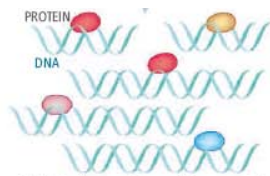
Protein purification  
Protein arrays  
Protein interactions

Protein localization  
Real time imaging  
Protein trafficking  
Protein turnover

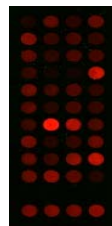
*In vivo* fluorescent imaging



**HaloTag®  
Purification**



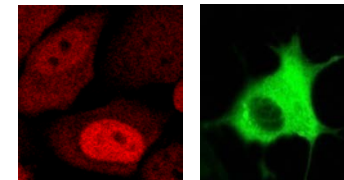
**HaloCHIP™  
Protein:DNA**



**HaloLink™  
Protein Arrays**

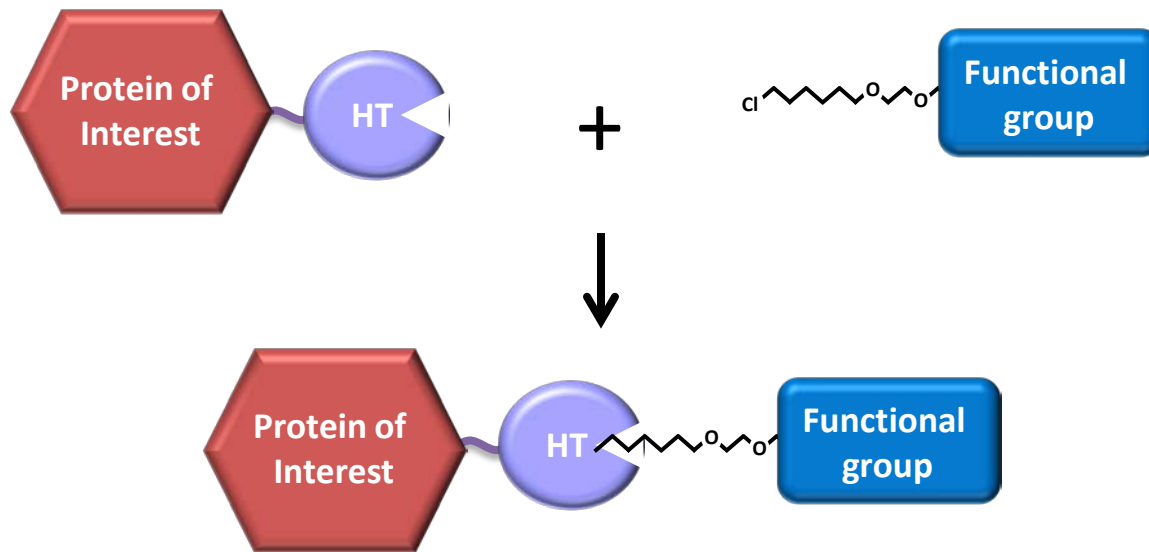


**HaloTag®  
Pull-Down**



**Fluorescent  
Ligands**

# *HaloTag is a Genetically Engineered Protein Fusion Tag*

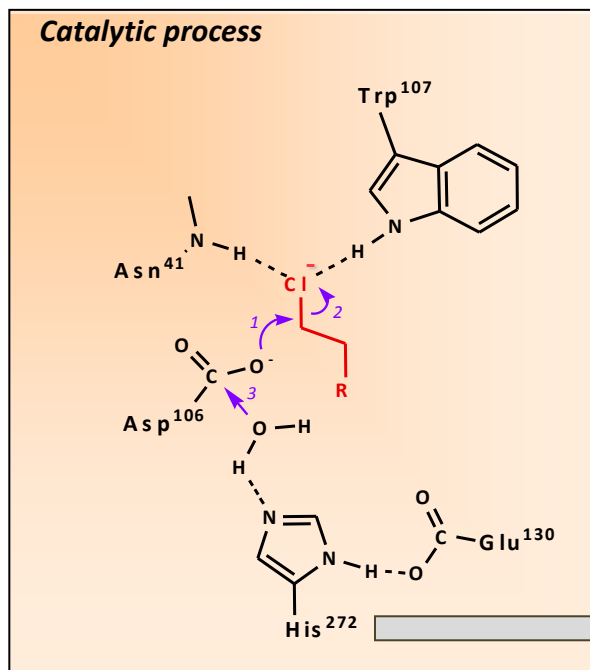


- A monomeric , 34 kDa, modified bacterial dehalogenase genetically engineered to covalently bind specific, synthetic HaloTag<sup>®</sup> ligands
- Irreversible, covalent attachment of chemical functionalities
- Suitable as either N- or C- terminal fusion

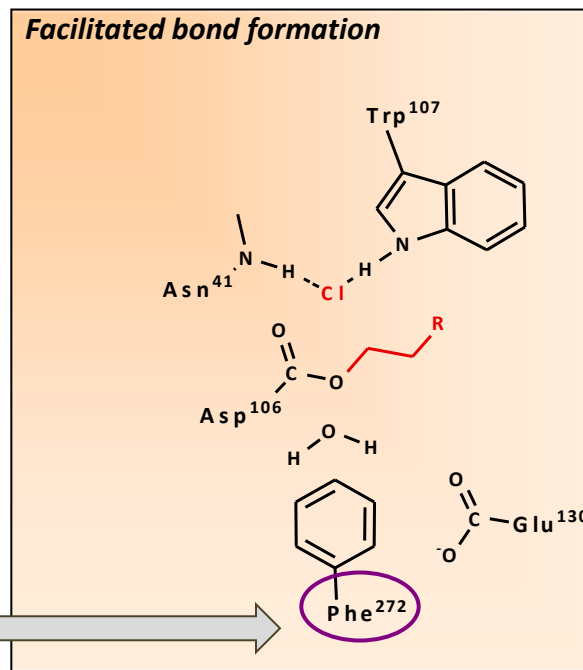
# Mutagenized HaloTag<sup>®</sup> Protein Enables Covalent HaloTag<sup>®</sup>-Ligand Complex



## Hydrolase (DhaA)



## HaloTag<sup>®</sup>



## HaloTag<sup>®</sup>:

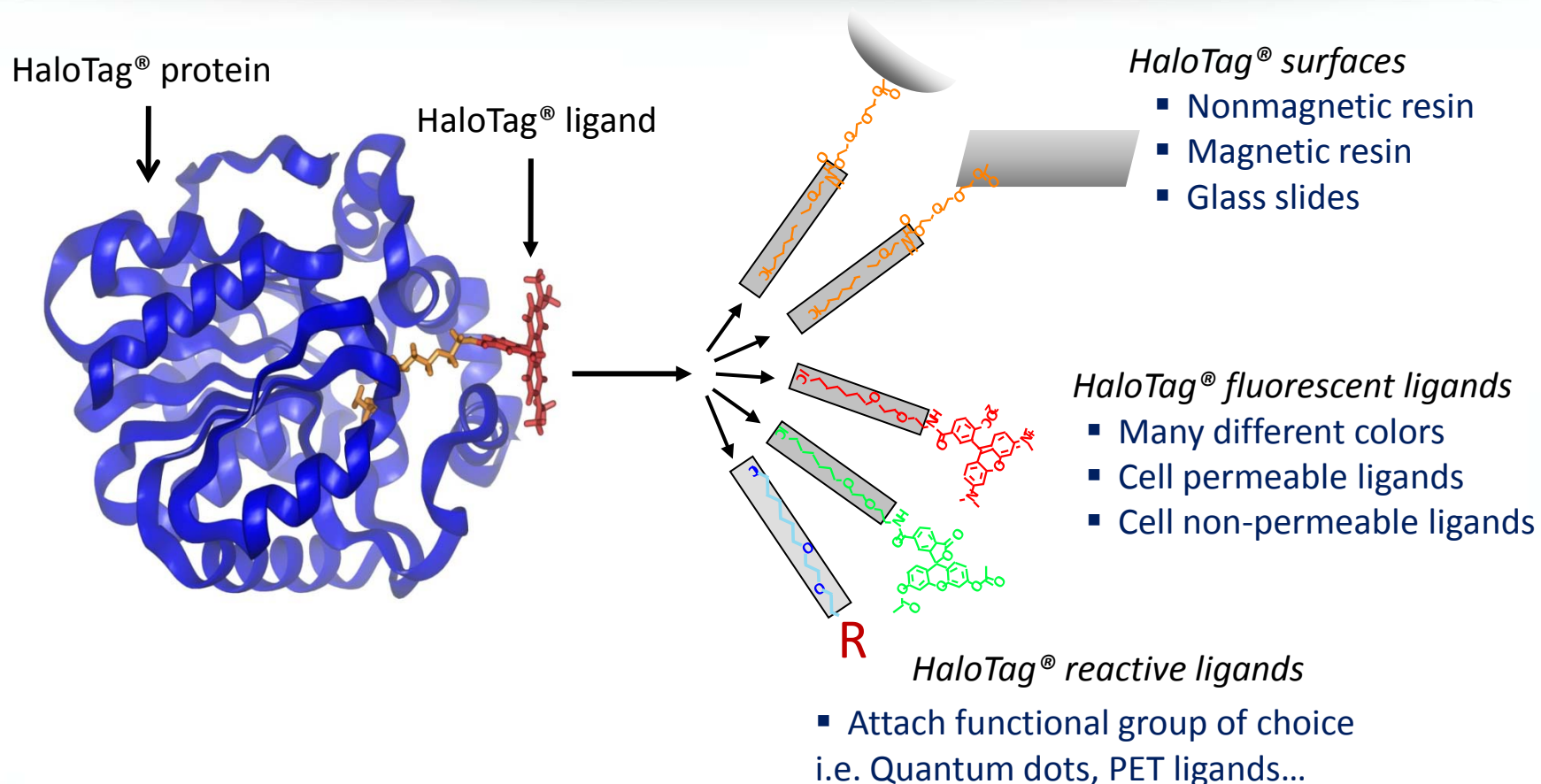
- 34kDa protein
- Monomeric
- Single change: His272Phe for covalent bond.

## Covalent bond:

- Stable after denaturation.
- Confirmed by Mass spec.

- DhaA is a rare, bacterial hydrolase.
- Binds to chloroalkane substrates.
- **Forms a covalent intermediate.**
- Activation of water by His drives hydrolysis.

# Ligands Impart Multi-functionality



- **Selectable functionalities:** a single fusion construct may be attached to a broad range of functional properties

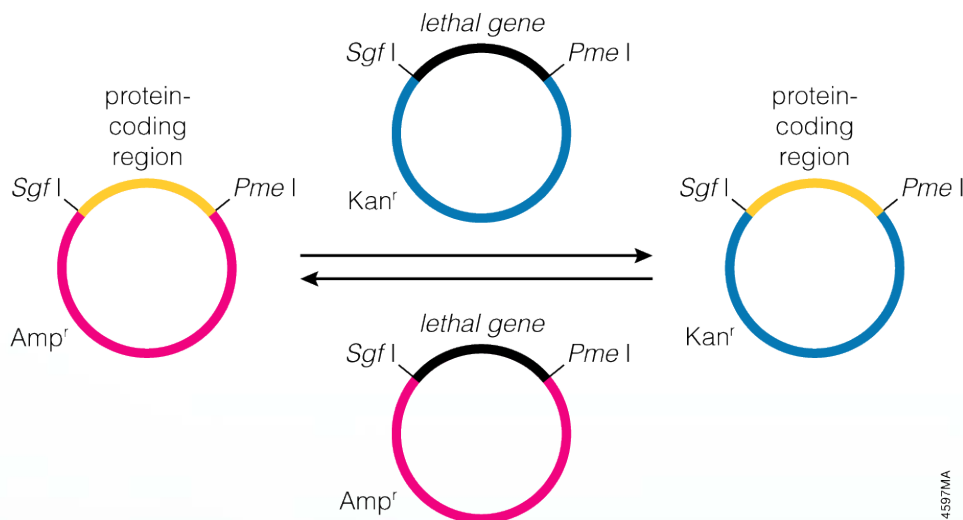
# Generation of a N- & C-Terminal HaloTag® Fusions



Flexi vectors for expression of HaloTag fusions in mammalian cells:

Expression	Promoter	N-terminal HaloTag®	C-terminal HaloTag®
Maximal expression	CMV	pFN21	pFC14
Promoter deletion series to optimize mammalian expression level	CMVd1	pFN22	pFC15
	CMVd2	pFN23	pFC16
	CMVd3	PFN24	pFC17

## Flexi® Vector cloning system



- Flexible system for directional cloning that utilizes REs that are infrequent in ORFs
- Efficient transfer to multiple vectors  
Sequence once, transfer to many



# No Cloning Necessary

## HaloTag<sup>®</sup>-Fusion Clones are Readily Available

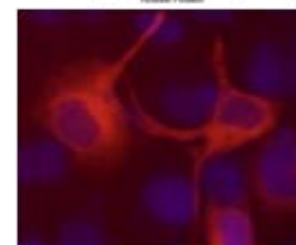
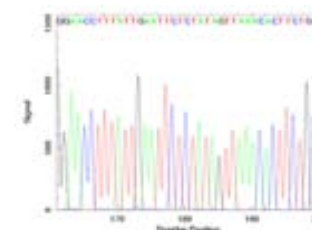
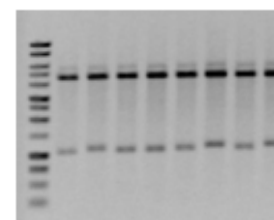


### Kazusa DNA Research Institute

Human ORFs N-terminal fusion constructs in Flexi<sup>®</sup> vector pFN21A for expression in mammalian cells

<http://www.kazusa.or.jp/kop/dsearch-e//>

Features	Flexi-HaloTag Collection PID beginning with FHC as of Nov, 2010
Size of Collection	7,100 clones
Fusion Tag	HaloTag <sup>®</sup> <small>for protein purification, imaging &amp; interactions</small>
<b>Validated Clones</b>	
> Sequence Validated	<b>Yes</b> (100% clones)
> Insert Validated	<b>Yes</b> (99.7% clones)
> Expression Validated	<b>Yes</b> (99.8% clones)
> Localization Validated	<b>Yes</b> (80.1% clones)
Format	DNA
Typical Delivery	2-4 weeks
Price (\$USD)	\$500.00 per clone

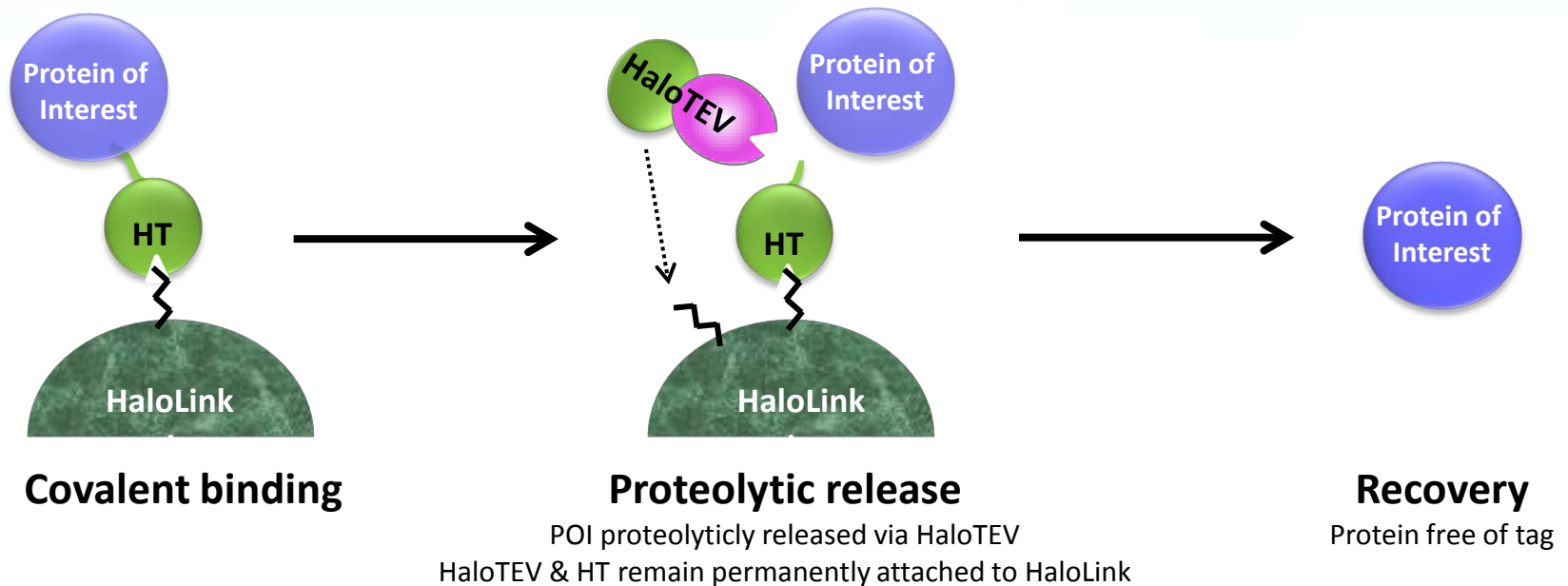


### GeneCopoeia

Human and mouse ORFs N- /C- terminal fusion constructs available in OmicsLink<sup>™</sup> vectors for expression in mammalian cells <http://www.genecopoeia.com/product/halo/>



# HaloTag<sup>®</sup>-based Protein Purification Scheme



## ❑ Covalent binding:

- Efficient capture regardless of expression level
- Stringent washes possible
- Minimal loss of bound protein

## ❑ Streamlined protocol:

- Proteolytic release coupled with protease & tag removal
- One physiological buffer and no need for buffer exchange

# *Protein Purification from Mammalian Cells*

## *Efficient, Sensitive & Gentle*



### **Mammalian cells: high quality proteins**

- Native environment
- Proper folding
- Protein processing
- Correct post-translation modifications

**Low expression level : Low yields; Low recovery; Impurities**

### **HaloTag<sup>®</sup>:**

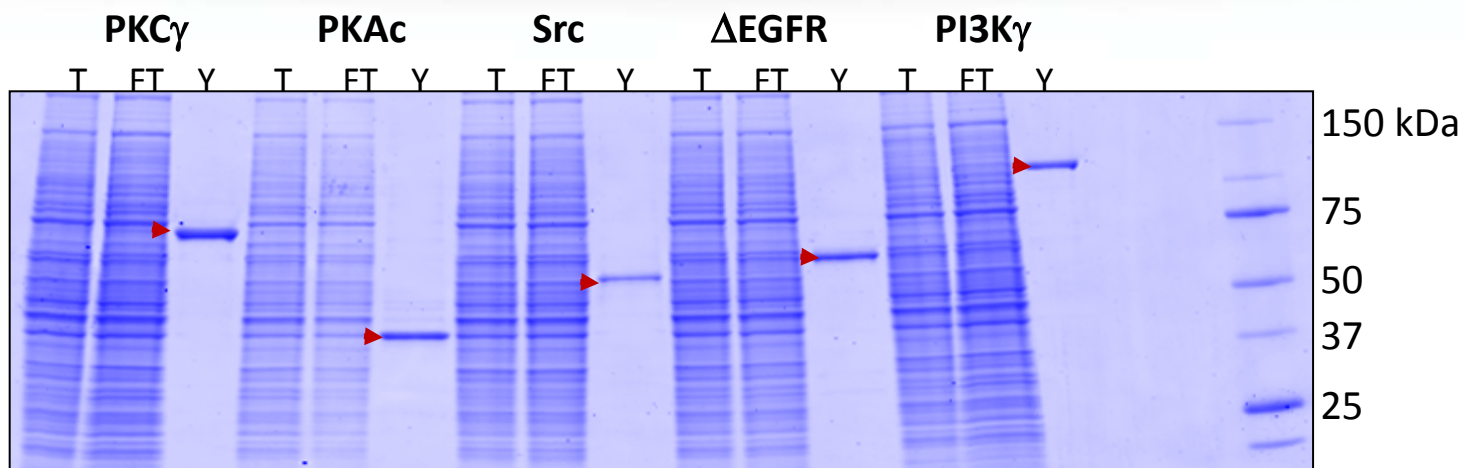
#### **Selective & Covalent capture**

- Efficient protein capture regardless of expression levels
- No loss of bound protein during washes

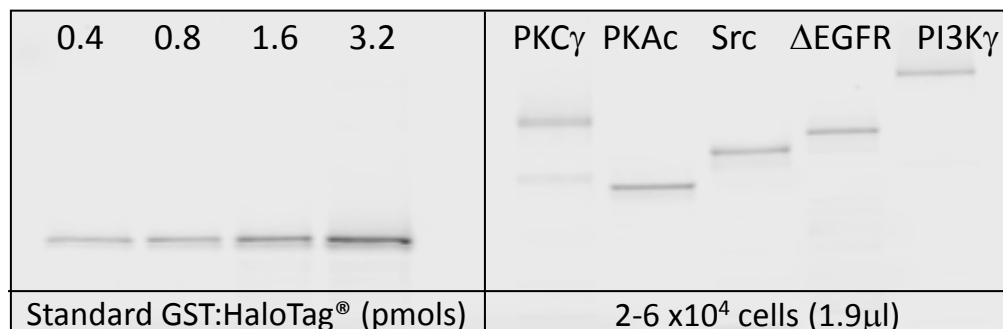
#### **Rapid sensitive detection**

- Optimization of expression levels

# Purification of Human Kinases from Transient Transfected HEK293T Cells

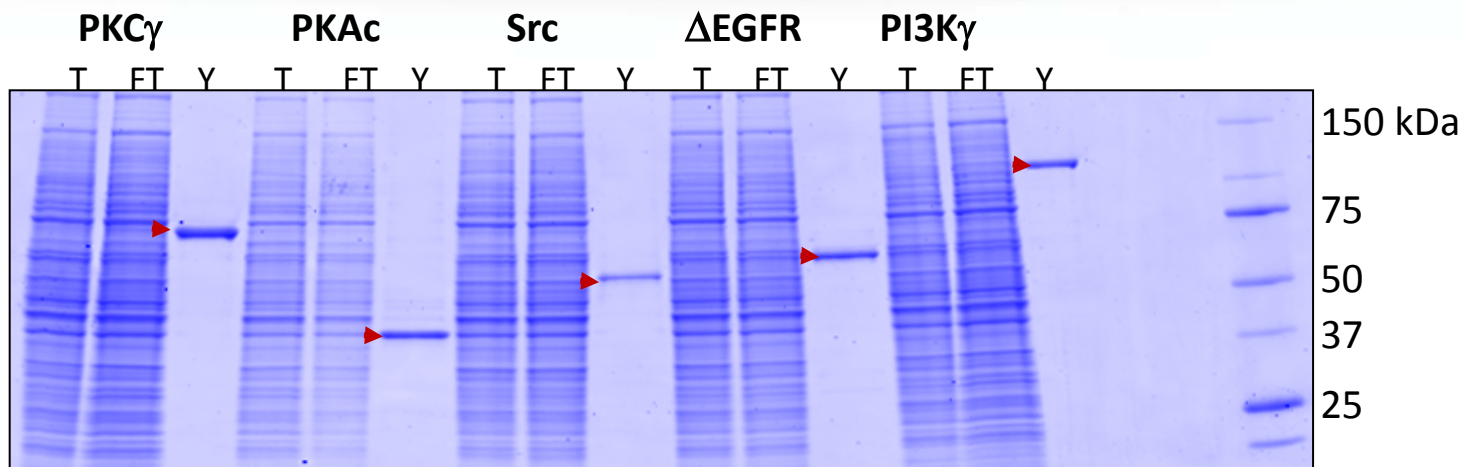


T: starting material    FT: unbound    Y: yield



Protein standard and fusions were labeled with HaloTag<sup>®</sup> TMRDirect ligand

# Purification of Human Kinases from Transient Transfected HEK293T Cells

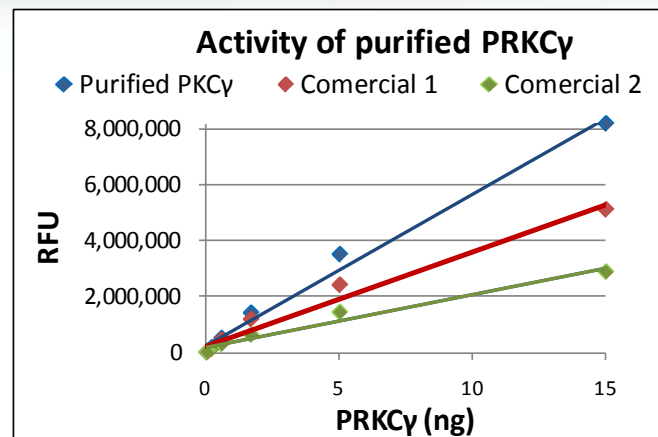
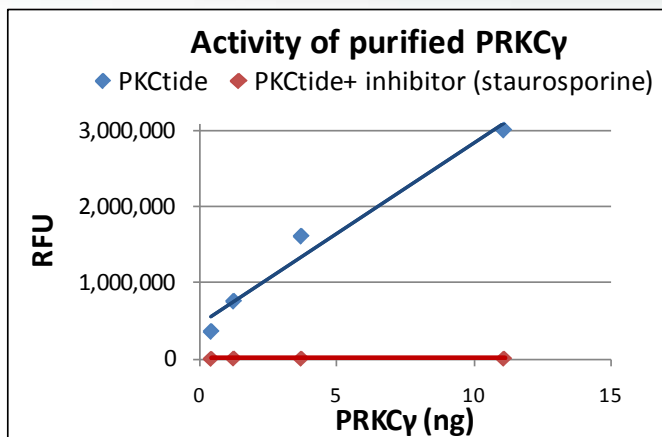


T: starting material FT: unbound Y: yield

Kinases	PKC $\gamma$	PKAc	SRC	$\Delta$ EGFR	PI3K $\gamma$
Estimated POI expression ( $\mu$ g)	289	159	181	177	247
Yield ( $\mu$ g)	244	140	137	167	221
% Recovery	84%	88%	76%	94%	89%

**Highly efficient protein capture and recovery**

# Purified Kinases are Highly Active

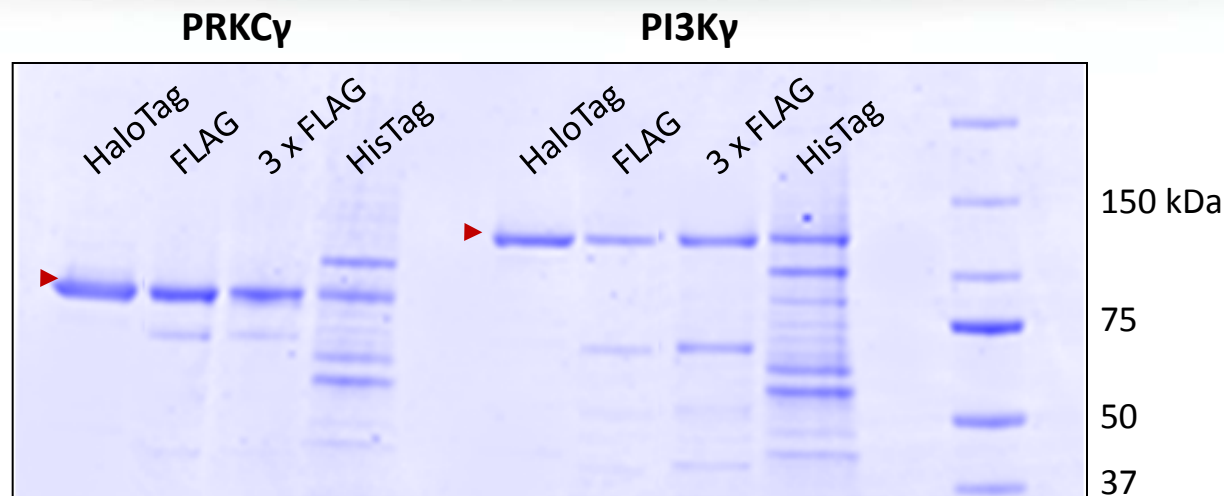


Kinase	Measured Specific activity (nmol/min/mg)	Reported Specific activity (nmol/min/mg)
PRKCy	16,551	2,260
PKAc	9,670	8,580
Src	1,624	1,032
$\Delta$ EGFR	196	101
PI3K $\gamma$	233	39

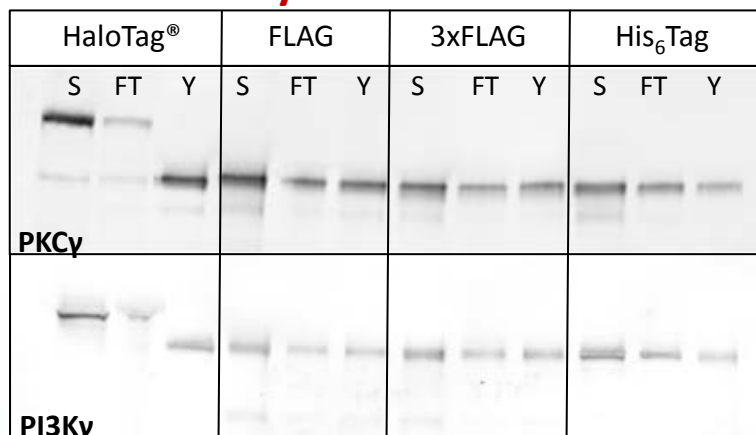
**Measured specific activities are in agreement with reported values**

Kinase activity was assayed using ADP-Glo™ assay

# Comparative Analysis with Other Affinity Tags



## Western analysis



T: starting material FT: unbound Y: yield

## Protein recovery (%)

Tag	PKC $\gamma$	PI3K $\gamma$
HaloTag <sup>®</sup>	86	88
FLAG	60	40
3xFLAG	52	58
His <sub>6</sub> Tag	31	33

**HaloTag<sup>®</sup>: Greater protein recovery  
Higher yields  
Higher purity**

## ***Summary: HaloTag<sup>®</sup>-Based Purification***

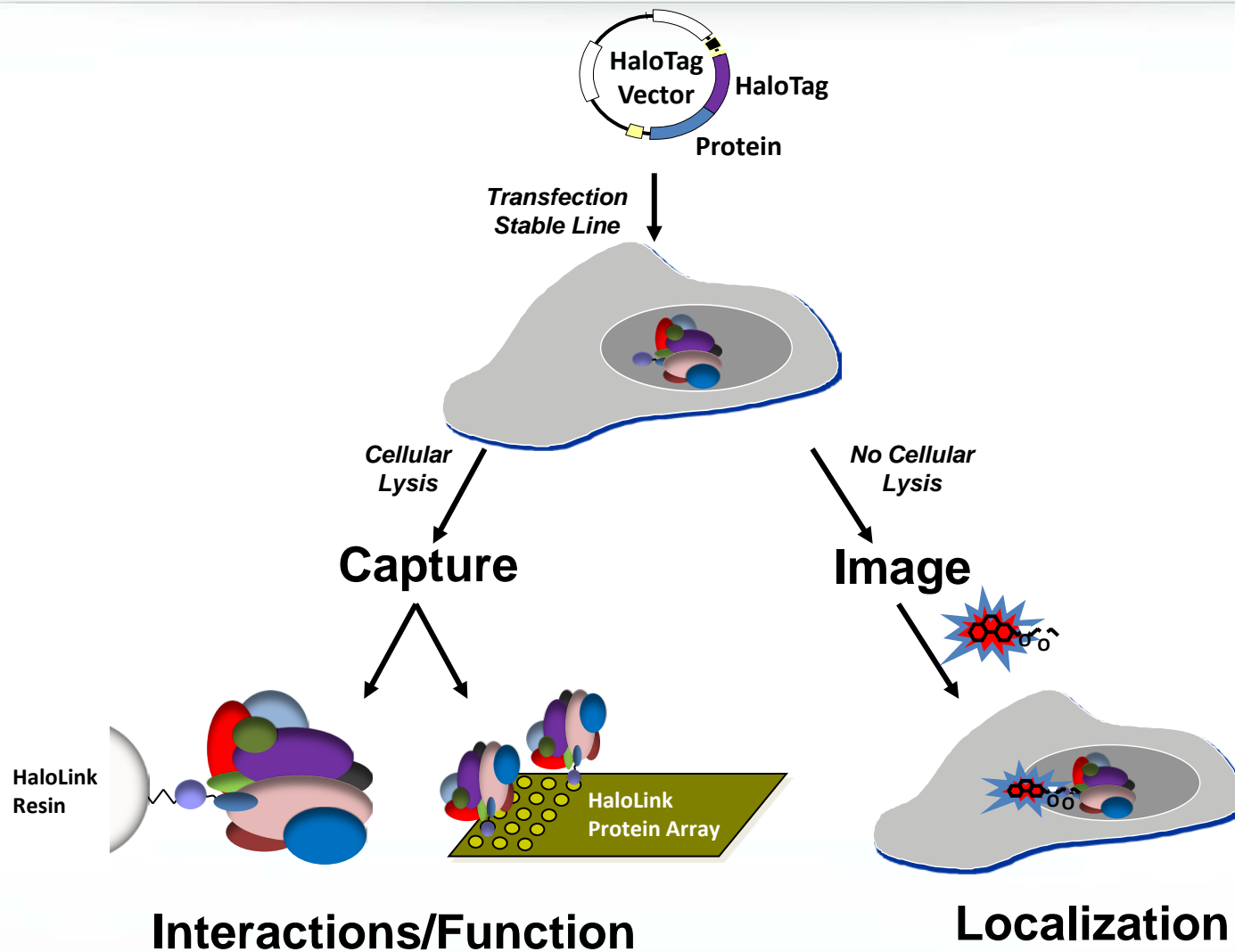
### **HaloTag<sup>®</sup>-based protein purification**

- Highly efficient purification regardless of expression levels
- Greater yields, purity and recovery than traditional affinity tags
- Streamlined protocol for proteolytic release coupled with protease & tag removal
- One physiological buffer and no need for buffer exchange

**Simple to use fluorescent detection for rapid optimization of expression levels**



# Studying Intracellular Protein Function

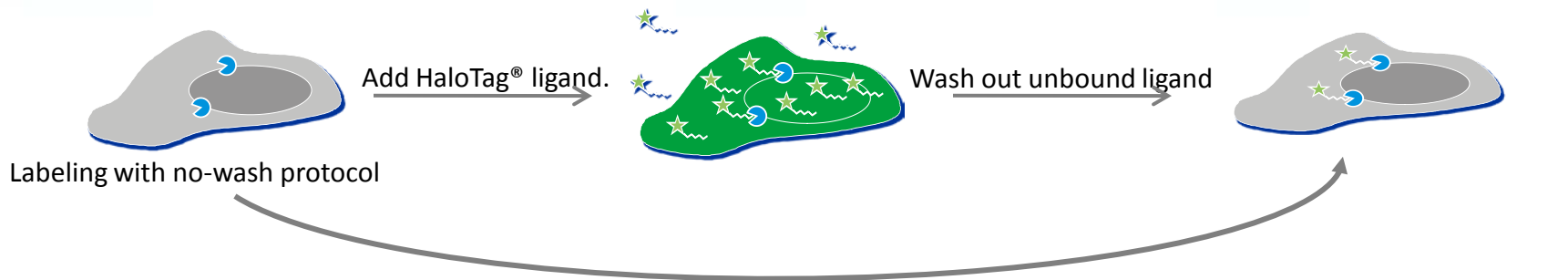


# Intracellular Protein Labeling and Imaging



Transfect HaloTag®-fusion construct into cells

Image; live or fixed cells

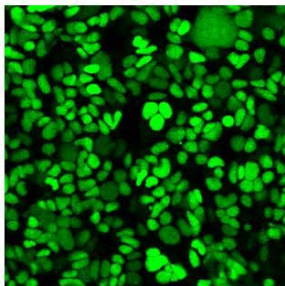


- No cytotoxicity
- Stable labeling
- Multiple colors
- Rapid and easy protocol

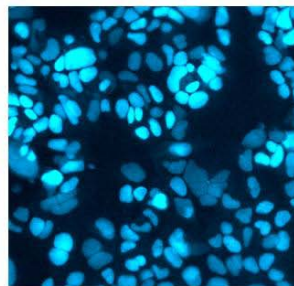
HaloTag®-NLS<sub>3</sub> fusion protein – HEK293 cells



TMR

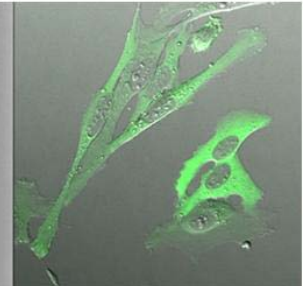
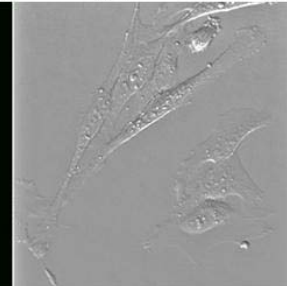
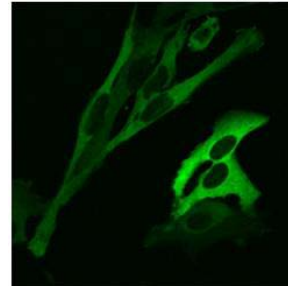


dAc-FAM



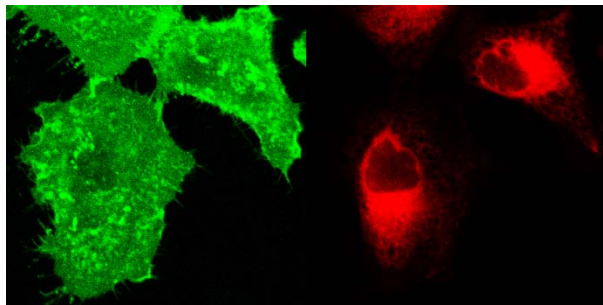
Coumarin

U2OS cells stably expressing p65-HaloTag.

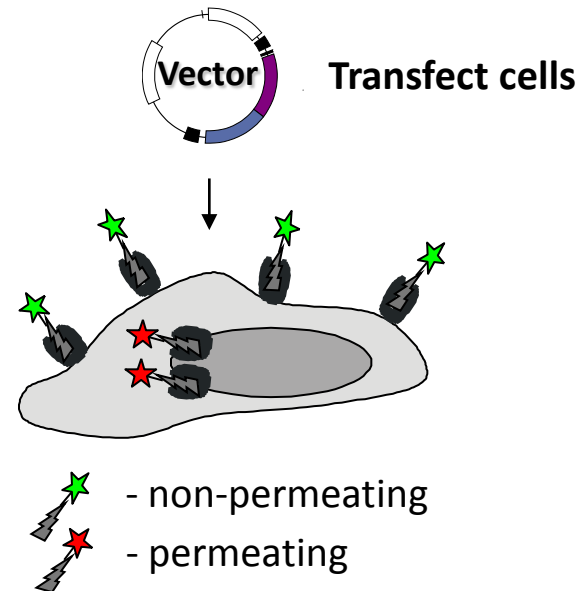
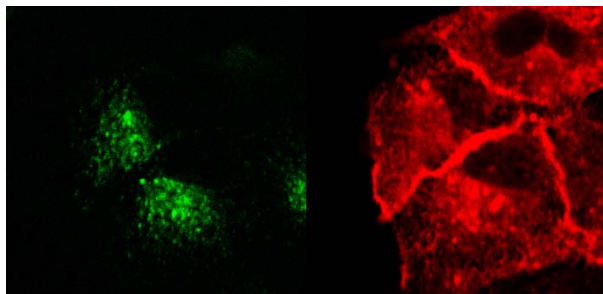


- All that is done with GFP can be done with HaloTag® and more...

# Analysis of Protein Trafficking



↓ 12hrs

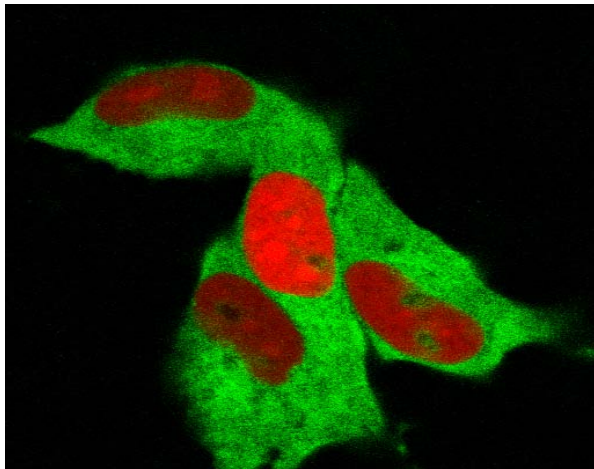


- Spatial control of labeling.
- Follow protein trafficking of distinct protein pools.

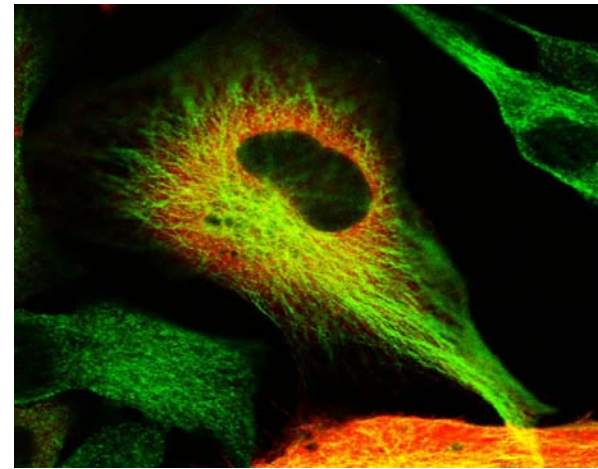
# Multiplexing with Other Labeling Technologies



hMGFP- $\alpha$ -tubulin  
HaloTag<sup>®</sup>-NLS<sub>3</sub>-TMR ligand



p65-HaloTag<sup>®</sup>-TMR Ligand  
Alexa 488 Ab



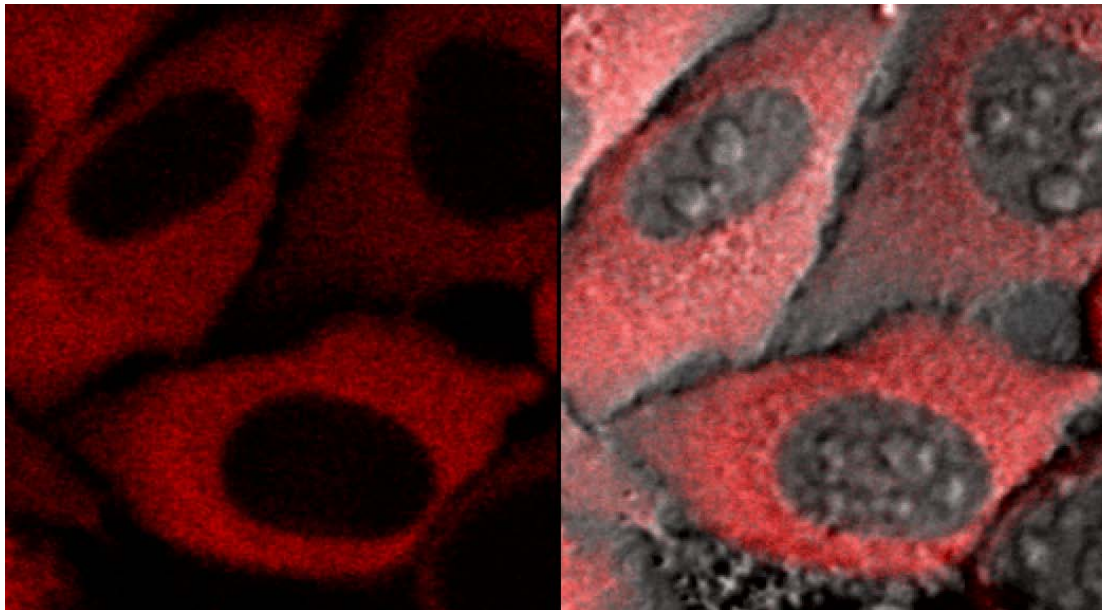
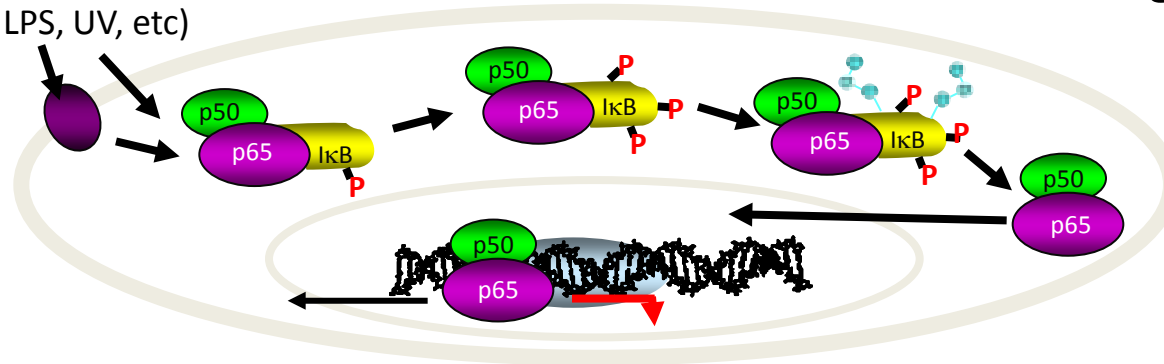
p65-HaloTag labeled with TMR Ligand  
fixed, then processed for ICC with anti- $\beta$ -tubulin Ab  
and Alexa-488-conjugated secondary antibody

- HaloTag<sup>®</sup> is compatible with fluorescent protein fusions
- HaloTag<sup>®</sup> is compatible with fixing and antibody staining
- Labeling simultaneous with fixation also possible

# Real-time Imaging

Signal (TNF $\alpha$ ,  
LPS, UV, etc)

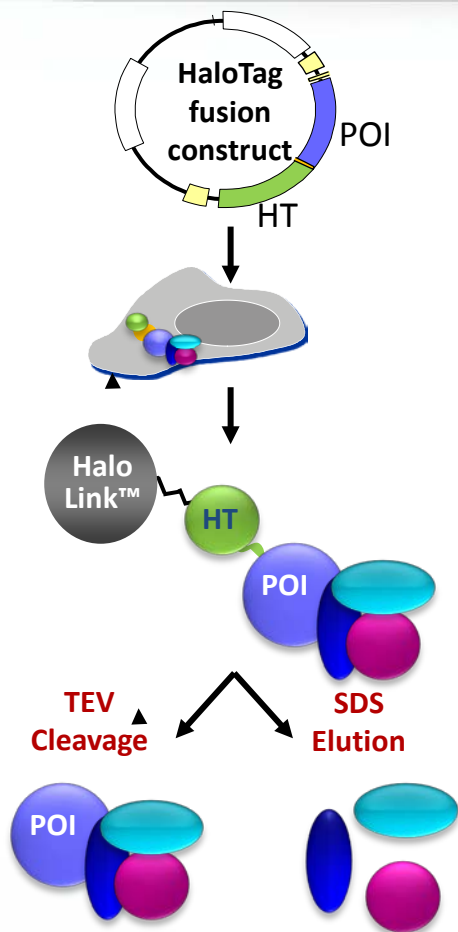
## NF $\kappa$ B Signaling Pathway



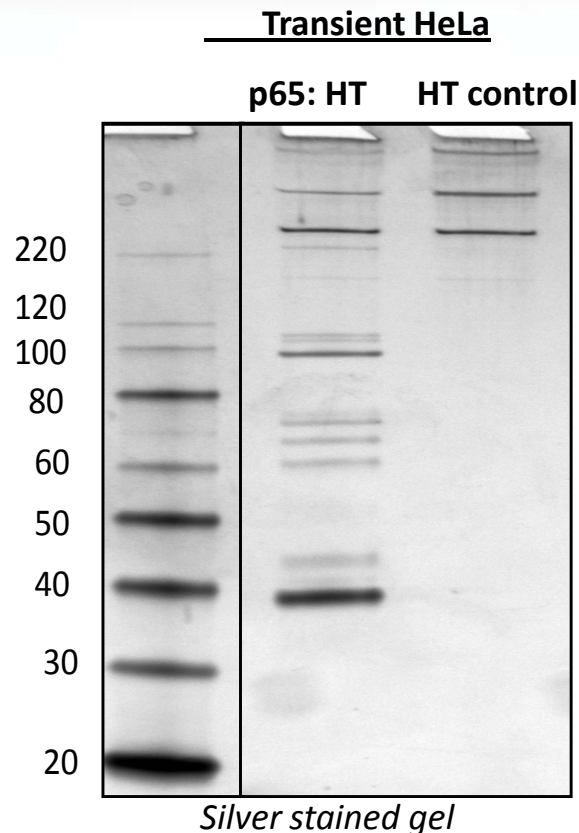
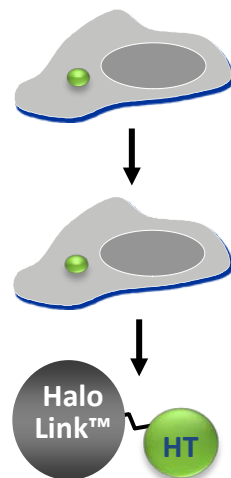
- HeLa cells expressing p65-HaloTag labeled with TMR Ligand
- Treated with TNF $\alpha$
- Imaged (5min/frame; 120min)



# Capture of Protein:Protein Complexes



Negative control:



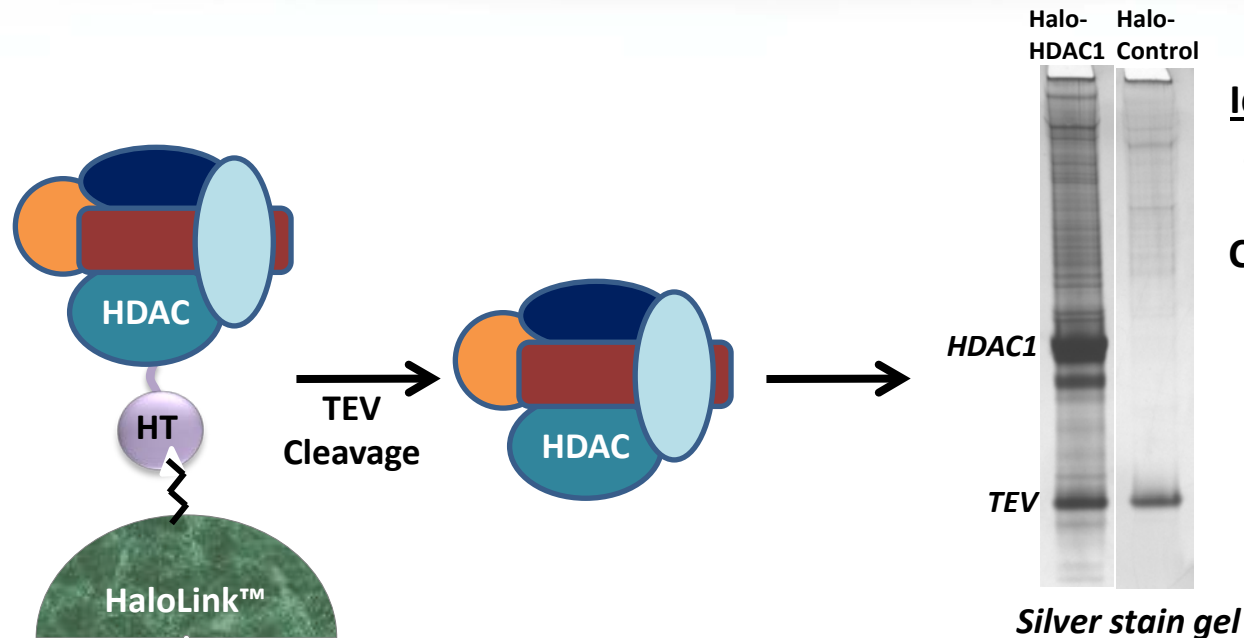
**Protein identified by MS analysis:**

p105 (p50)  
p100  
Rel A (p65)  
p50  
p52  
Rel B  
C-Rel  
I $\kappa$ B $\alpha$ , I $\kappa$ B $\beta$ , I $\kappa$ B $\epsilon$

\*Same results from TEV cleavage and in-solution digestion

•p65-HaloTag specifically pull-down expected protein partners of the NF $\kappa$ B pathway

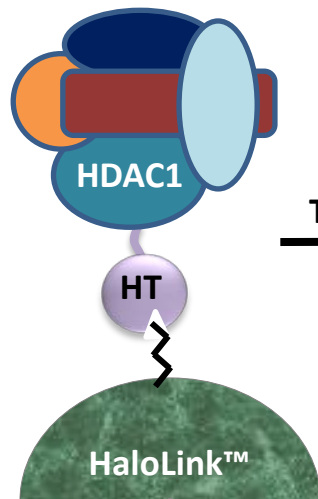
# HDAC1 Complex Purification



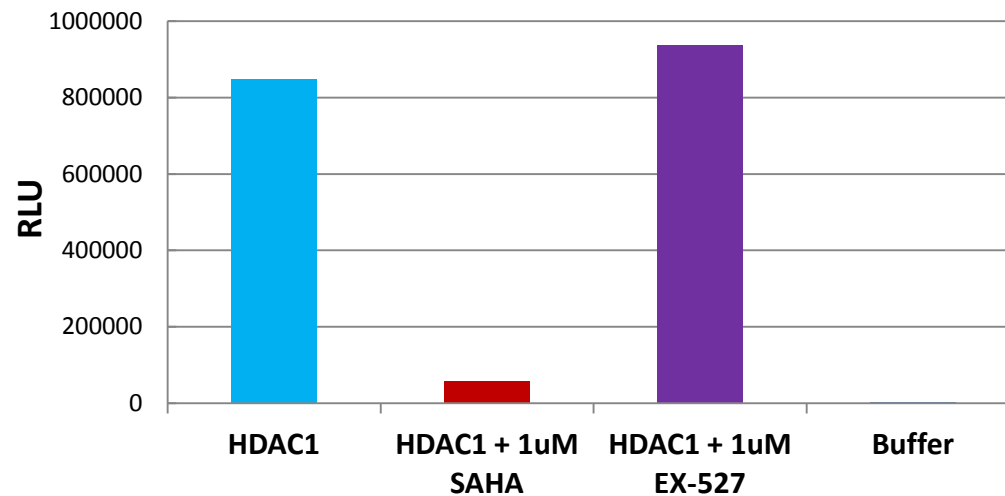
- Expected HDAC complex capture as determined by Mass Spec
- TEV cleavage allows for HDAC complexes to be released in tact.
- Compatible with downstream functional analysis



# Isolated HDAC1 Complexes Show Specific Activity



## *HDAC-Glo™ +/- HDAC (SAHA) and SIRT (EX-527) inhibitors*

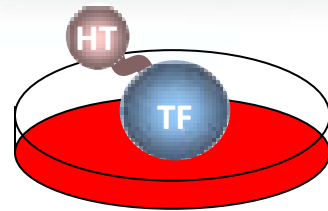


- Enrichment of specific HDAC activity from HDAC complex purification
- Able to screen effects of inhibitors on purified physiological complexes
- Overall technology extended to other epigenetic complex

# Intracellular Protein:DNA Interactions -HaloCHIP™

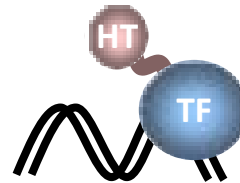


1) Expression of HaloTag® (HT) Trxn. factor (TF) fusion protein.

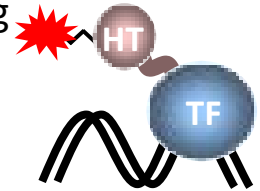


**Controls**  
Untransfected Cells

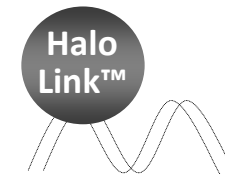
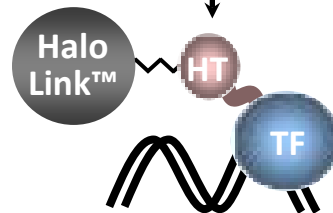
2) Crosslinking, lysis, and sonication.



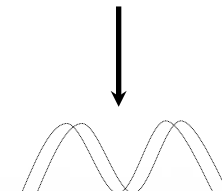
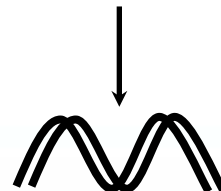
**OR**  
Block HaloTag® binding



3) Covalent capture on HaloLink™ resin followed by stringent washing.



4) Release of DNA by reversal of crosslinks.



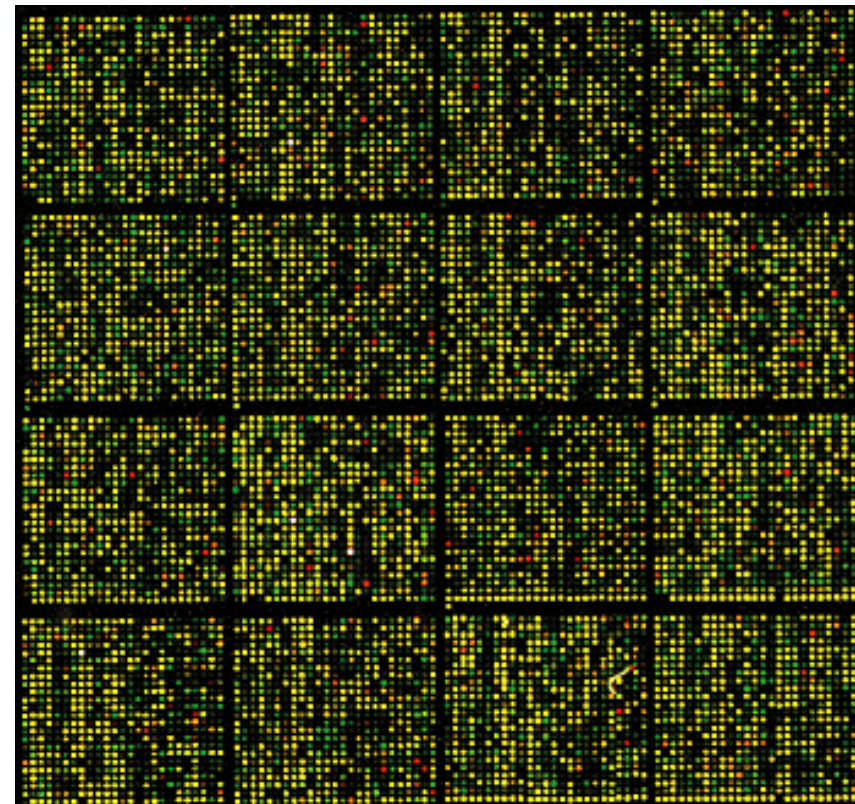
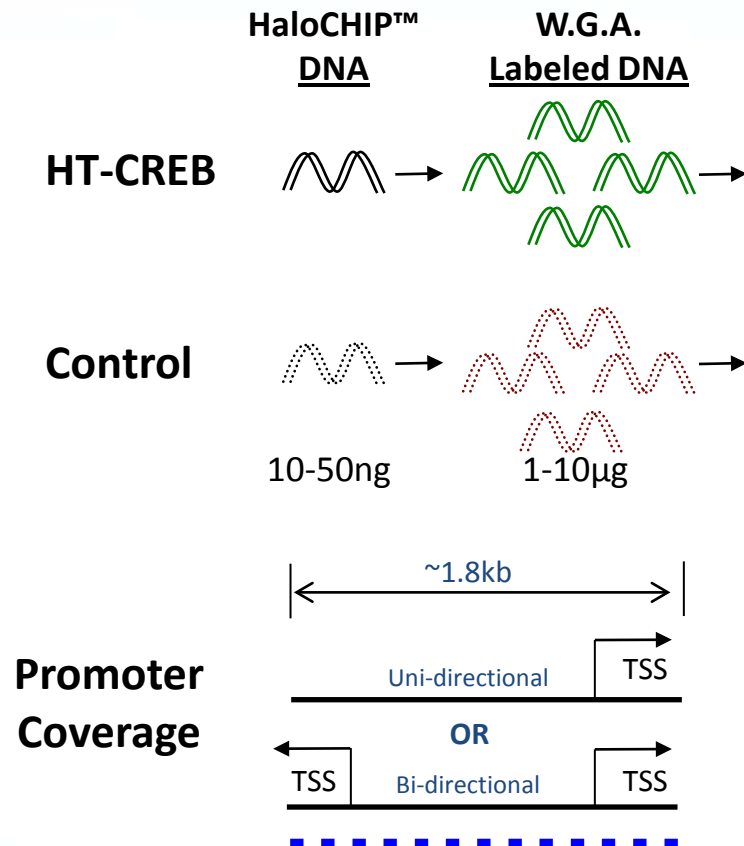
**Sample DNA**

**Background DNA**

# CREB HaloCHIP-chip genome wide analysis



## CREB HaloCHIP™-chip array



27,661 promoters - 385,000 probes

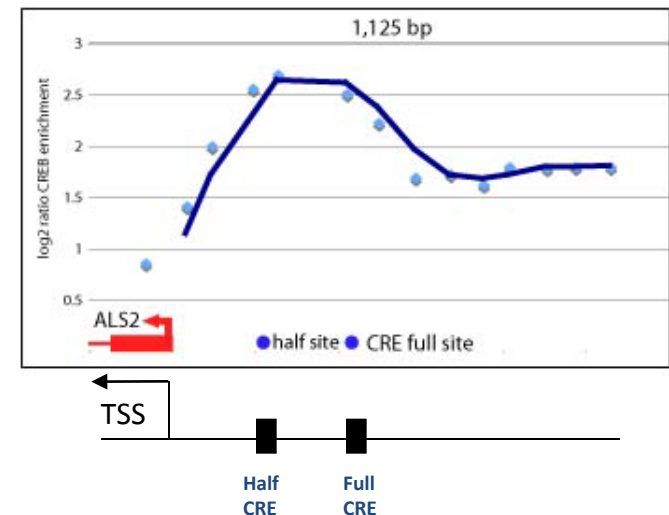
# Gene Ontology (GO) and Promoter Analysis



## CREB HaloCHIP™-ChIP Top 1% Promoters

<u>Cellular Functions</u>	<u># of Promoters</u>	<u>p-value</u>
Histone Assembly	12/65	1.26E-06
Chromatin architecture	20/261	7.63E-07
Ribonucleic Complexes	26/392	7.06E-07
RNA processing	26/395	8.01E-07
DNA metabolism	38/638	2.93E-08
Nucleic acid binding	110/2764	2.19E-09

## Promoter Binding Profile

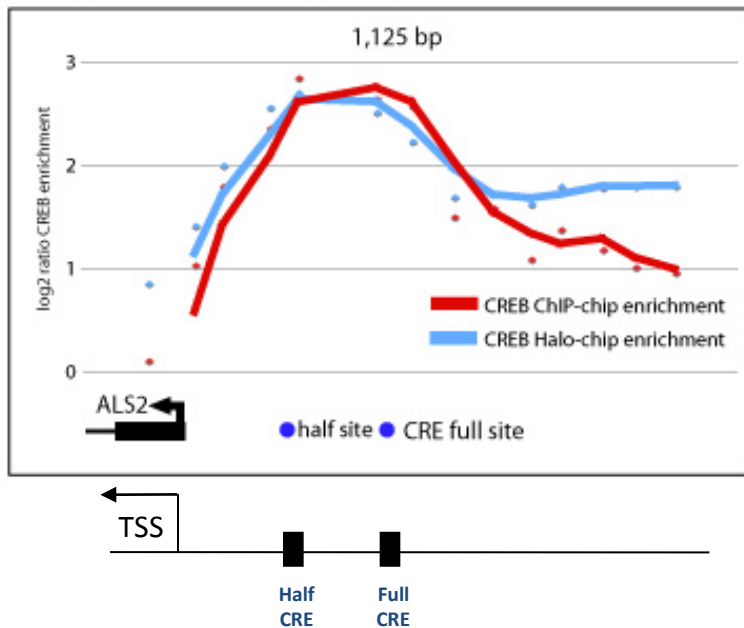


- List of promoters all involved in processes CREB is known to regulate.
- Binding profile shows peaks binding above CRE consensus sites

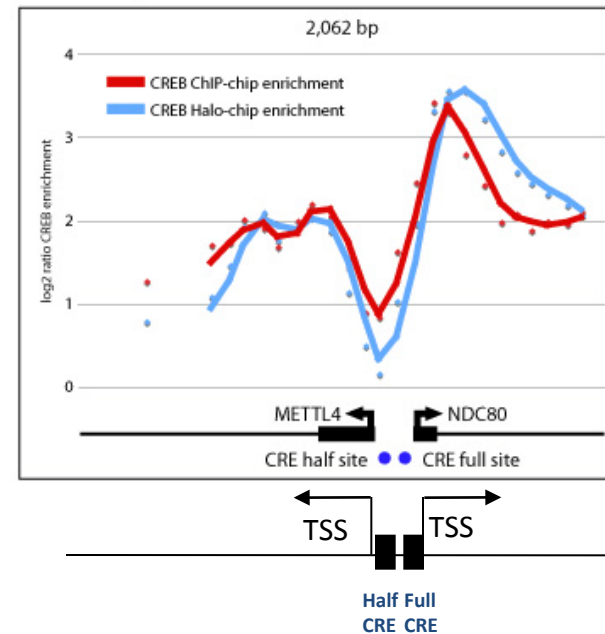
# HaloCHIP™ and ChIP Binding Patterns



## Uni-directional Promoter

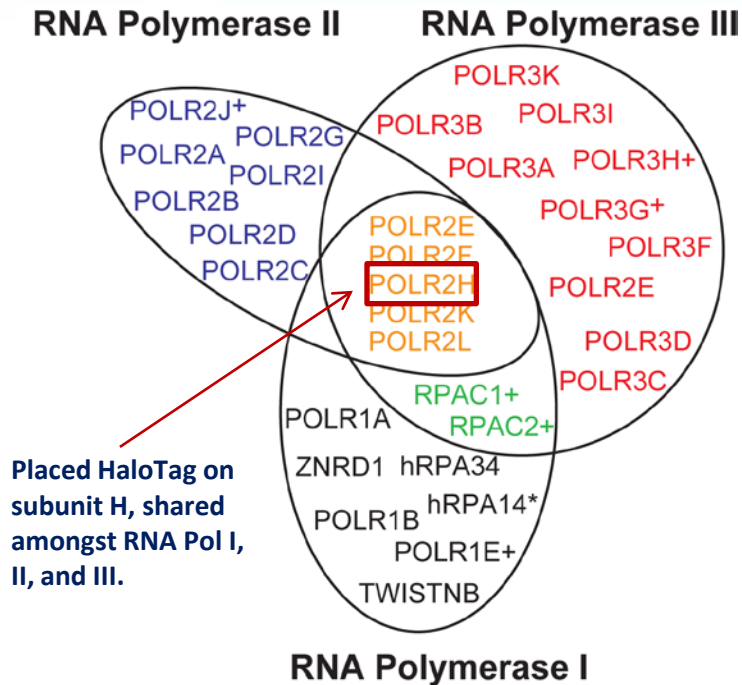


## Bi-directional Promoter

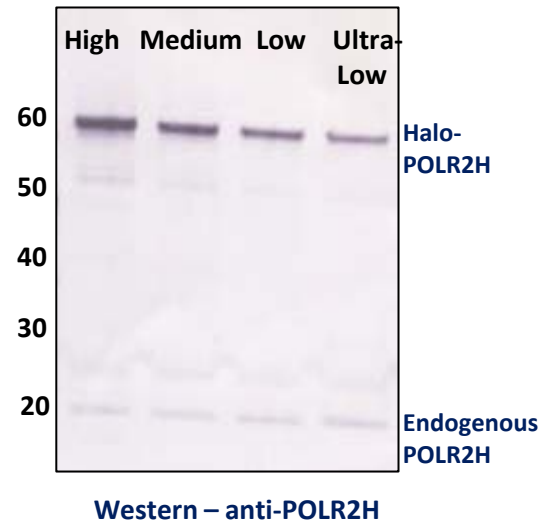


- Overlapping genomic binding patterns between endogenous CREB and Halo-CREB
- High percentage of bi-directional promoters showing downstream CREB binding

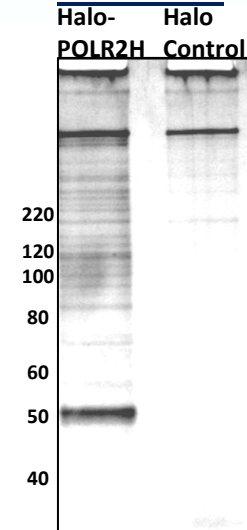
# Expression Studies and Isolation of Eukaryotic RNA Polymerases



Expression levels of Halo-POLR2H  
CMV promoter deletion series



Pull-down



- A CMV deletion series yielded HT-POLR2H expression over a 50-fold range.
- Pull-down performed for each in triplicate and analyzed by MudPIT mass spectrometry.



# Qualitative Data from Triplicate Experiments – RNAP Subunits

	HT Control	HT High	HT Medium	HT Low	HT Ultra
POLR1A		XXX	XXX	XXX	XXX
POLR1B		XXX	XXX	XX	XXX
RPAC1		XXX	XXX	XX	XXX
POLR1D		XXX	XXX	XX	XX
POLR1E		X	XX		X
hRPA34		X	XXX	X	XXX
TWISTNB		X	X	X	
ZNRD1		XX	XX	X	X
POLR2A		XXX	XXX	XXX	XXX
POLR2B		XXX	XXX	XX	XXX
POLR2C		XXX	XXX	XX	XXX
POLR2D		XXX	XXX	XX	XXX
POLR2E		XXX	XXX	XX	XXX
POLR2F		XX			
POLR2G		XXX	XXX	X	XX
POLR2H		XXX	XXX	XXX	XXX
POLR2I		XXX	XX	X	XX
POLR2J		XXX	XXX	XX	XX
POLR2K		X	X		X
POLR2L		XXX	XXX	X	XX
POLR3A		XXX	XXX	XX	XXX
POLR3B		XXX	XXX	X	XXX
POLR3C		XXX	XXX		X
POLR3D		XXX	XXX	XX	XX
POLR3E		XXX	XXX	X	XX
POLR3F		XXX	XXX	X	XXX
POLR3G		XXX	XXX	X	X
POLR3G-LIKE		XX	XX		X
POLR3H		X	XXX		
POLR3I		XX	XXX		X
POLR3K		XXX	XXX	X	XXX

Number of hits out of 3 replicates

	0 of 3
X	1 of 3
XX	2 of 3
XXX	3 of 3

- Excellent subunit recovery across CMV deletion series

Overall recovery out of all possible subunits. ←



# Comparison with FLAG-POLR2H

Acronym	HT-POLR2H	HT Control	FLAG-POLR2H	FLAG Control
POLR1A	XXX		XX	
POLR1B	XXX		X	
RPAC1	XXX		XXX	XX
POLR1D	XXX			
POLR1E	X			
hRPA34	X		XX	
TWISTNB	X			
ZNRD1	XX			
POLR2A	XXX		XXX	X
POLR2B	XXX		XXX	
POLR2C	XXX		XX	
POLR2D	XXX		X	
POLR2E	XXX		X	
POLR2F	XX			
POLR2G	XXX			
POLR2H	XXX		XXX	
POLR2I	XXX			
POLR2J	XXX			
POLR2K	X			
POLR2L	XXX			
POLR3A	XXX		X	
POLR3B	XXX			
POLR3C	XXX		X	
POLR3D	XXX		XX	
POLR3E	XXX			
POLR3F	XXX			
POLR3G	XXX			
POLR3G-LIKE	XX			
POLR3H	X			
POLR3I	XX			
POLR3K	XXX			

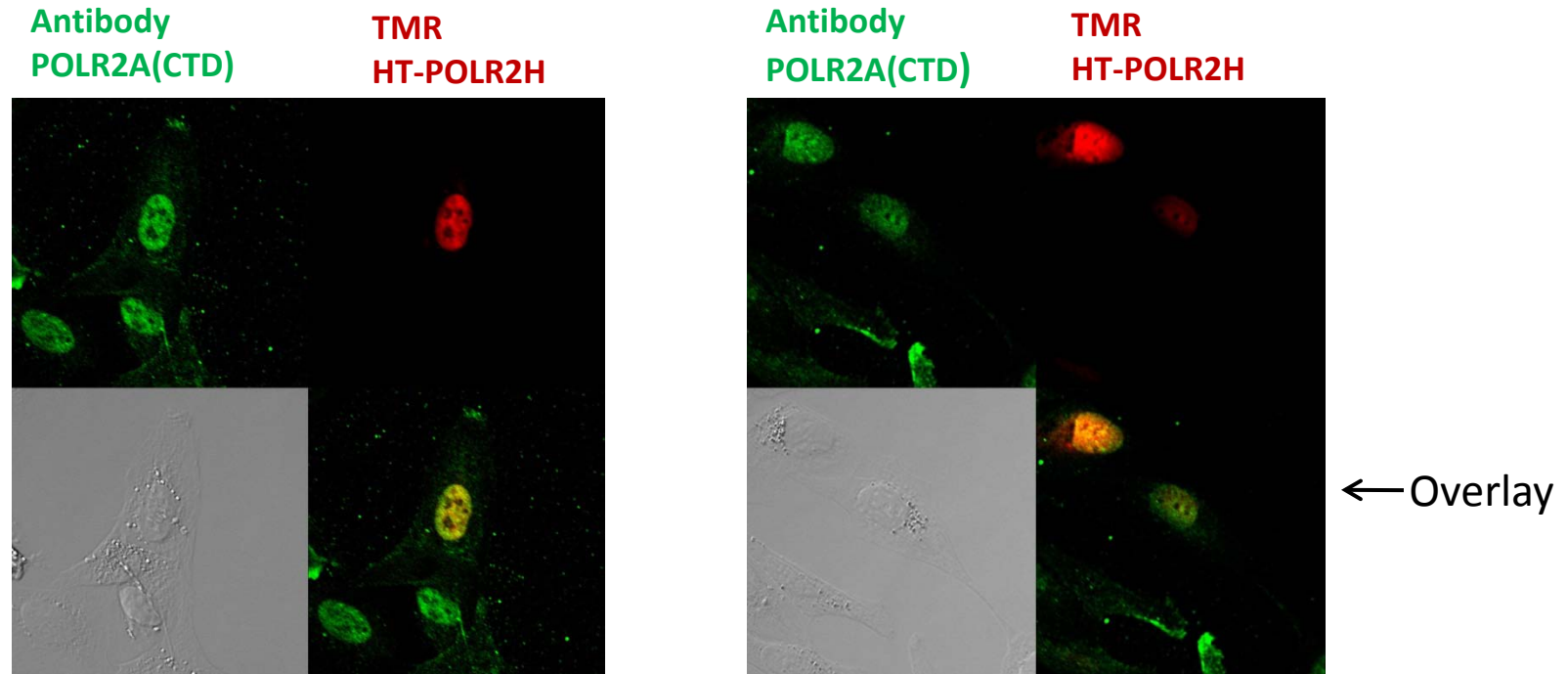
Number of times core RNAP subunits identified by MudPIT analysis in triplicate experiments.

	0 of 3
X	1 of 3
XX	2 of 3
XXX	3 of 3

- Improved recovery with HaloTag
- Higher reproducibility
- Lower background

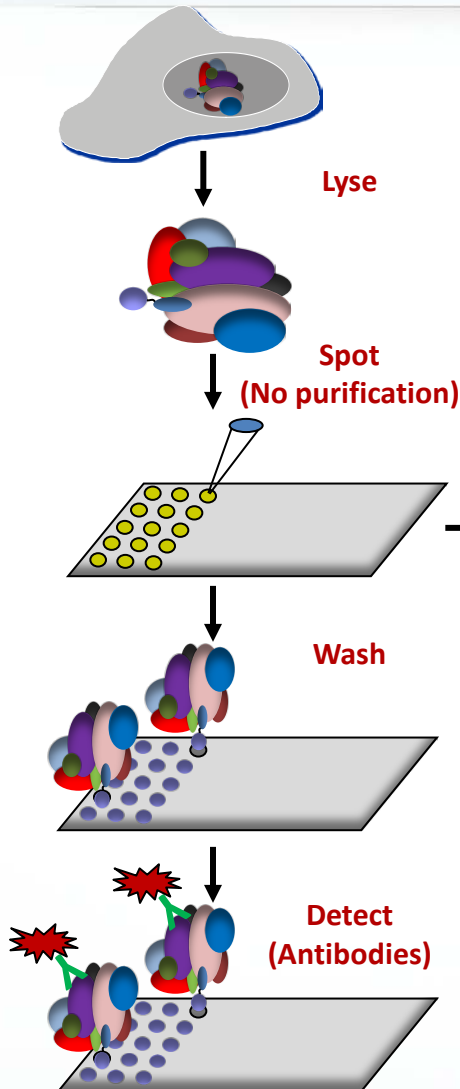
# Imaging of HaloTag<sup>®</sup>-POLR2H

## Co-localization with Endogenous POLR2A

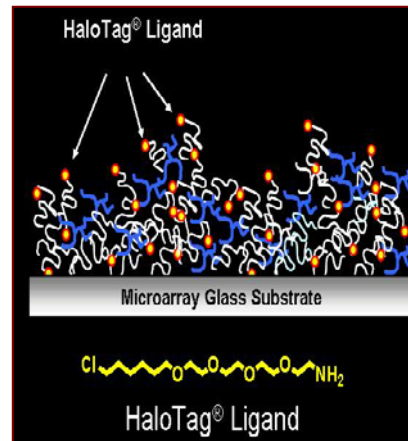


- HaloTag-POLR2H is properly localized to the nucleus.
- Co-localization with antibody specific to the CTD of POLR2A.

# Oriented Capture of HaloTag-POLR2H on HaloLink™ Arrays



HaloLink™ Slide Surface  
(PEG linker with Chloroalkane)



## HaloTag-POLR2H Capture on Slide

### Antibodies

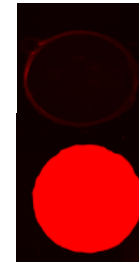
POLR2A  
(CTD)

HaloTag

Experimental  
(Halo-POLR2H)



Control  
(HaloTag)

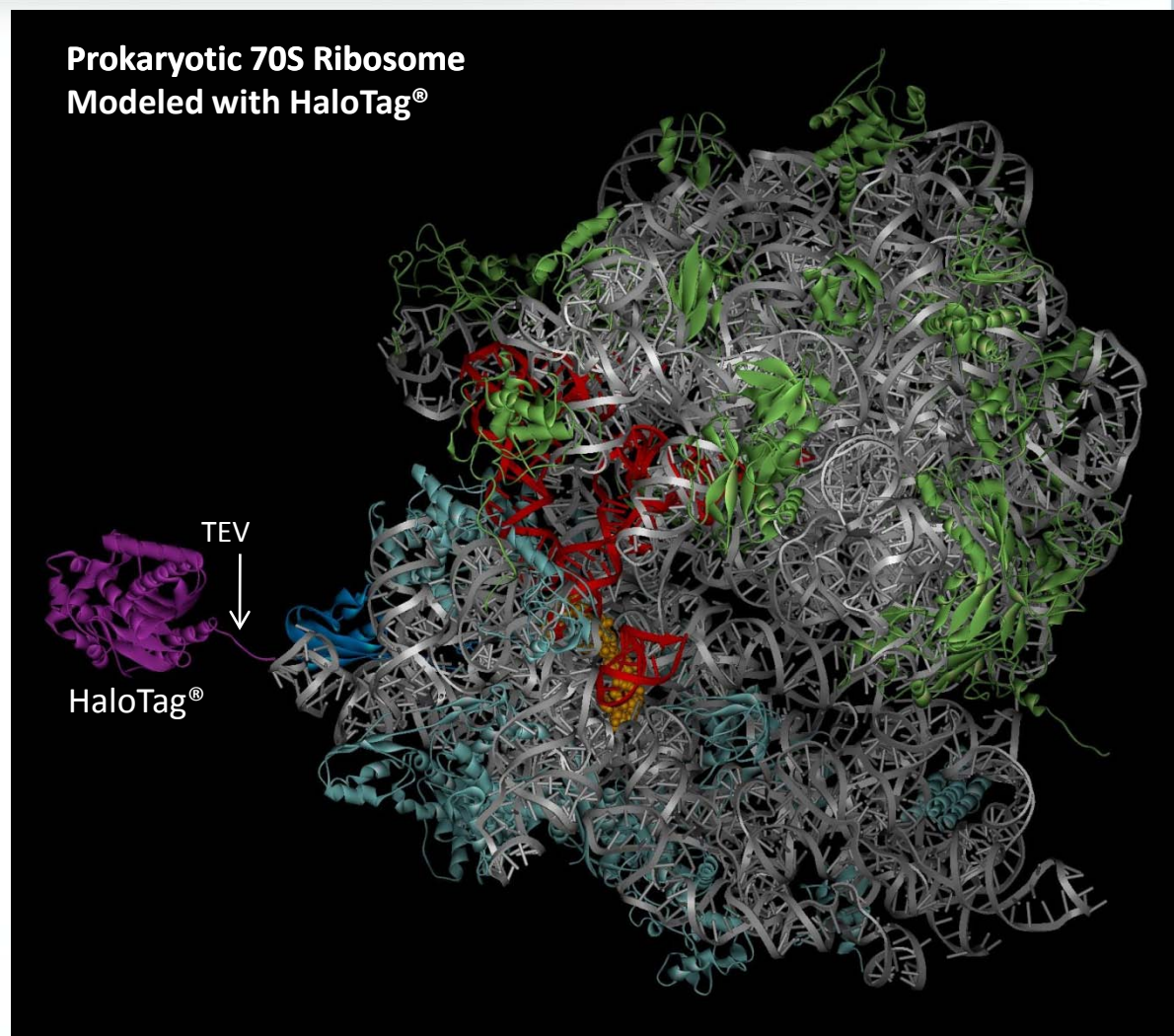


❖ HaloTag-POLR2H isolated on slide directly from lysate.

❖ Detection of binding partner suggests RNA Polymerase complex capture on slides.

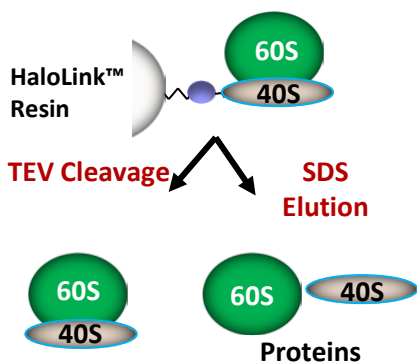
## Isolation of the Ribosome

- One of the largest macromolecular machines.
- Highly abundant.
- Many interacting partners, mRNAs.
- Difficult to capture using a single protein fusion tag.
- Placed HaloTag<sup>®</sup> on 40S subunit protein, Human RPS9.



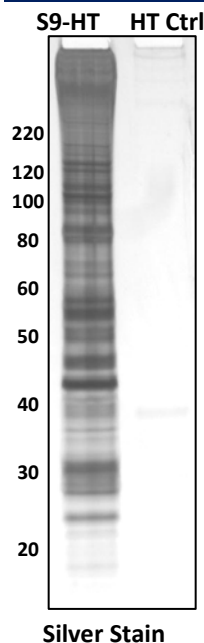
# Capture of Eukaryotic Ribosomes Using S9-HT

## S9-HT Pull-down



- HaloTag® remains on resin
- Rapid binding -15min.
- No diffusion off resin
- Capture low levels of protein

## TEV Cleavage

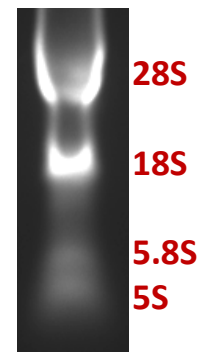


## Identified by MS (LC-MS/MS)

- 31 of 33 40S proteins
- 42 of 50 60S proteins
- 2 Poly-A binding proteins
- 1 GNF exchange protein
- 9 Nuclear ribonucleoproteins
- 2 Initiation factors
- 2 Elongation factors
- 2 Splicing factors

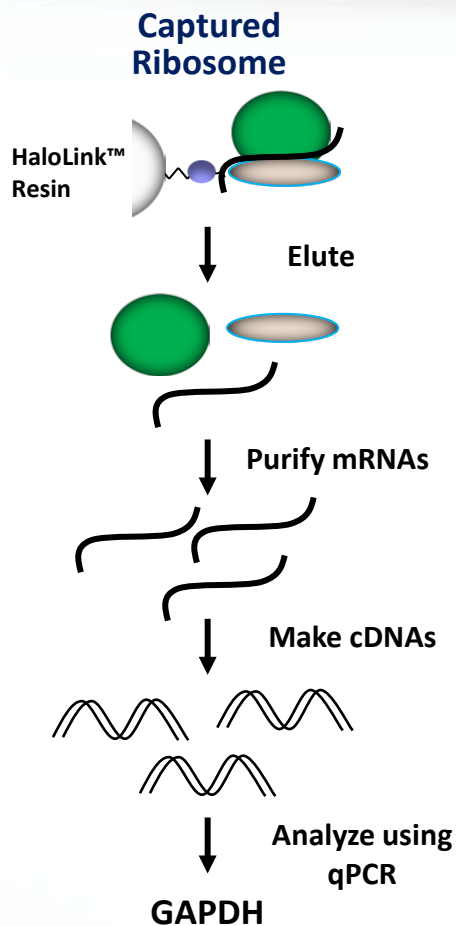
## RNA Capture

Purified RNAs

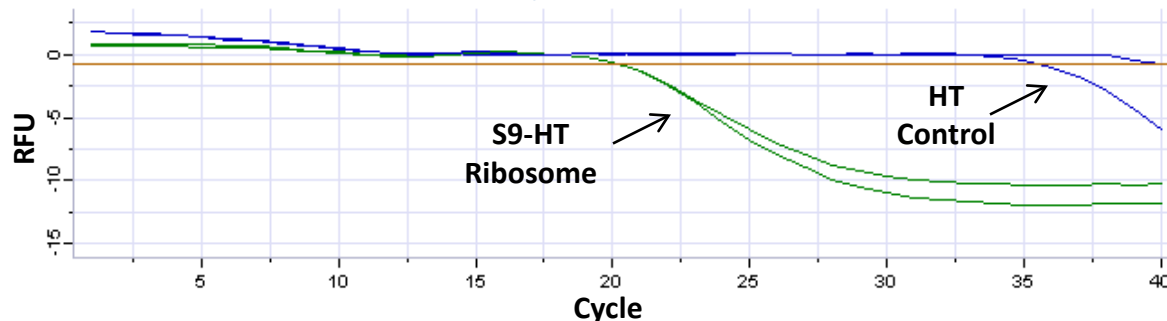


- Purified the 80S ribosome with single step from HeLa cells.
- Efficiently cleave complex from resin with TEV protease.

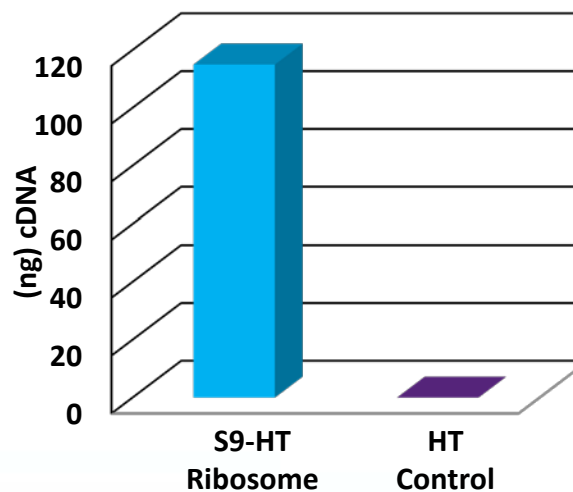
# Isolated Ribosomes are Bound to mRNAs



## Plexor qPCR GAPDH cRNA Amplification



## Total GAPDH cDNA

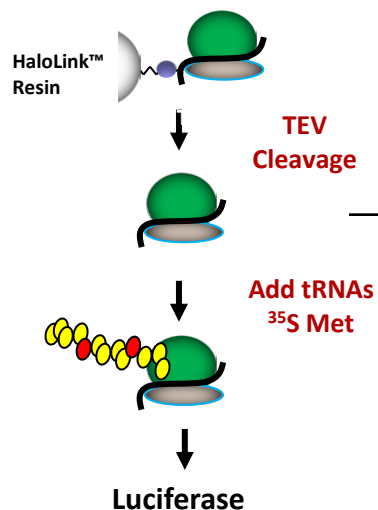


- GAPDH mRNA bound to isolated ribosomes
- Potential for complete mRNA analysis.

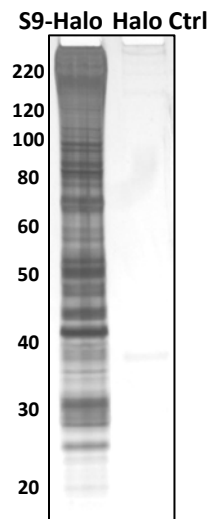


# Isolated Ribosomes are Active for Translation

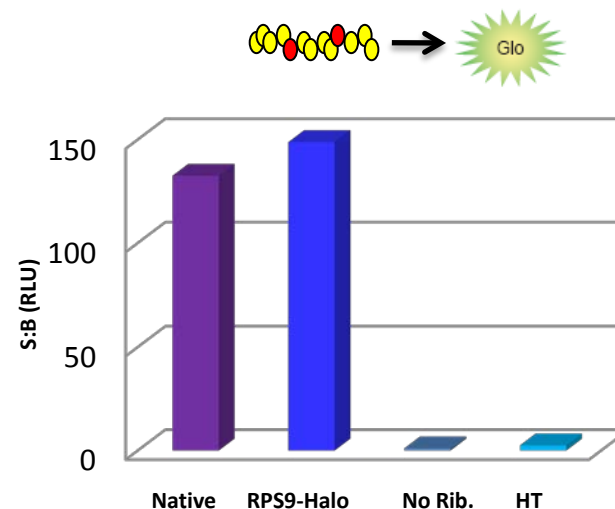
## Pulldown of Luc-mRNA RPS9-HT in Luc-HEK293



## Ribosome Pulldown



## Specific luciferase synthesis



- Ribosomes which have incorporated RPS9-Halo are active for elongation.
- Bound to Luc-mRNA and can translate functional luciferase protein.



# Monitoring Ribosomal Trafficking and Populations

**Serum Starve**  
**RPS9-HT**

**Pulse**  
**TMR Label**

**Serum Recovery**

**Chase**  
**Green Label**

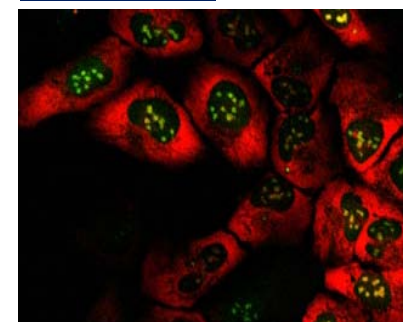
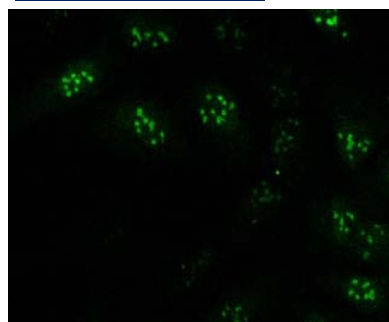
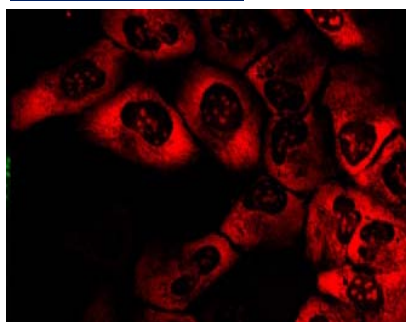
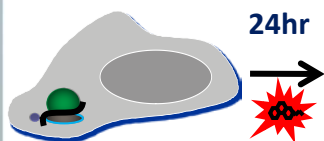
**Wash**

**Image Overlay**

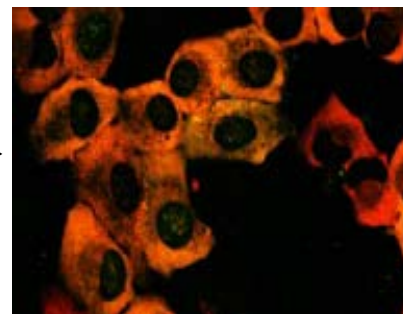
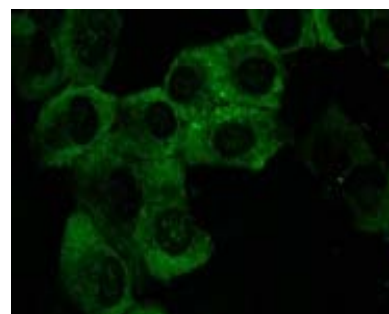
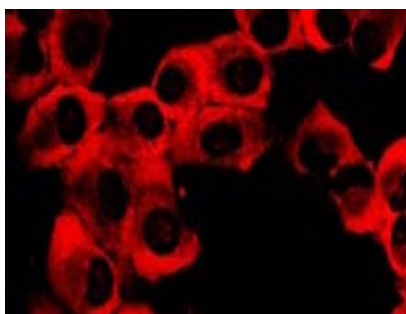
**Old population**

**New population**

**Old and new**



OR



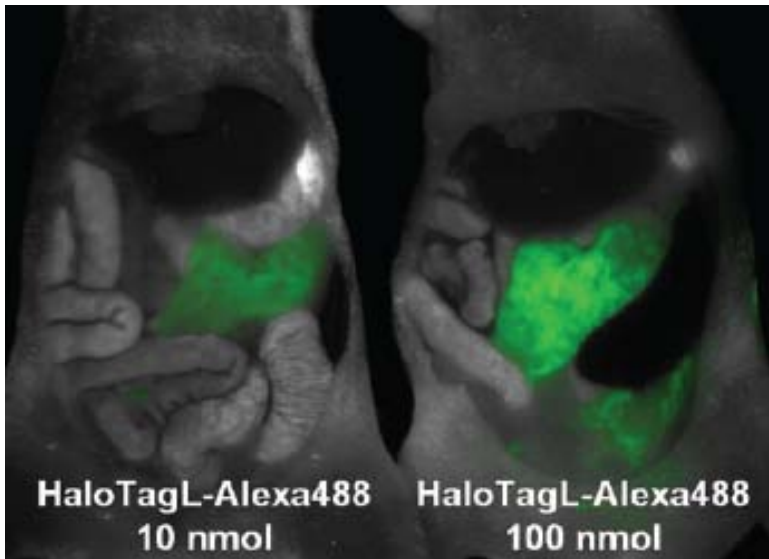
- Newly synthesized ribosomal proteins localized to nucleoli 3hrs post-stress.
- After 24hrs of recovery, ribosomal populations re-localized to cytoplasm.

# HaloTag<sup>®</sup> Platform



- In vivo fluorescent imaging
- Future directions

# *In Vivo* Imaging

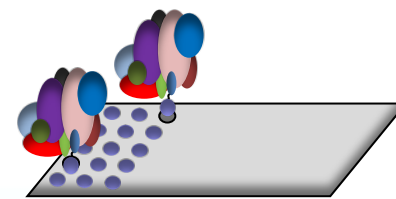
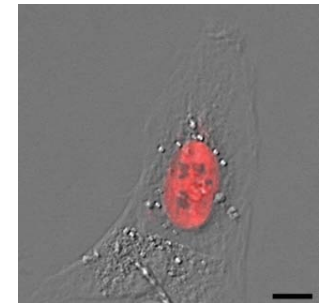
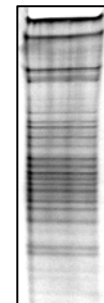
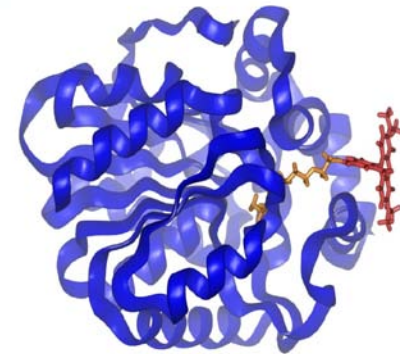


Fluorescent *in vivo* imaging using HaloTag<sup>®</sup> technology allows for development of imaging ligands

- PET
- near IR

# Summary

- Evolved HaloTag<sup>®</sup> protein for specific, covalent, and rapid binding.
- HaloTag<sup>®</sup> technology shows strength and advantages for a variety of mammalian applications:
  - *Protein purification*
  - *Capture on surfaces*
  - *Protein:DNA interactions*
  - *Protein complex isolation*
  - *Cellular and in vivo imaging*



Glass Slides