

# ***In Vitro and In Vivo Methods to Study Protein:Protein Interactions***

Rob Brazas, Ph.D.

November, 2011



# ***Presentation Overview***

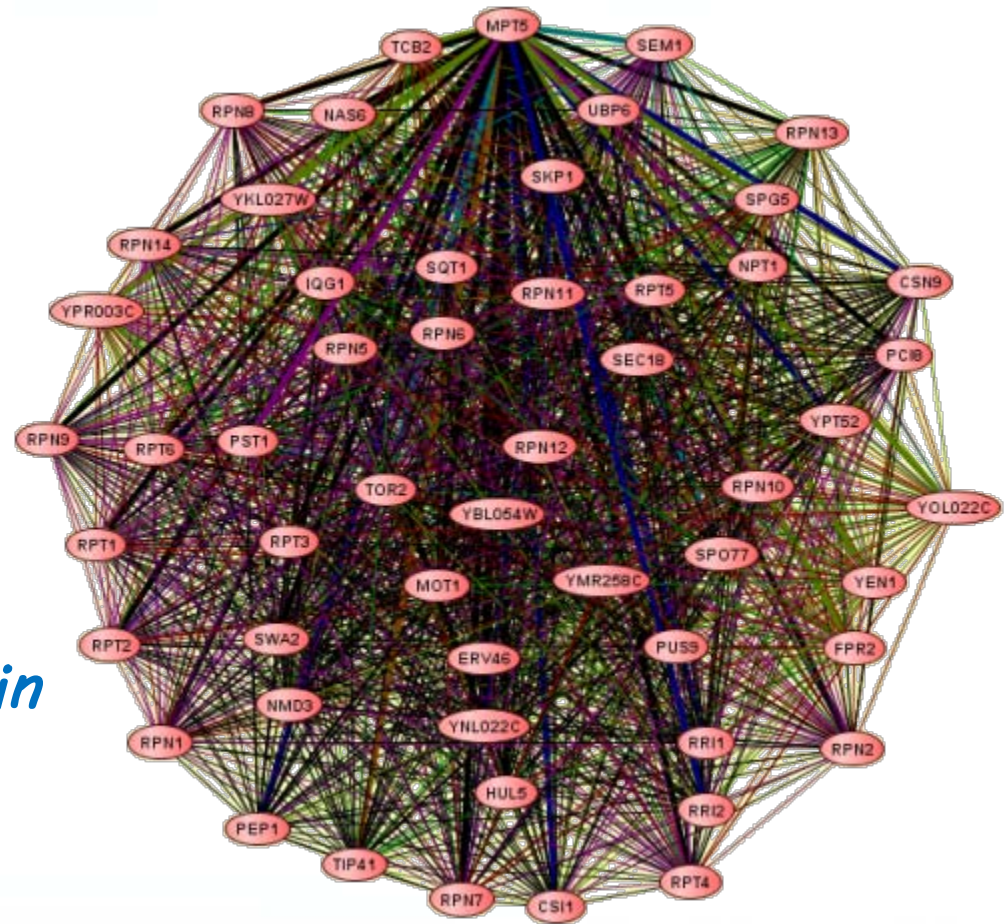
- Review of Methods to Study Protein:Protein Interactions
- In Vitro (Cell-Free) Protein Expression and Application to Interaction Studies
- HaloTag® Fusion Protein Technology and Protein Interaction Experiments
- Applying These Technologies to the Study of Protein-Protein Interactions
  - Case Studies Illustrating Both Discovery and Verification Applications
- Mammalian Two Hybrid Assays
  - Adapting the Yeast Two Hybrid Assay to Mammalian Cells
  - Case Study: Cdk3 Interaction with ATF1
- Other Applications of In Vitro (Cell-Free) Expression and HaloTag® Technology
- Summary

# Protein-Protein Interactions are Critical to All Cellular Processes



- Replication
- Transcription
- Translation
- Signal transduction & more

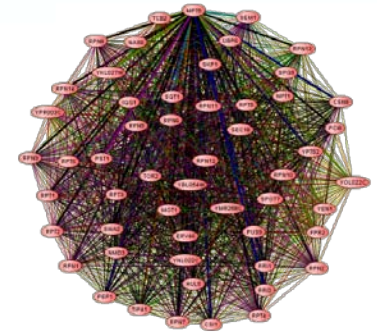
*How are protein:protein interactions studied?*



# ***Common Methods Used to Study Protein:Protein Interactions***



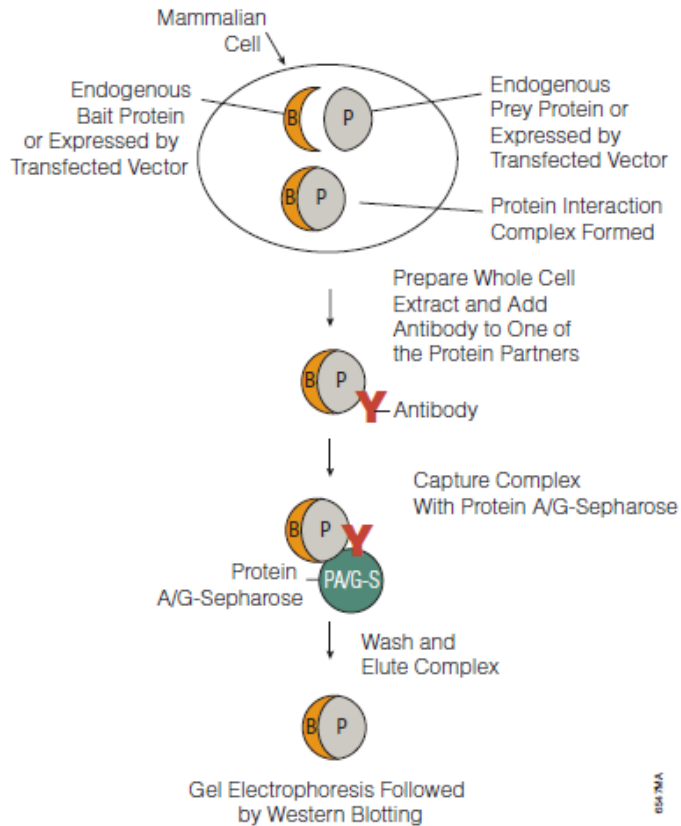
- Co-immunoprecipitation
- Protein Affinity Purification
- Far Western Blotting
- Two Hybrid Assays (yeast, mammalian)
- In Vivo Förster Resonance Energy Transfer (FRET)
- and more...



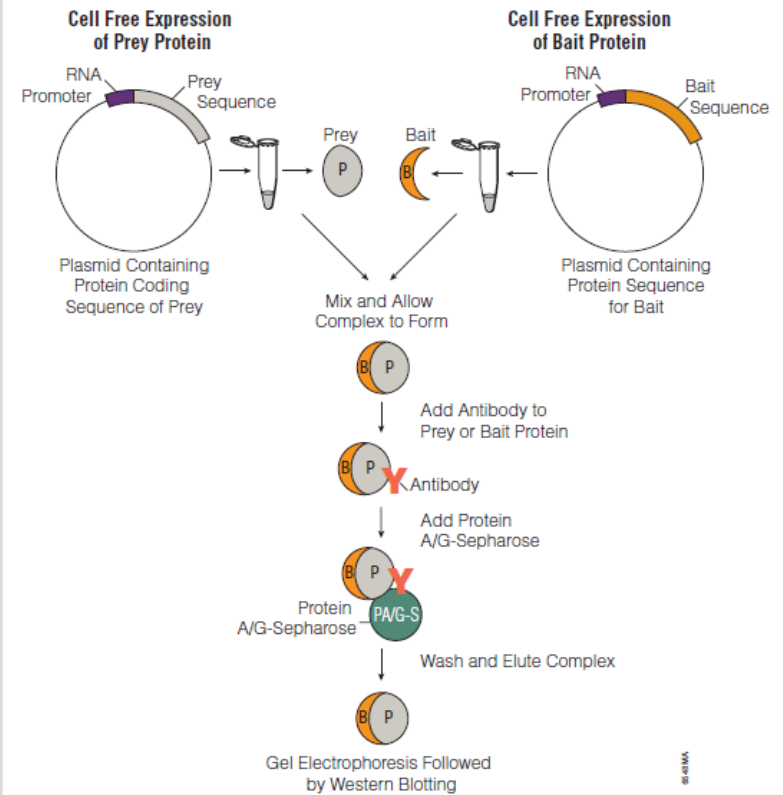
# Co-Immunoprecipitation (Co-IP)

## A Classic Method to Study Interactions

### In Vivo



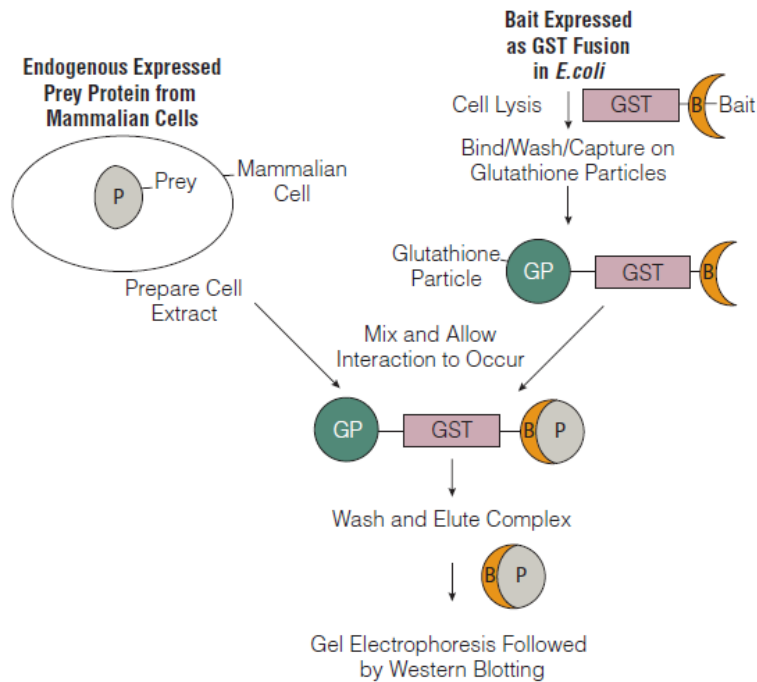
### In Vitro



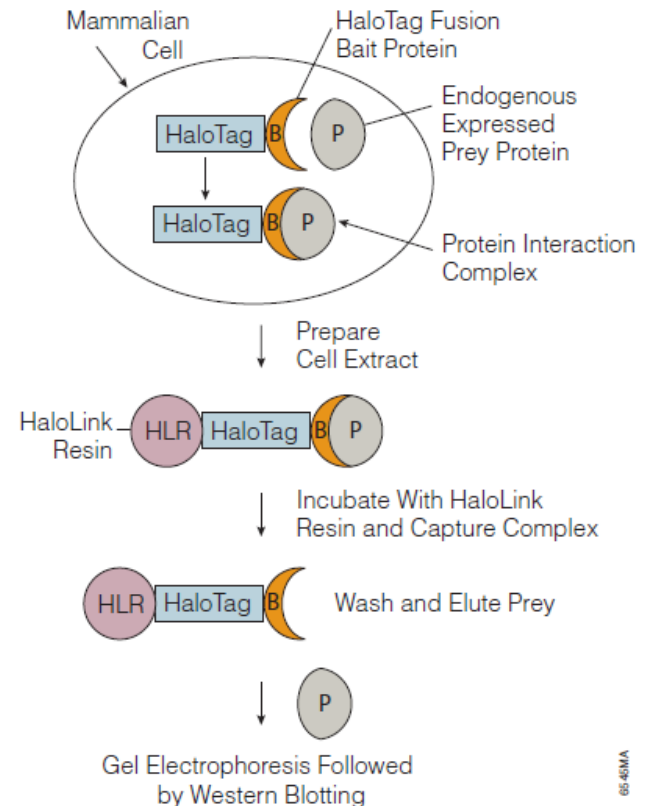
# Protein Affinity Purification Methods

## Identification of Novel Interacting Partners

### Purified Bait + Cell Extracts



### In Vivo with Tagged Bait

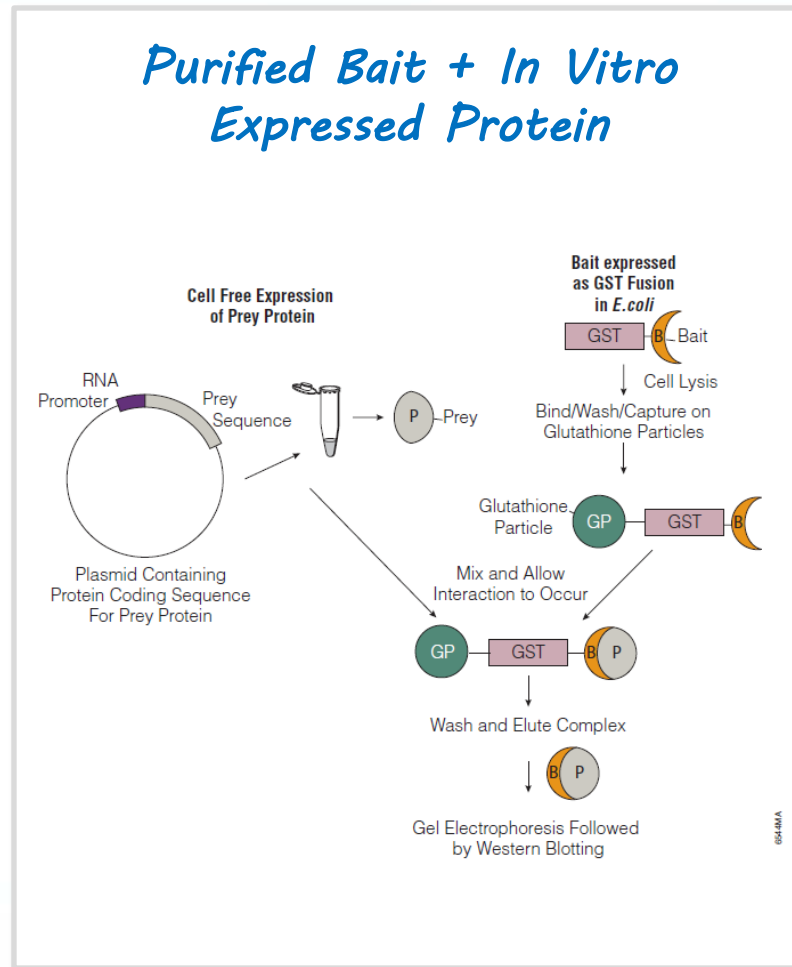


# Protein Affinity Purification Methods

## A Quick Method for Verifying an Interaction

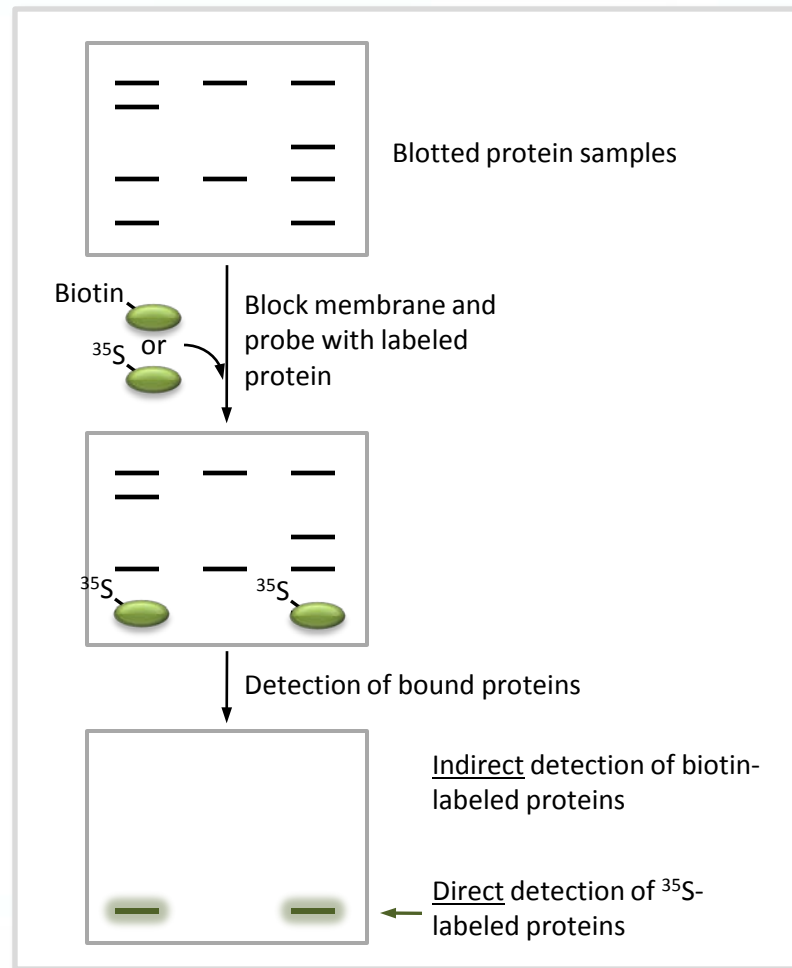


### Purified Bait + In Vitro Expressed Protein



# Far Western Blots

## A Rapid Method for Testing Interactions

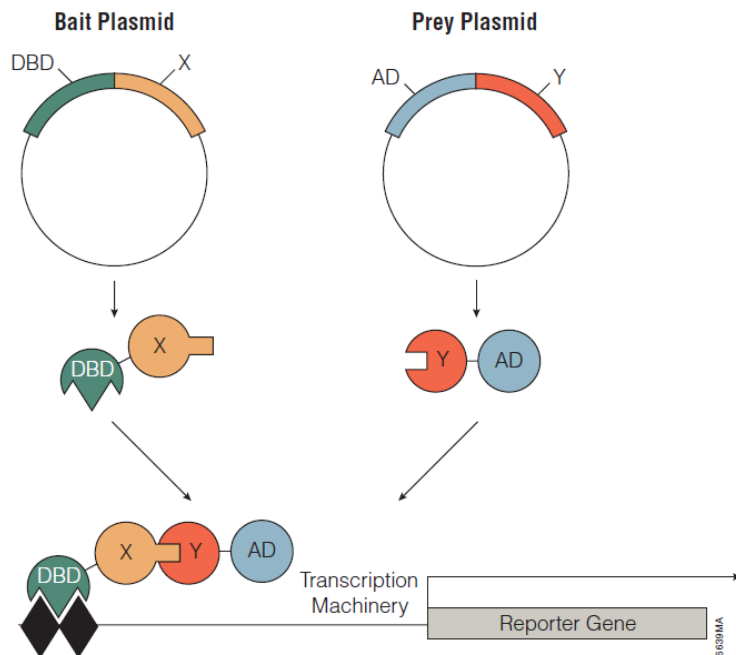




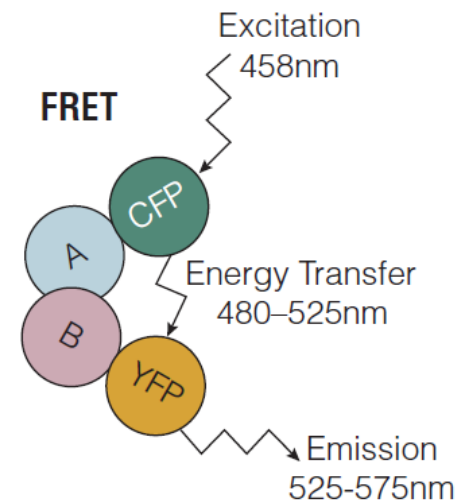
# Two Hybrid and FRET Assays

## Newer Protein:Protein Interaction Assays

### Two Hybrid Assays



### Förster Resonance Energy Transfer (FRET)

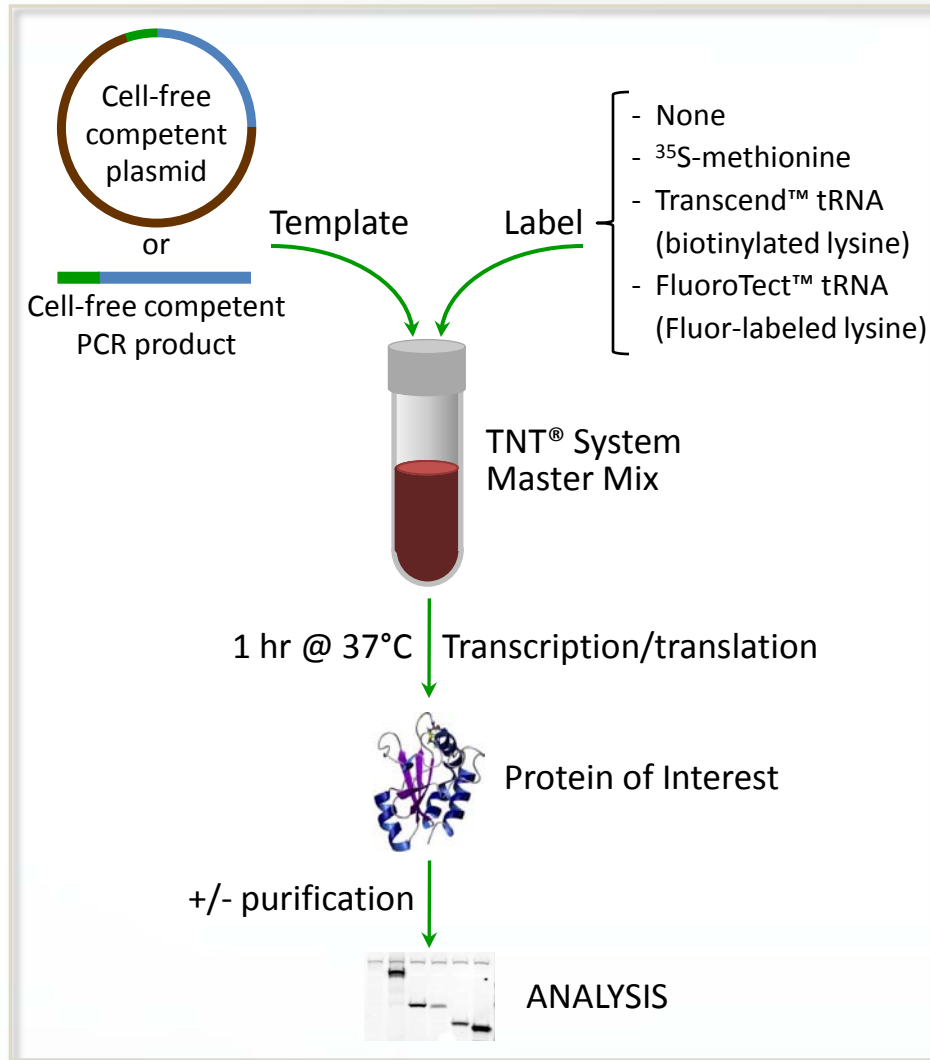


# ***In Vitro (Cell-Free) Expression – the Rapid, Easy-to-Use Solution***

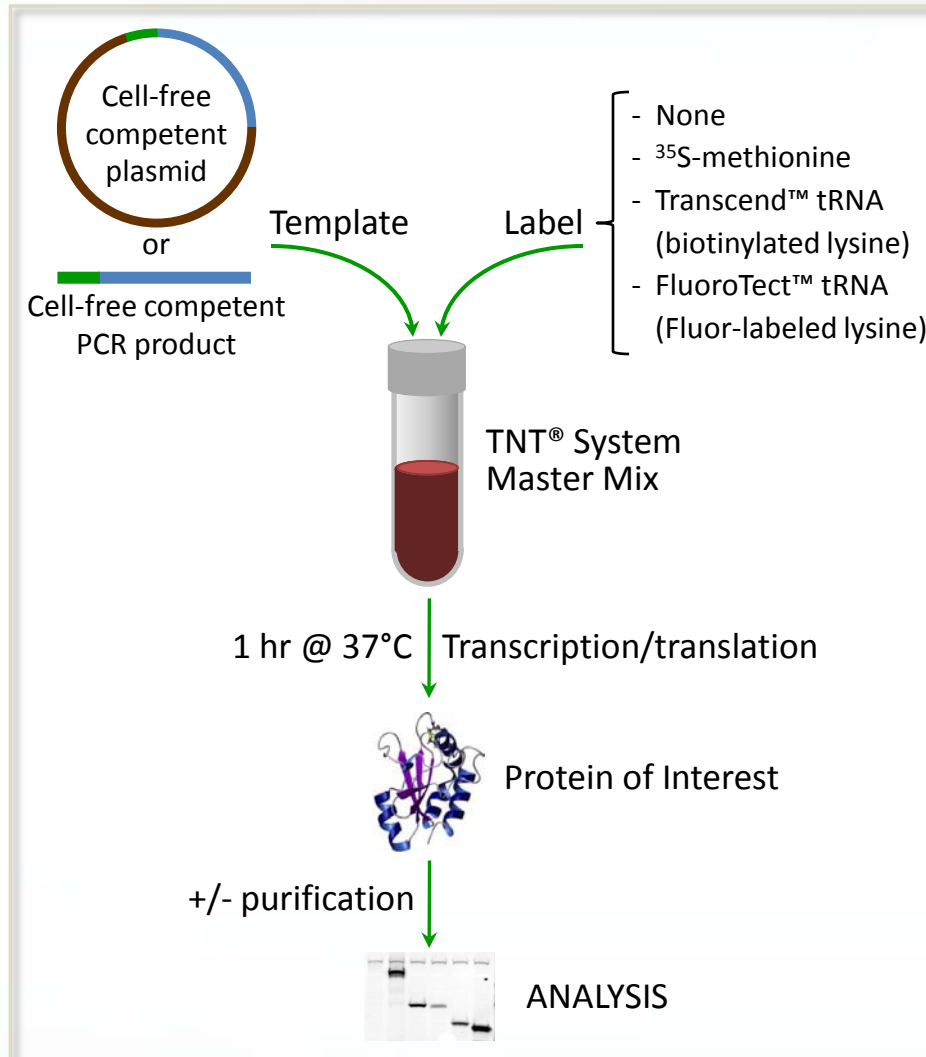
In Vitro Production of Target Proteins Using Cellular Extracts for Protein:Protein Interaction Experiments and More



# Easiest/Fastest Method to Go from DNA to Protein



# Easiest/Fastest Method to Go from DNA to Protein



## • Saves valuable time

- Produce protein in 1-2 hours vs. days to weeks in *E. coli* or mammalian cells
- Use PCR product templates and bypass cloning ORF into expression vector

## • Produces sufficient protein for many applications including:

- Protein-protein interactions
- Co-immunoprecipitations
- Gel-shift assays
- Enzymatic assays

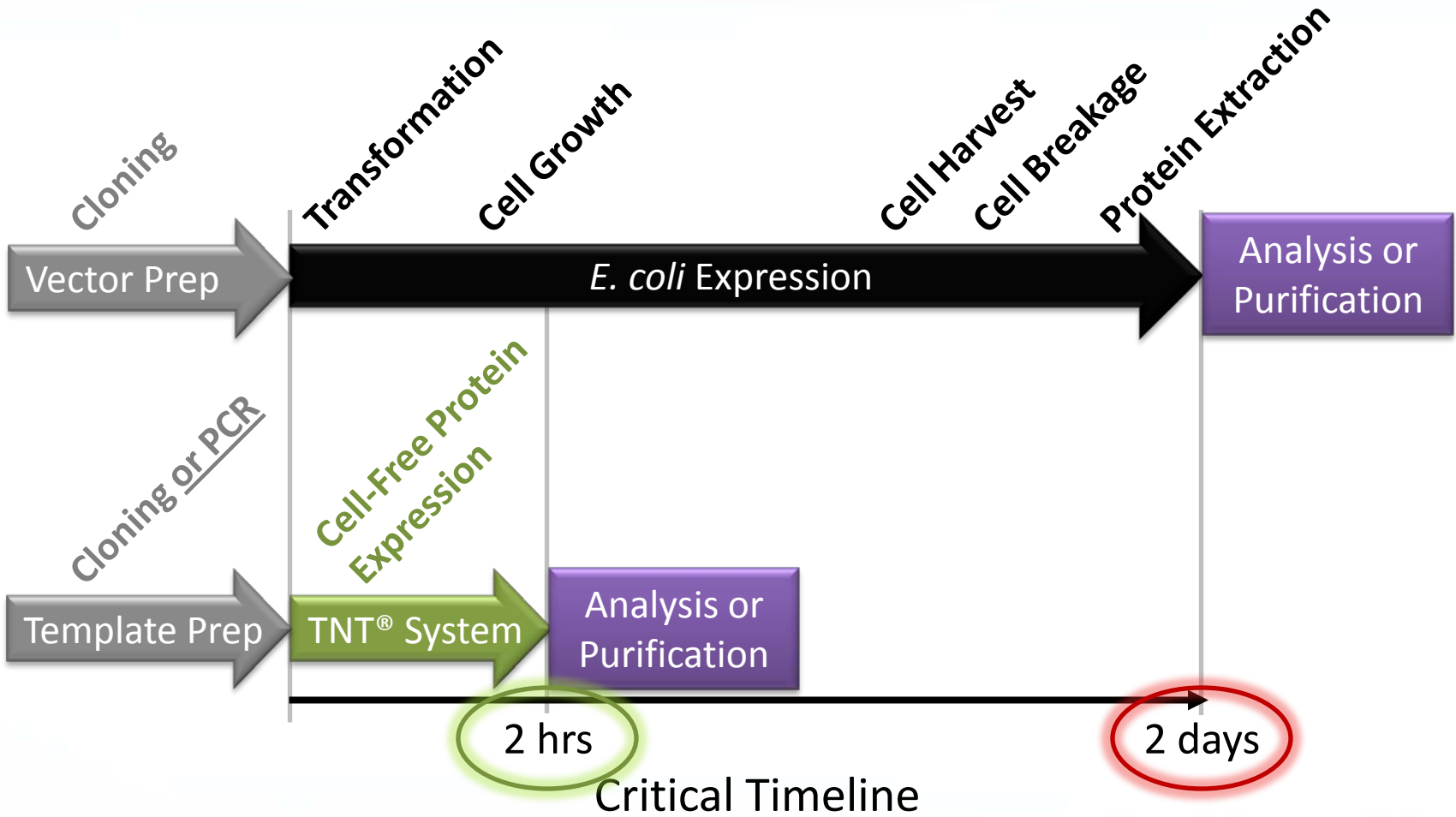
## • Enables analysis of difficult proteins

- Normally toxic to cells
- Insoluble in *E. coli*

## • Simplified detection

- Directly label protein during synthesis
  - Fluorophore, <sup>35</sup>S, biotin

# Faster Protein Production than *E. coli* Systems



# Choices to Match Your Research Needs



>75 citations  
in 2011  
Jan-July

Prokaryotic

*Bacteria*



*E. coli*

Eukaryotic

Plant



Wheat germ

Insect



*Spodoptera frugiperda*

Mammalian



Rabbit reticulocyte

S30 T7 High-Yield  
Protein Expression  
System

Highest Yield

TNT® SP6 High-Yield  
Wheat Germ  
System

Maximal  
soluble protein

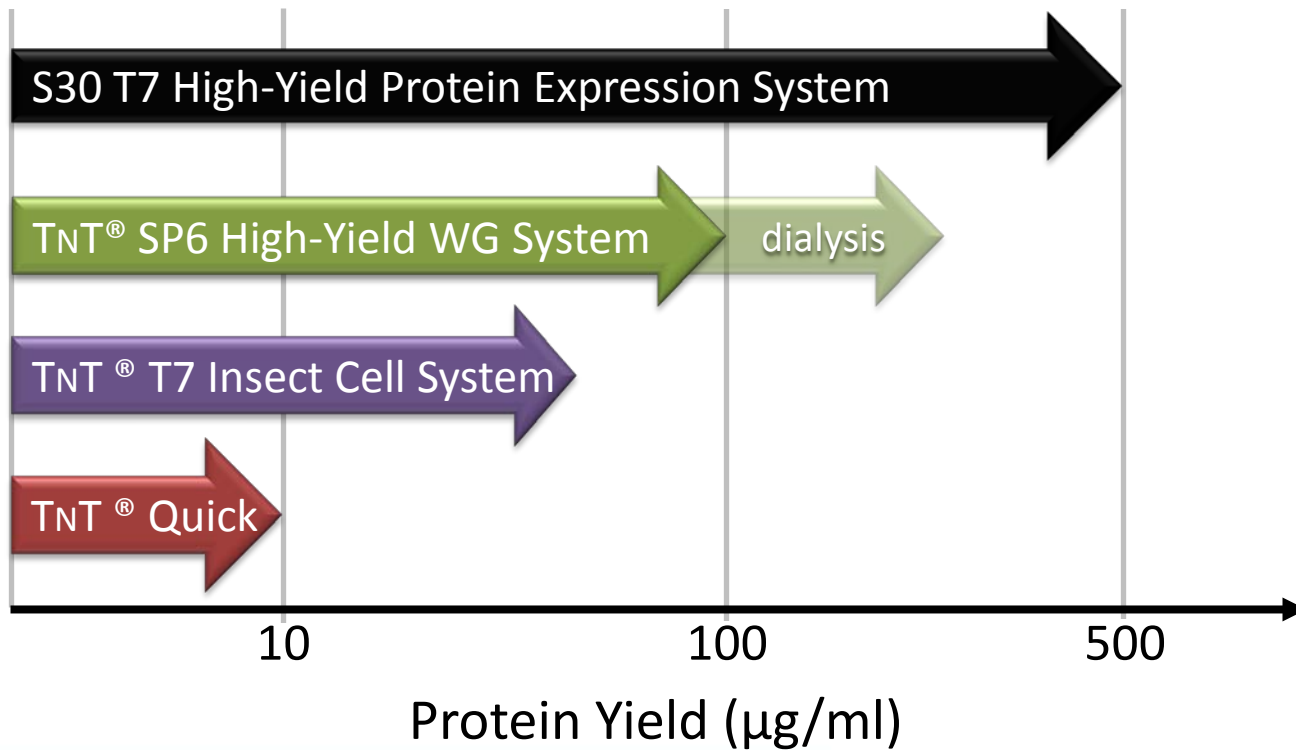
TNT® T7 Insect Cell  
Extract System

Most active  
protein

TNT® Quick Coupled  
T7 and SP6 Systems

Native mammalian  
system

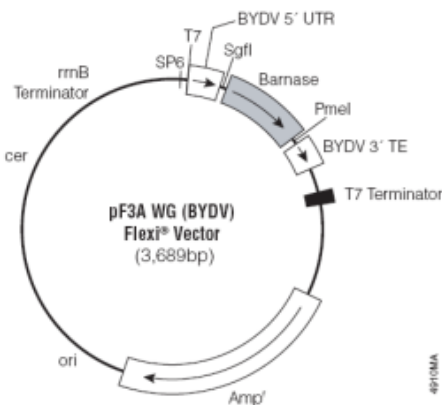
# Maximal Yields from Each System



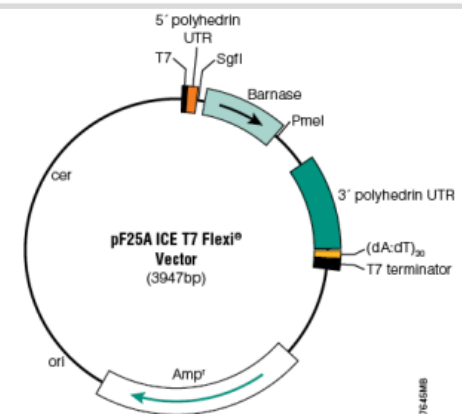
# Characteristics of In Vitro Expression Systems

System	Time	Yield	Requirements
TNT® T7 or SP6 Quick Coupled System	1 hour	≤0.5µg/50µl	<b>Any vector</b> containing a T7 or SP6 promoter upstream of coding sequence.
S30 High Yield (Bacterial)	1 hour	≤25µg/50µl	T7 promoter-driven bacterial expression vector. Can also use very active bacterial promoters (T5)
TNT® SP6 High-Yield Wheat Germ System	2 hours	≤5-12.5µg/50µl	Highest yield with specialized vector containing plant viral sequences (≤5.0µg/rxn). Greatest yield using dialysis method (≤12.5µg/rxn).
TNT® T7 Insect Cell System	4 hours	≤4.0µg/50µl	Require use of a baculovirus expression vector with T7 promoter upstream.

*Optimized Wheat Germ Vector*

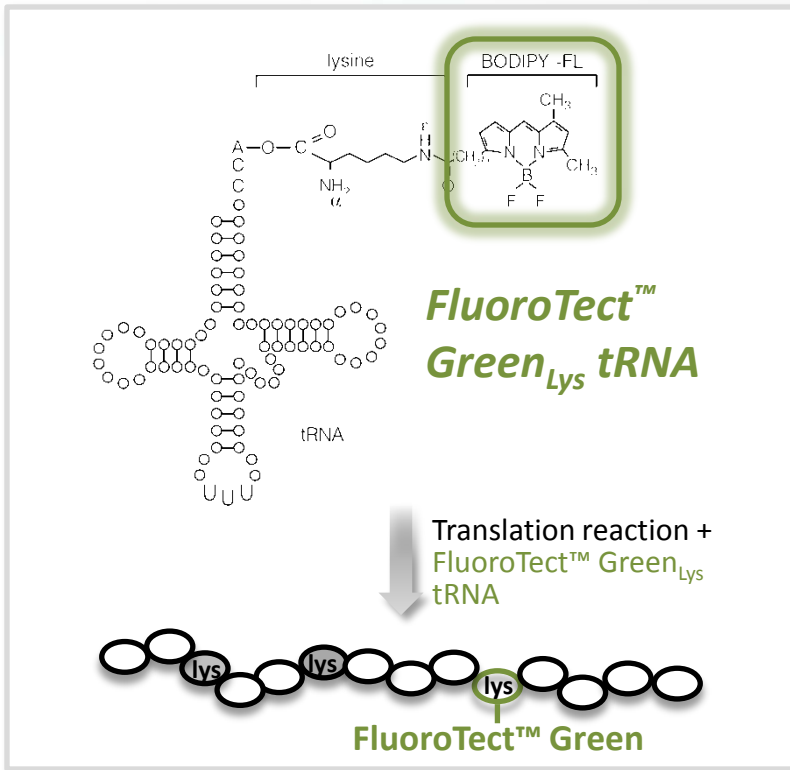


*Insect Cell Vector Example*

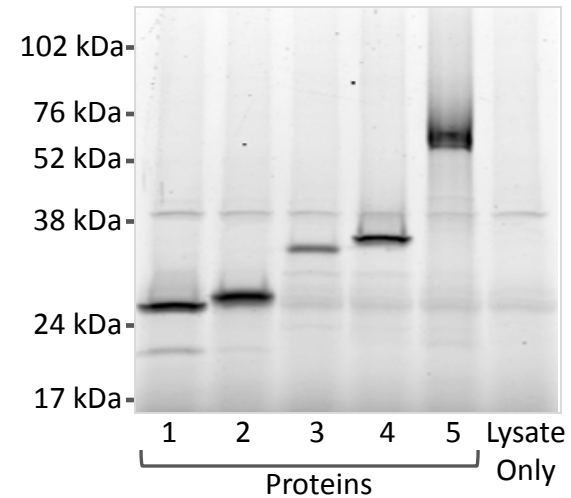




# Fluorescent Detection of In Vitro Expressed Proteins Non-Radioactive Co-translational Labeling



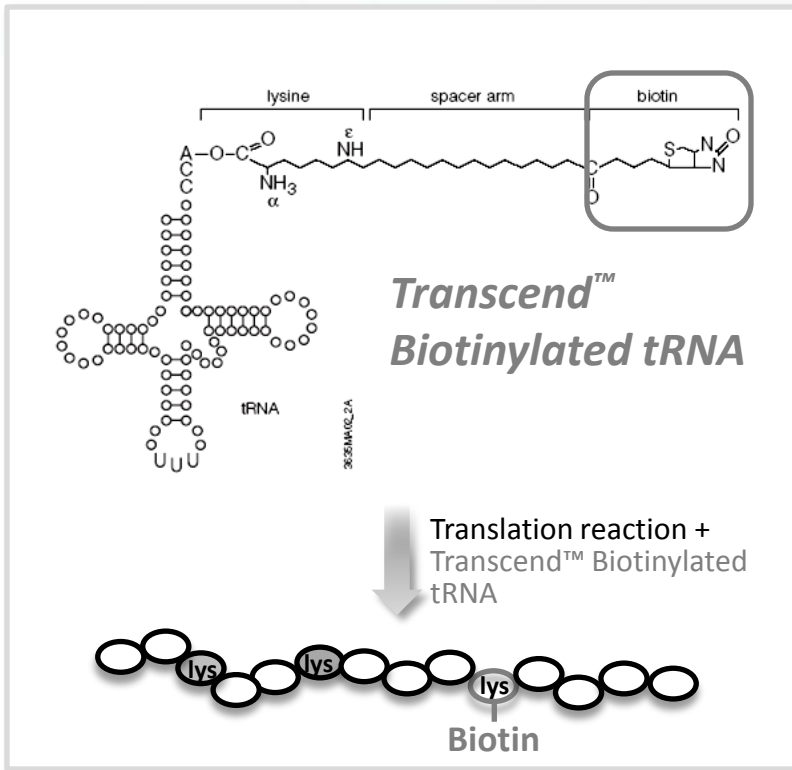
## TNT® T7 Quick Coupled System



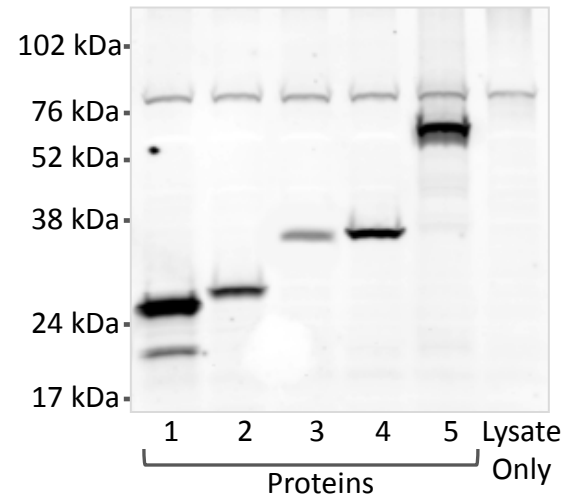
- ✓ Produce active/detectable proteins without radioactivity
- ✓ Direct detection of fluorescently labeled proteins in gels
- ✓ Use in many applications including pulldowns, co-immunoprecipitations, mobility shift assays...

# Indirect Detection of In Vitro Expressed Proteins

## Transcend™ Biotin Co-translational Labeling



### TNT® SP6 Wheat Germ High Yield System



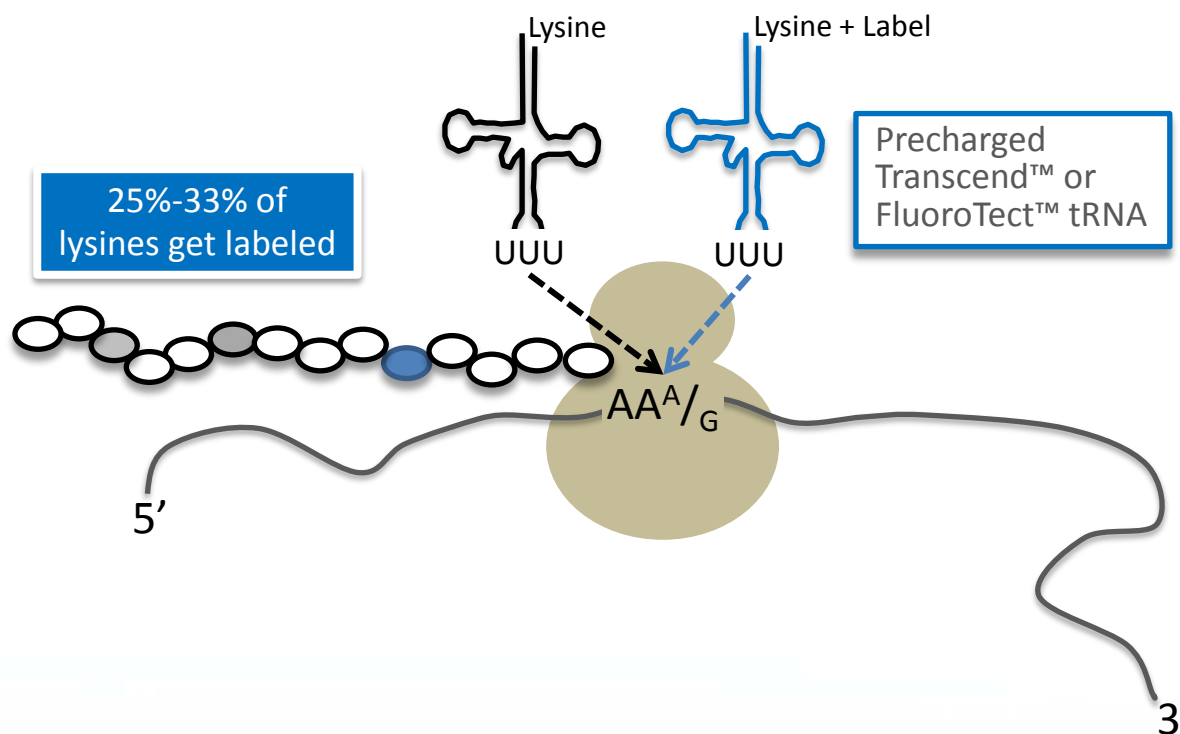
- ✓ Produce active/detectable proteins without radioactivity
- ✓ Indirect detection using streptavidin conjugates (HRP, AlkPhos) of biotin labeled proteins
- ✓ Use in many applications including pull downs, co-immunoprecipitations, mobility shift assays...

# What if There is a Lysine in the Active Site

## No Problem – Only 25-33% of Lysines are Labeled



*FluoroTect™ or Transcend™ tRNA compete with natural lysyl tRNA for incorporation into growing peptide chain*



# ***HaloTag<sup>®</sup> Fusion Protein***

A Unique, Multifunctional Fusion Tag Well-Suited to Protein:Protein Interaction Studies



# What is HaloTag® Technology?

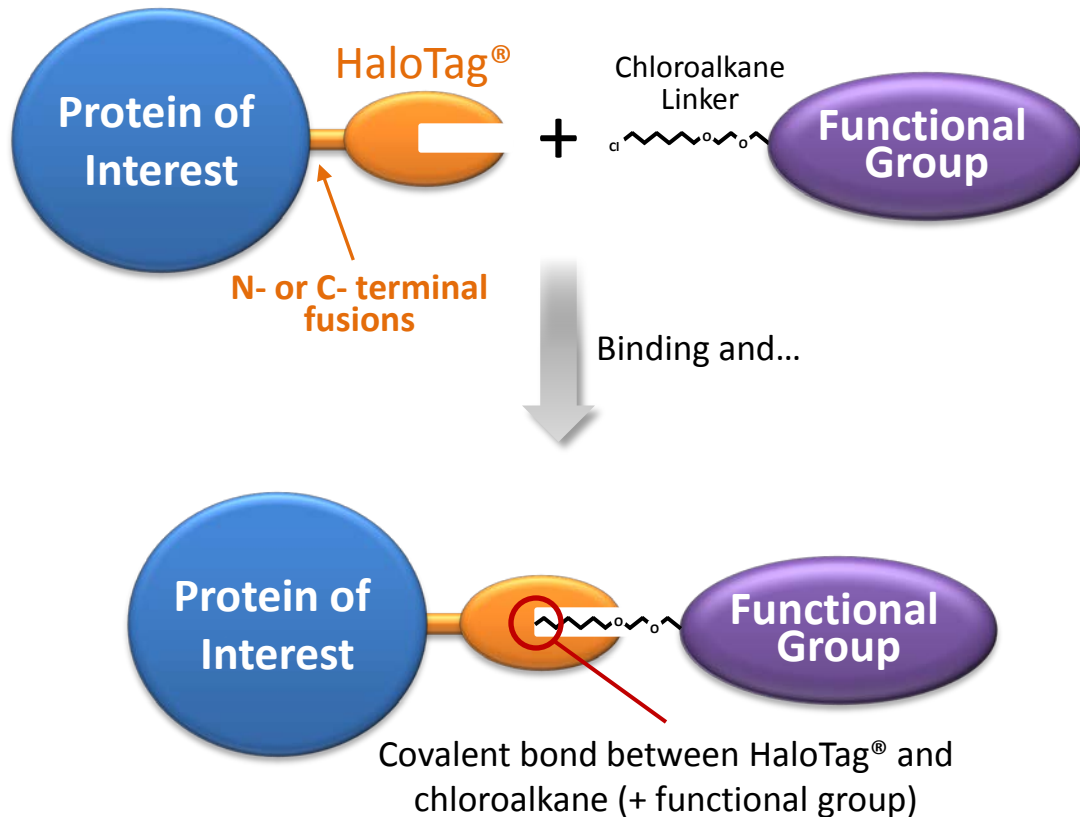
## A Unique, Multifunctional Protein Fusion Tag



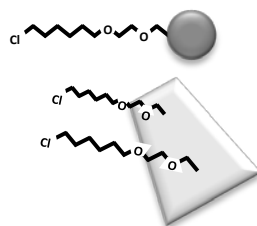
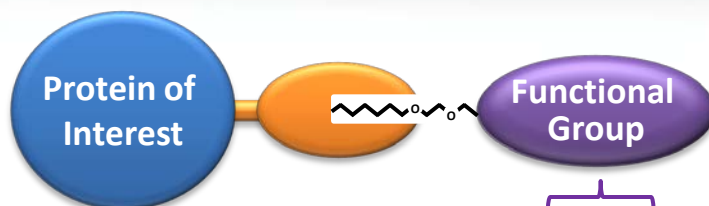
### HaloTag®:

- Engineered 34.1kDa halophilic bacterial hydrolase
- Binds to chloralkane substrate and locks with covalent attachment
- Faster kinetics than the biotin:streptavidin interaction
- No homolog in mammalian cells = no background

Read more about the development of this powerful fusion tag:  
Ohana, R.F., *et al.*  
(2009) *Prot. Exp. Purif.*  
**68**, 110-120.



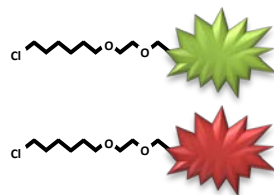
# Many Functional Groups are Available to Match Your Research Application(s)



## HaloTag® Surfaces/Resins

### *Capture and Display*

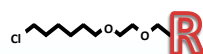
- ✓ Protein arrays
- ✓ Purification
- ✓ Interaction analysis



## HaloTag® Fluorescent Ligands

### *Labeling and Detection*

- ✓ Cellular imaging
- ✓ Gel analysis
- ✓ Quantitation

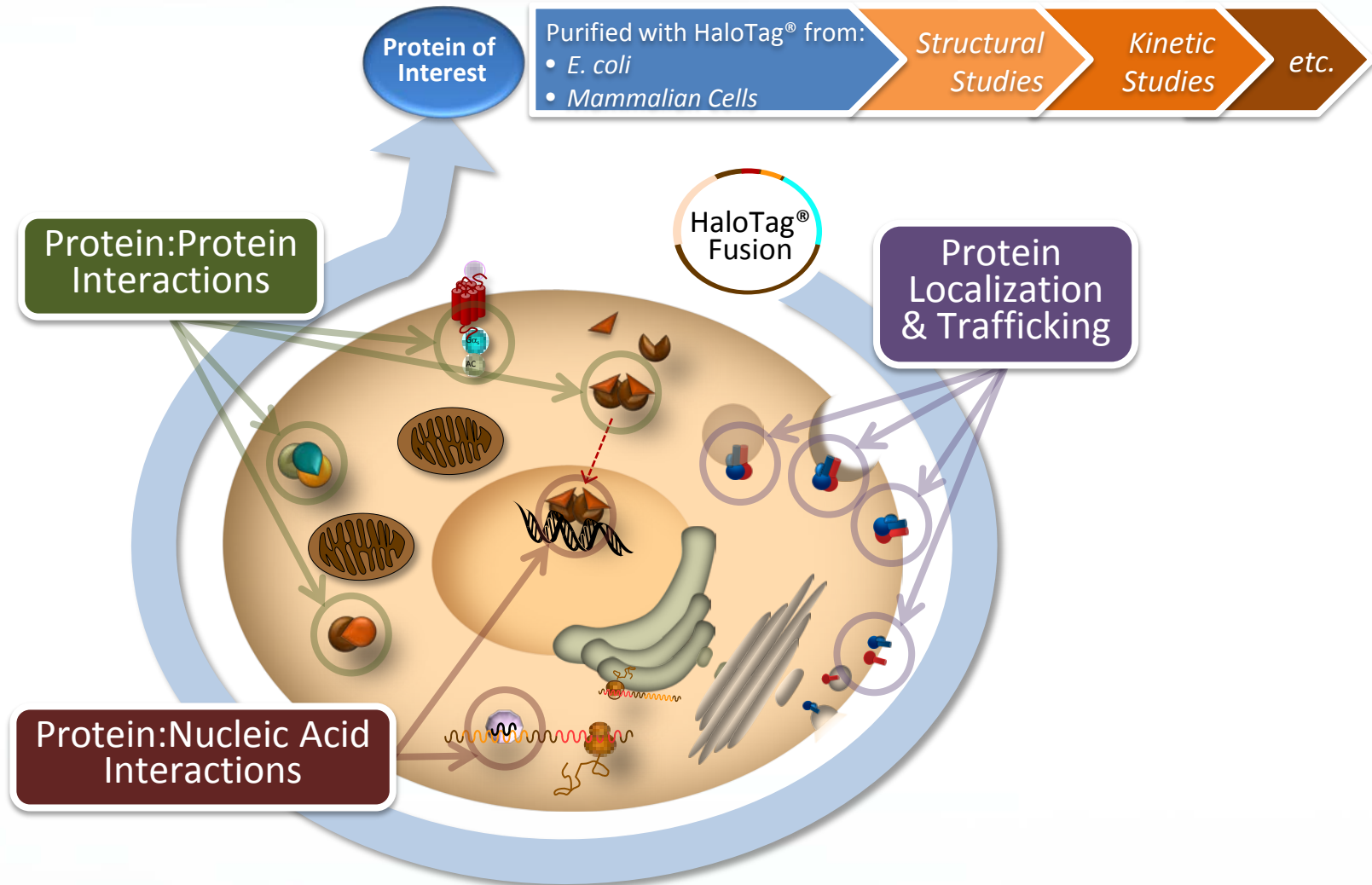


## HaloTag® Reactive Ligands

### *Custom Modifications*

- ✓ Attach to particles, surfaces
- ✓ Attach special ligands

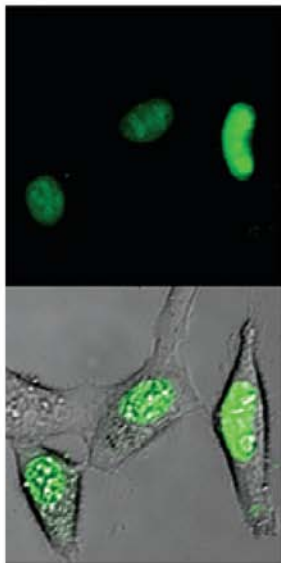
# One HaloTag® Fusion Protein = Global Protein Characterization Capabilities



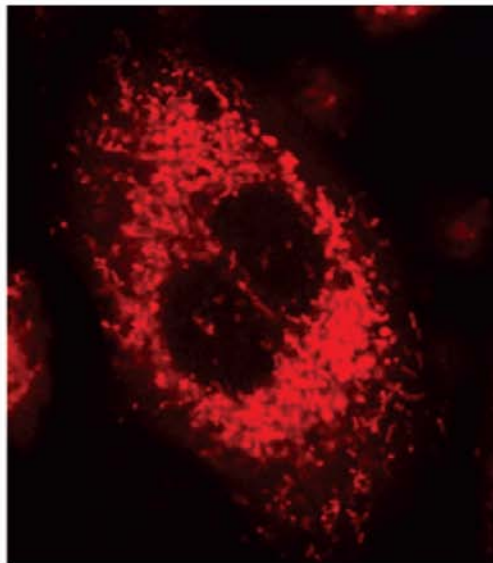
# ***HaloTag® Fusions Go & Are Detectable Anywhere*** *Examples Using Various Fluorophore Ligands*



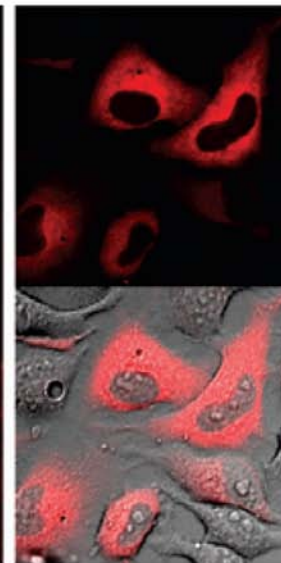
Nucleus  
HaloTag®-NLS<sub>3</sub>



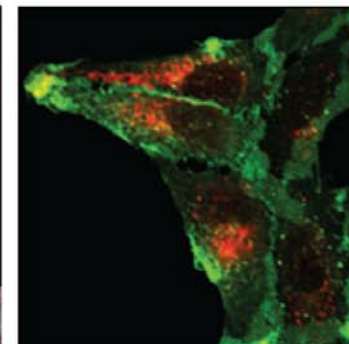
Mitochondria  
Mito-HaloTag  
Ligand



Cytosol  
p65-HaloTag  
Ligand



Membrane  
HaloTag-ECS



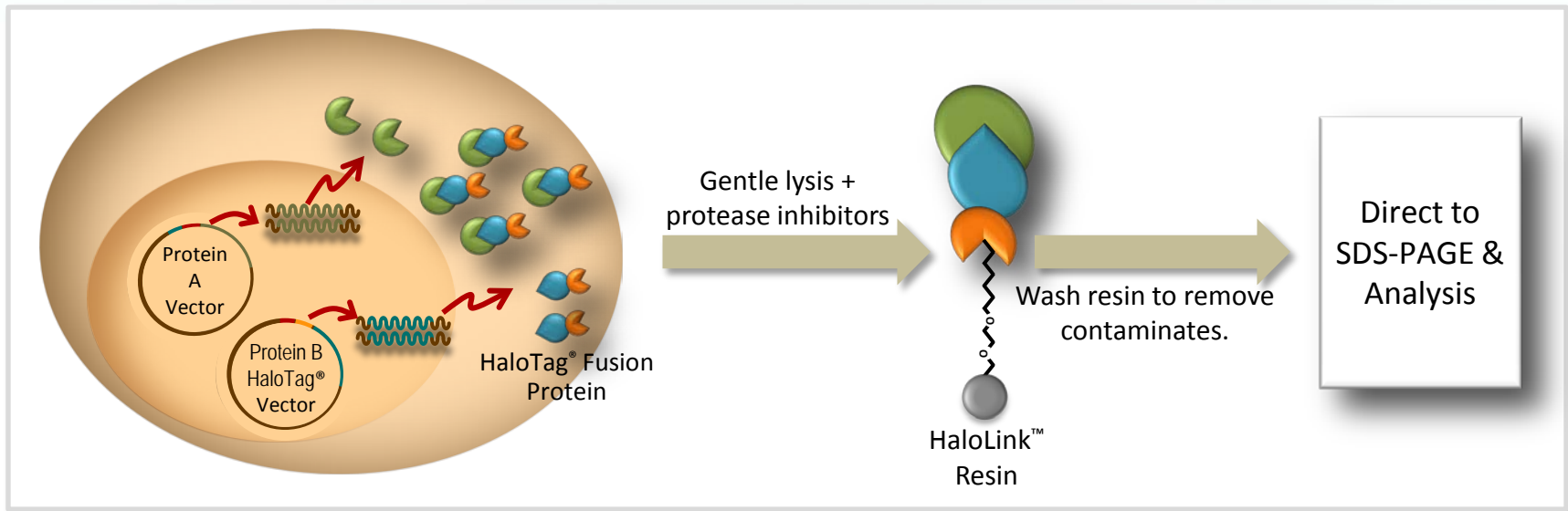
9771TA

4975TA

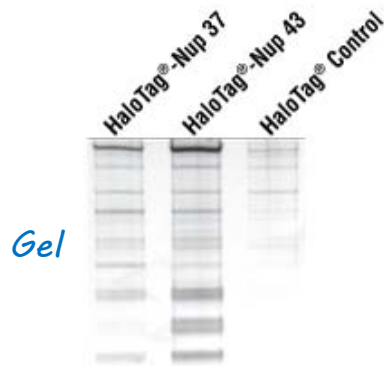


# Simple, Effective Mammalian Pulldown Assays

## Affinity Purification of the Nup 107-160 Complex



### Capture of Proteins Associated with HaloTag®-Nup 37 and -Nup 43 Fusions



#### Proteins Identified by LC/MS/MS Analysis

HaloTag®-Nup 37	HaloTag®-Nup 43	HaloTag® Control
Nup 160	Nup 160	No Nup subunits detected
Nup 133	Nup 133	
Nup 107	Nup 107	
Nup 98/96	Nup 98/96	
Nup 85/75	Nup 85/75	
Nup 43	Nup 43	
Nup 37	Nup 37	
Nup TPR	Nup TPR	
NUDC		

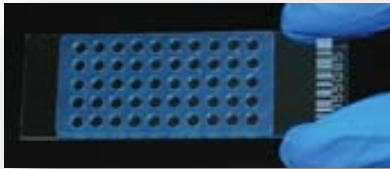
Méndez, J., et al. (2010) [Internet] [cited: 2011; July 20]. Available from:

<http://www.promega.com/resources/articles/pubhub/efficient-isolation-identification-and-labeling-of-intracellular-mammalian-protein-complexes/>

# Easy to Build and Customize Protein Arrays with HaloLink™ Array Slides



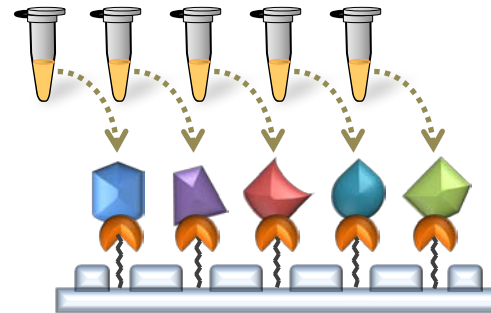
## HaloLink™ Array Slide



- Immobilize ≤50 proteins per slide
- HaloTag® fusion dictates orientation
- Easy generation/customization of bait proteins using cell-free expression

Hurst, R., *et al.* (2009) *Analytical Biochemistry* **392**, 45-53.

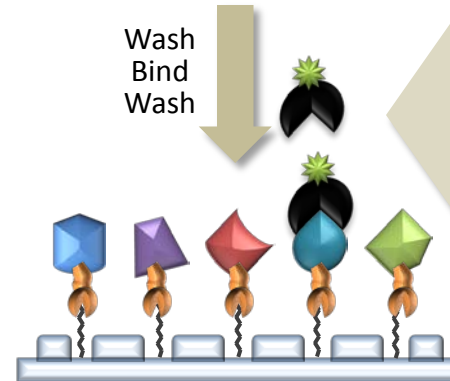
TNT® SP6 High-Yield WGE rxn



Synthesize HaloTag® fusion bait proteins using TNT® SP6 WG System

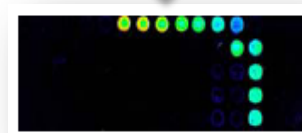
Immobilize  
1 bait/well

Wash  
Bind  
Wash



- FluoroTect™ label
- Transcend™ label
- HaloTag® Ligand-labeled
- Radioactive label
- No label: Ab detection

Detection/Visualization

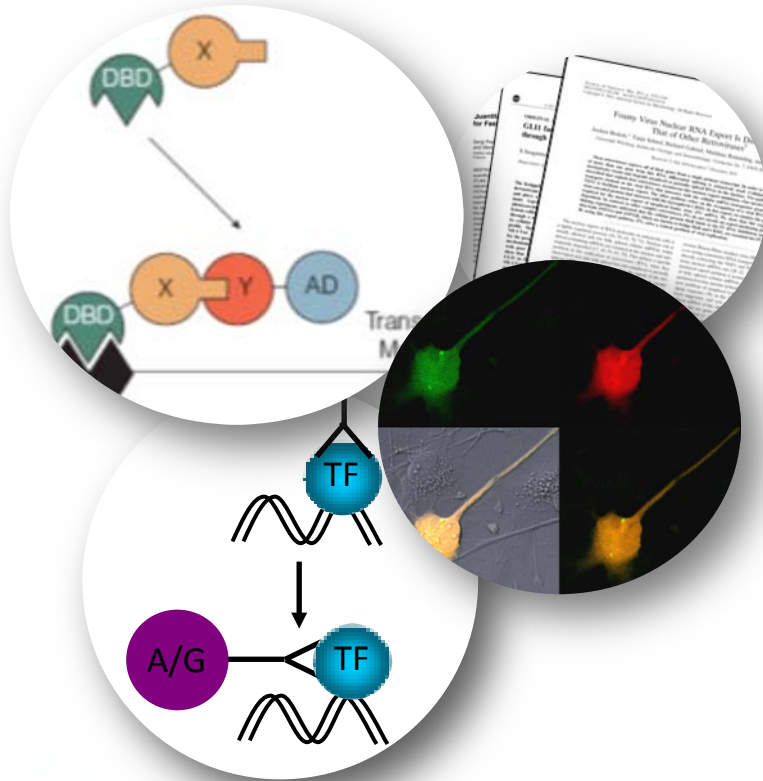


***Applying In Vitro (Cell-Free) Expression  
& HaloTag® Fusions to the Study  
Protein:Protein Interactions***



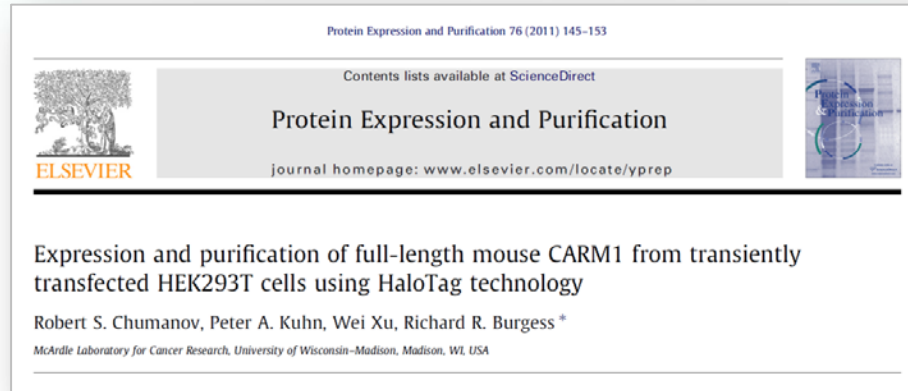
# Discovery Studies

## Identification of New In Vivo Protein Partners



- Protein affinity purification
  - All in vivo
  - In vitro with purified bait + cellular extracts
- Yeast Two-hybrid assays
  - Library screening
- etc.

# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



Chumanov, R., et al. (2011) *Prot. Exp. Purif.* **76**, 145-53

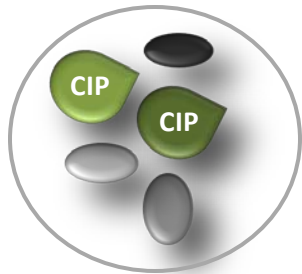
- **CARM1** (PRMT4)
  - Protein arginine methyl transferase
  - Transfers methyl group from SAM to Arg residues
  - Coactivator of transcriptional activation
- CARM1 studies have been hampered by inability to purify full-length protein
- **Goal:** Identify target substrates of CARM1 using affinity purification of interacting proteins followed by in vitro methylation experiments

# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



## Experimental Design - In Vitro Protein Affinity Purification

Nuclear Extract



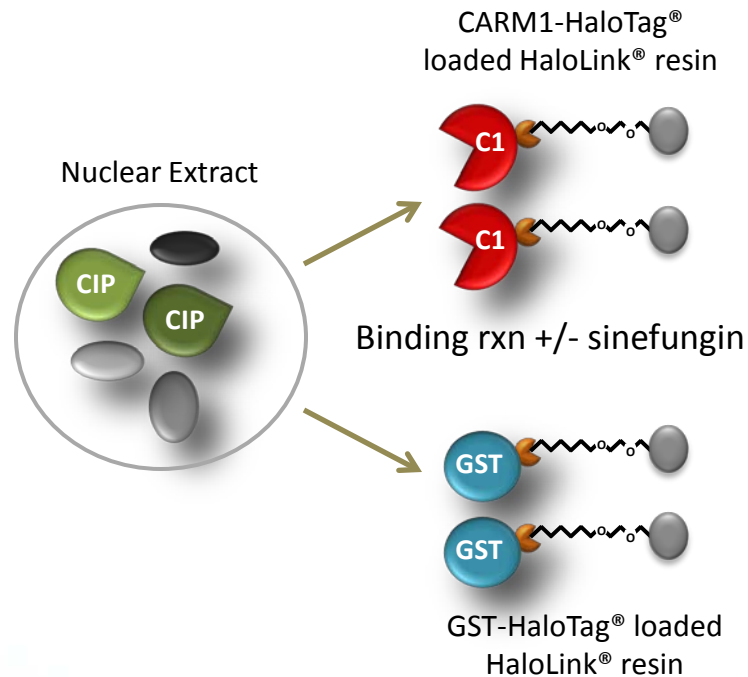
**C1** = CARM1

**CIP** = CARM1 Interacting Protein

# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



## Experimental Design - In Vitro Protein Affinity Purification



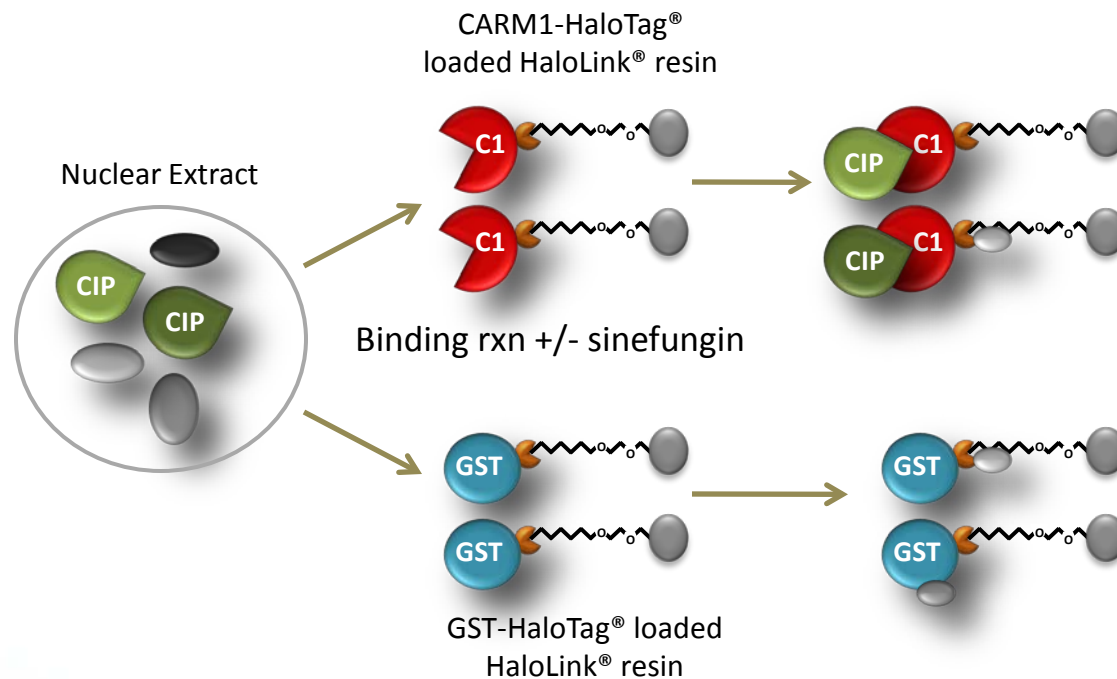
**C1** = CARM1

**CIP** = CARM1 Interacting Protein

# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



## Experimental Design - In Vitro Protein Affinity Purification



**C1** = CARM1

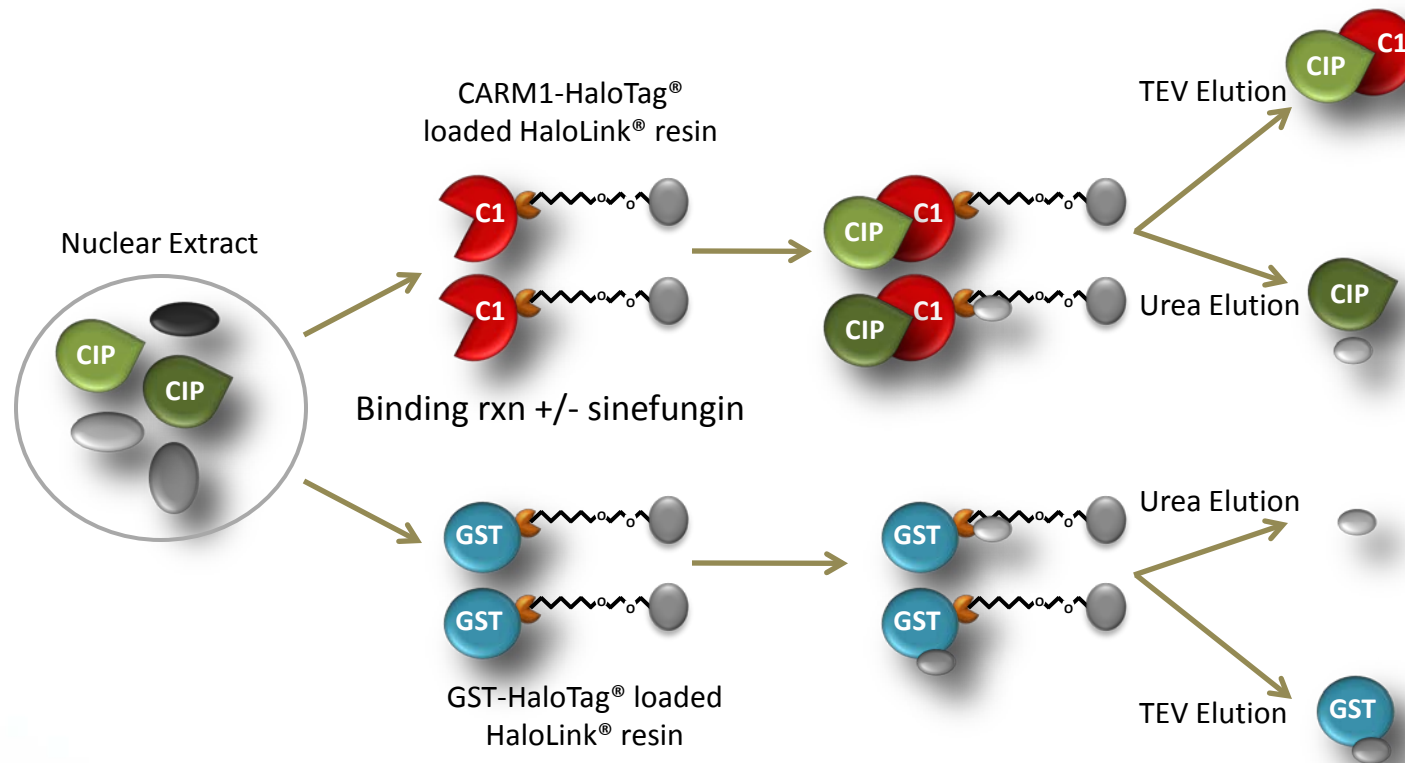
**CIP** = CARM1 Interacting Protein



# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



## Experimental Design - In Vitro Protein Affinity Purification



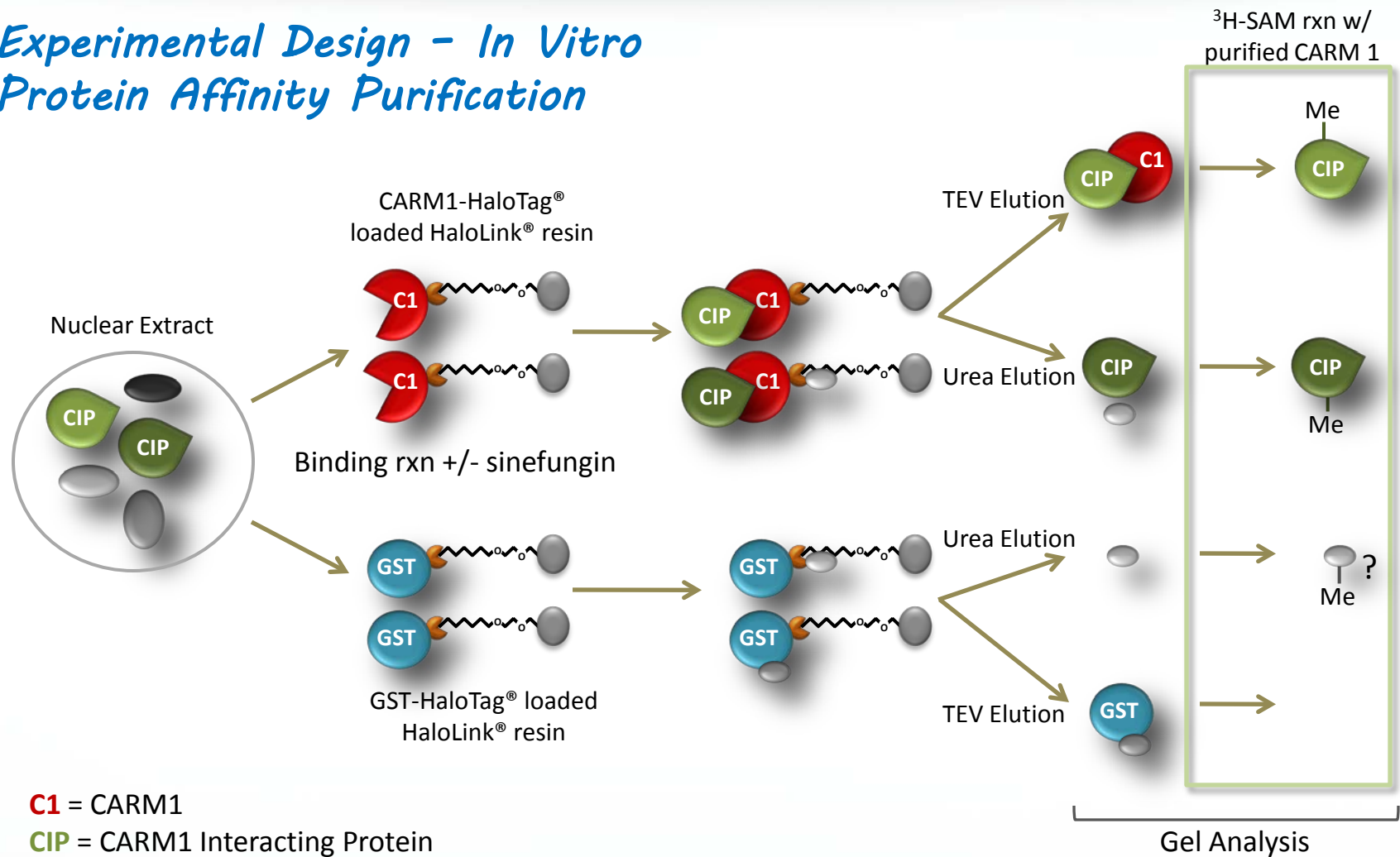
**C1** = CARM1

**CIP** = CARM1 Interacting Protein

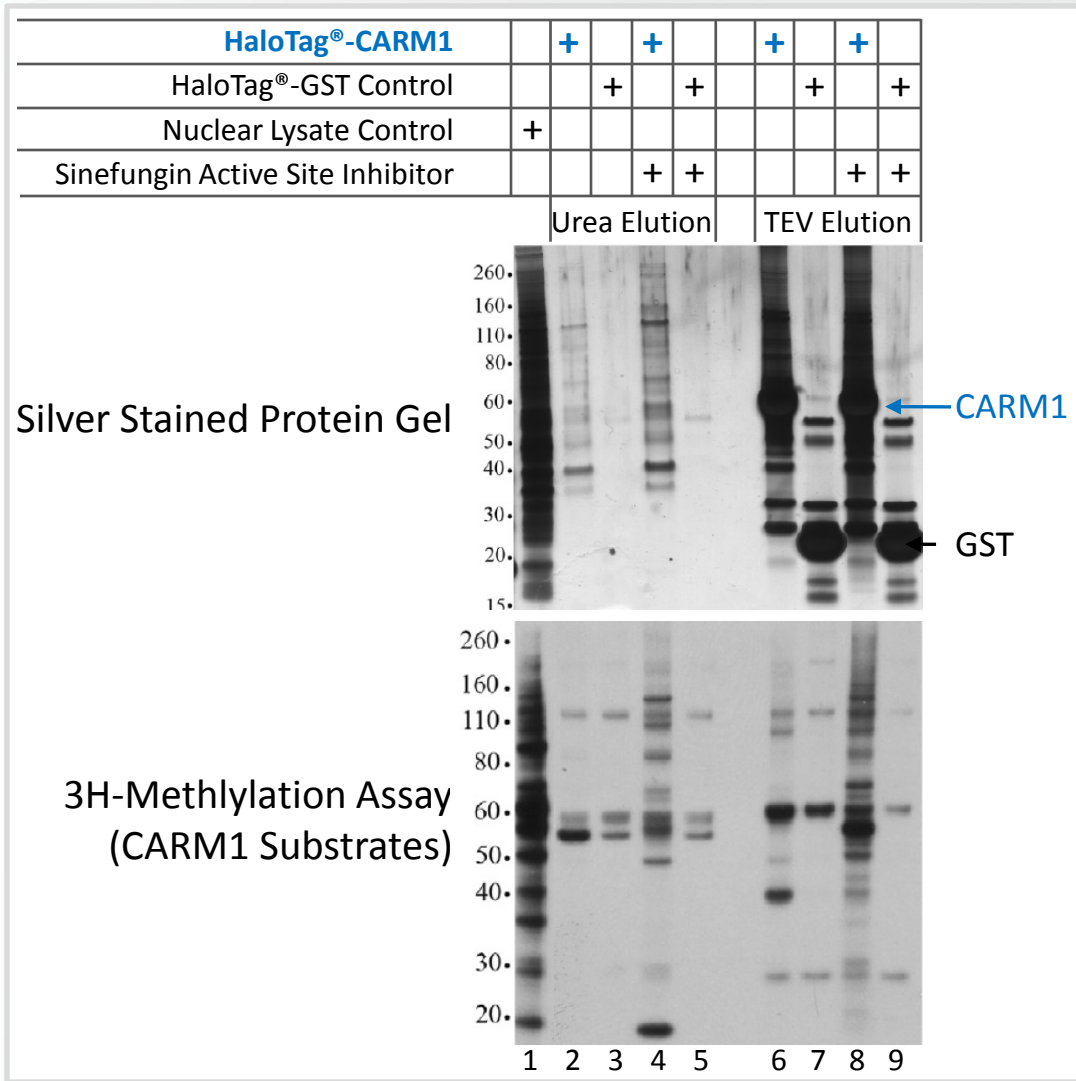
# Case Study: Identification of New CARM1 Interacting Proteins/Substrates



## Experimental Design - In Vitro Protein Affinity Purification



# Case Study: Identification of New CARM1 Interacting Proteins/Substrates

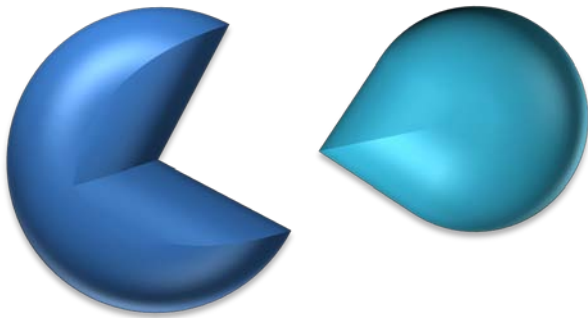


## Results

1. Identified multiple CARM1 Interacting Proteins
2. Many CIPs are also CARM1 substrates
3. Sinefungin increases capture efficiency of CIPs/substrates

# **Verification Studies**




## *Confirming Two (or more) Proteins Interact*



- In Vitro Confirmation
  - Protein affinity purification (pulldown assays)
  - Co-Immunoprecipitations
- In Vivo (cells) Confirmation
  - Protein affinity purification (pulldown assays)
  - Co-immunoprecipitations
  - Mammalian two-hybrid assays

# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



**BJ** [www.biochemj.org](http://www.biochemj.org)  This is a data-enriched, interactive PDF that provides the gateway to a world of information when opened in Utopia Documents. [Download FREE software now](#)   **387**

Biochem. J. (2011) **436**, 387–397 (Printed in Great Britain) doi:10.1042/BJ20101748

---

**The novel Nrf2-interacting factor KAP1 regulates susceptibility to oxidative stress by promoting the Nrf2-mediated cytoprotective response**

Atsushi MARUYAMA\*, Keizo NISHIKAWA†, Yukie KAWATANI‡, Junsei MIMURA\*, Tomonori HOSOYA\*, Nobuhiko HARADA\*, Masayuki YAMAMATO‡ and Ken ITOH\*<sup>1</sup>

\*Department of Stress Response Science, Center for Advanced Medical Science, Hirosaki University Graduate School of Medicine, 5 Zaifu-cho, Hirosaki, Aomori 036-8562, Japan, †Laboratory of Stem Cell Regulation, National Institute of Biomedical Innovation, 7-6-8 Saito, Asagi, Ibaraki, Osaka 567-0085, Japan, and ‡Department of Medical Biochemistry, Tohoku University Graduate School of Medicine, 2-1 Seiryō-cho, Aoba-ku, Sendai 980-8575, Japan

Maruyama A., et al. (2011) *Biochem. J.* **436**, 387-97

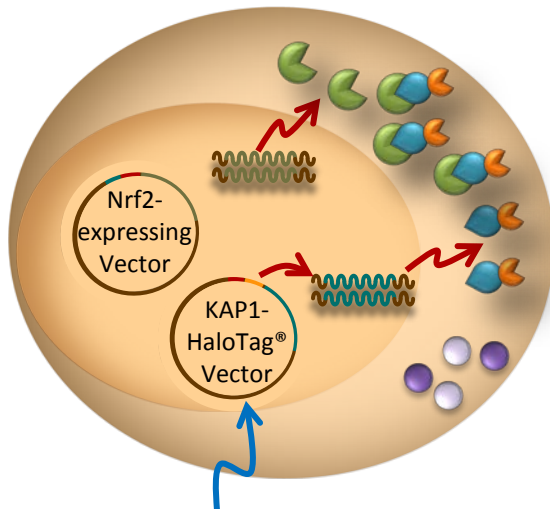
- **Nrf2** (Nuclear factor-erythroid 2-related factor)
  - Transcription factor
  - Co-ordinately regulates ARE (antioxidant-response element)-mediated induction of cytoprotective genes in response to oxidative stress & electrophiles
- Mechanism of action is not well understood
- **Goal:** After identification of KAP1 as an Nrf2 interacting protein, confirm and characterize the interaction

# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



## Experimental Design - Confirming the Interaction with a HaloTag® Pulldown Assay

HEK293



Validated HaloTag® clones  
available from

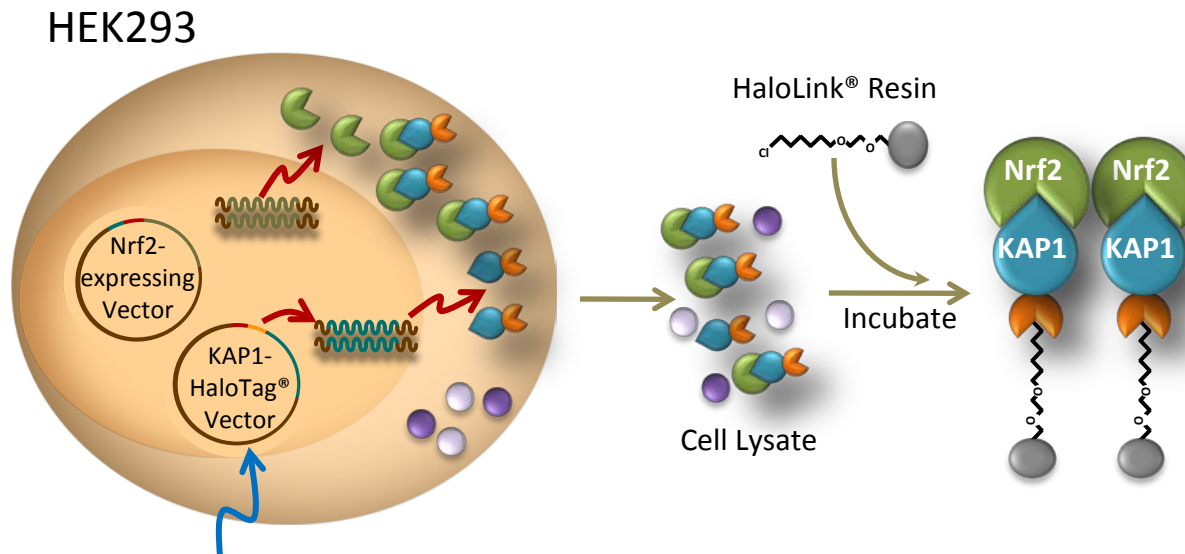
[www.promega.com/kazusa](http://www.promega.com/kazusa)

A partnership between Kazusa  
DNA Research Institute and  
Promega.

# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



## Experimental Design - Confirming the Interaction with a HaloTag® Pulldown Assay



Validated HaloTag® clones  
available from

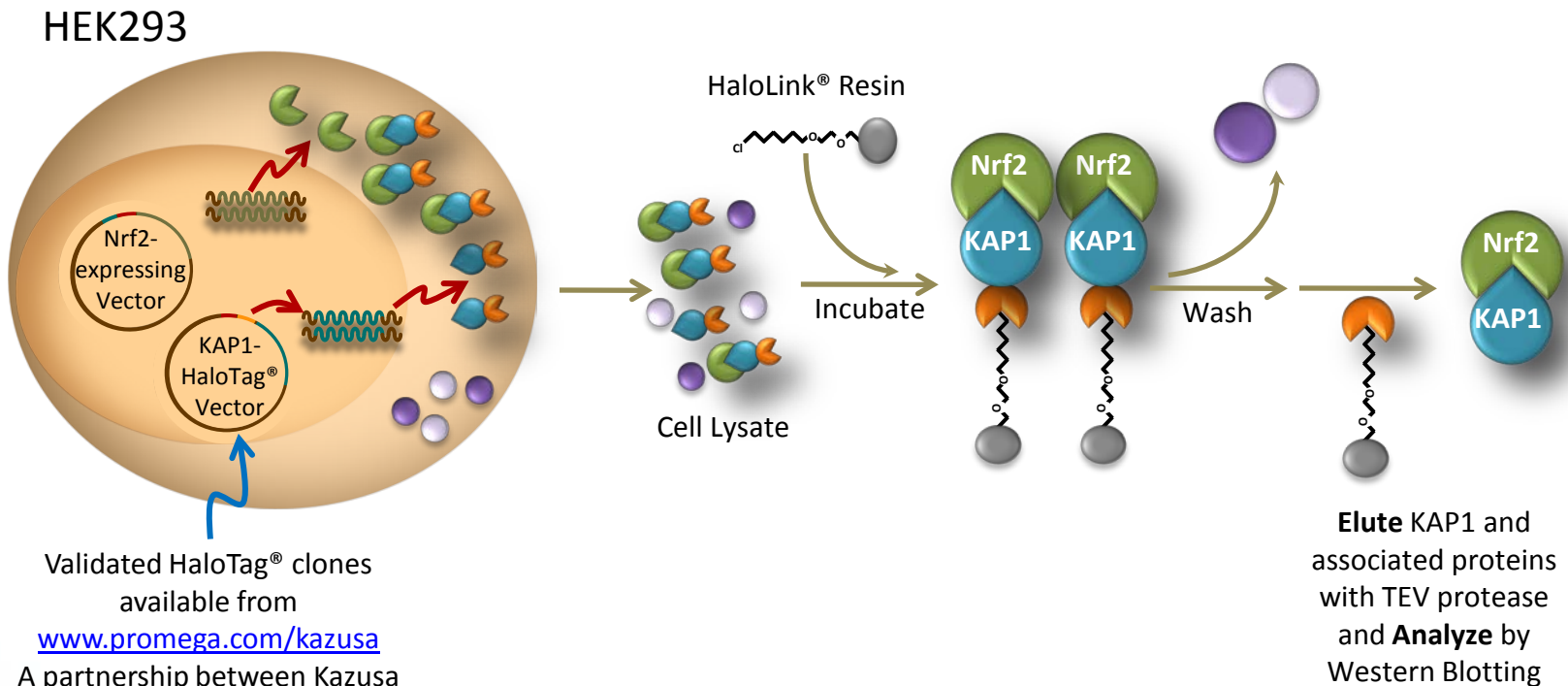
[www.promega.com/kazusa](http://www.promega.com/kazusa)

A partnership between Kazusa  
DNA Research Institute and  
Promega.

# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction

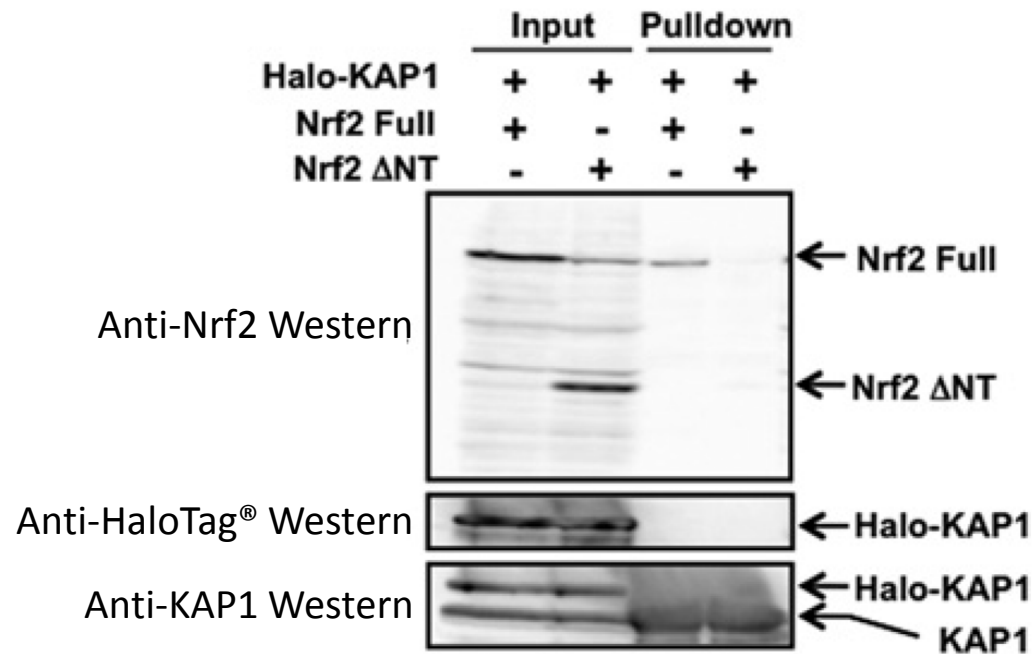


## Experimental Design - Confirming the Interaction with a HaloTag® Pulldown Assay





# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



Data from Maruyama A., et al. (2011) *Biochem. J.* **436**, 387-97

## Results

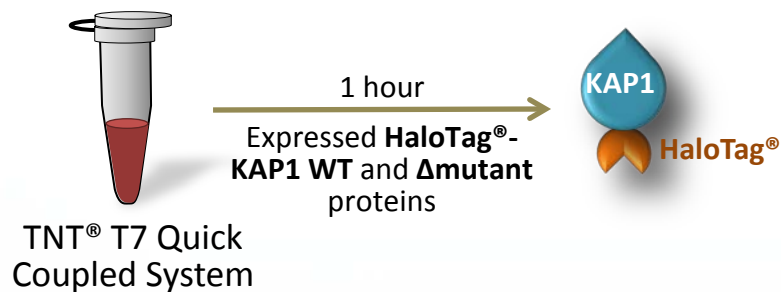
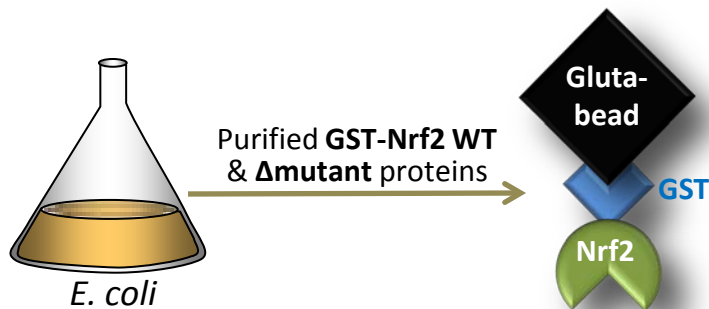
1. Verified KAP1:Nrf2 interaction *in vivo*

2. N-terminus of Nrf2 is required for interaction with KAP1

# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



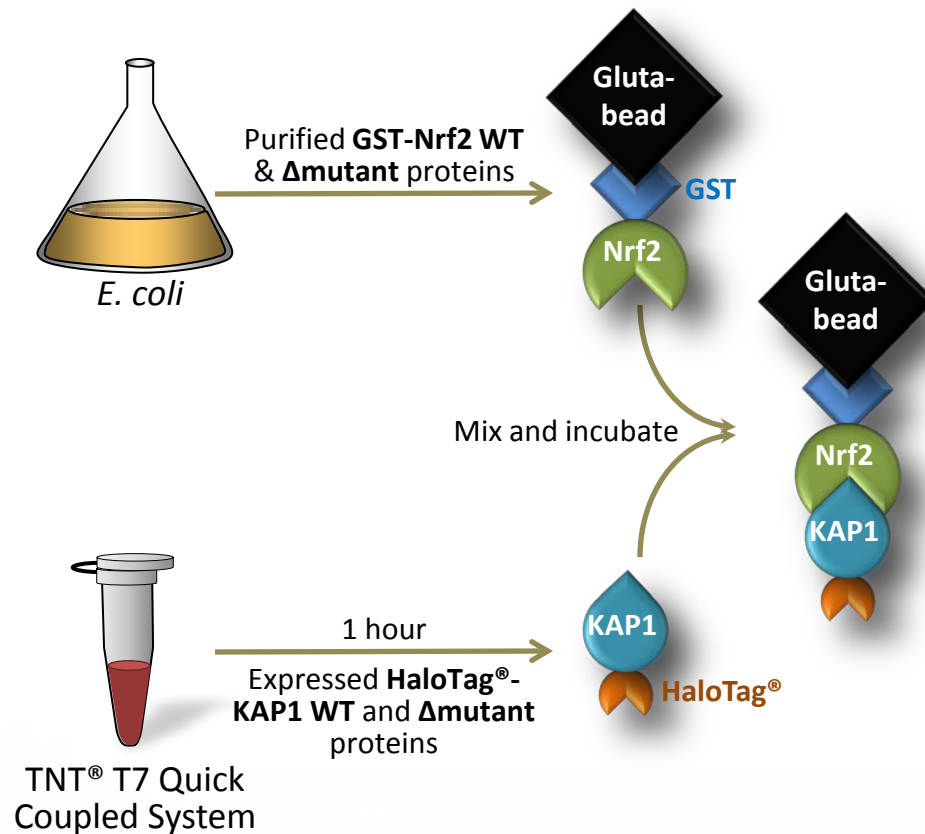
## Experimental Design - Characterizing the Interaction with GST Pulldowns



# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



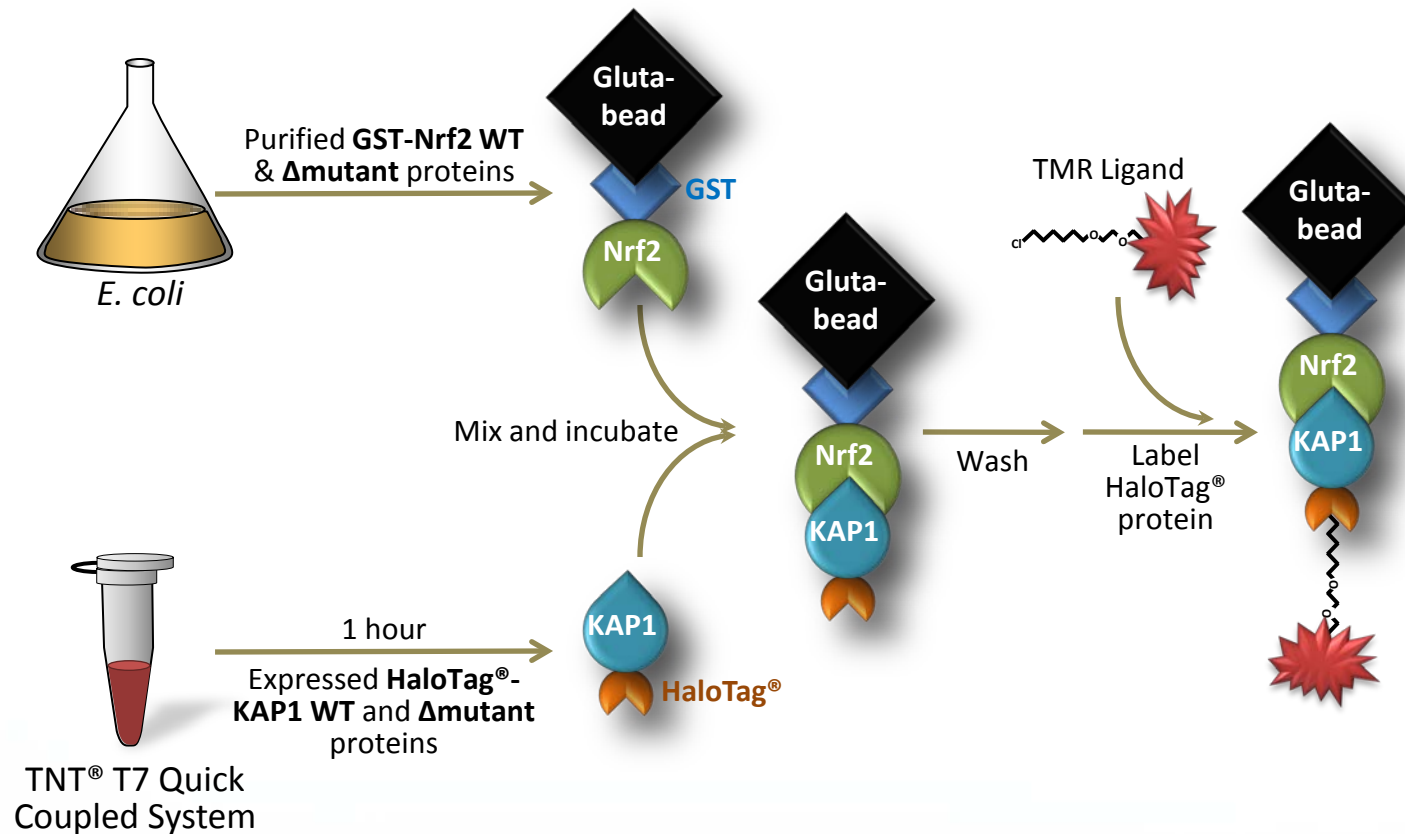
## Experimental Design - Characterizing the Interaction with GST Pulldowns



# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



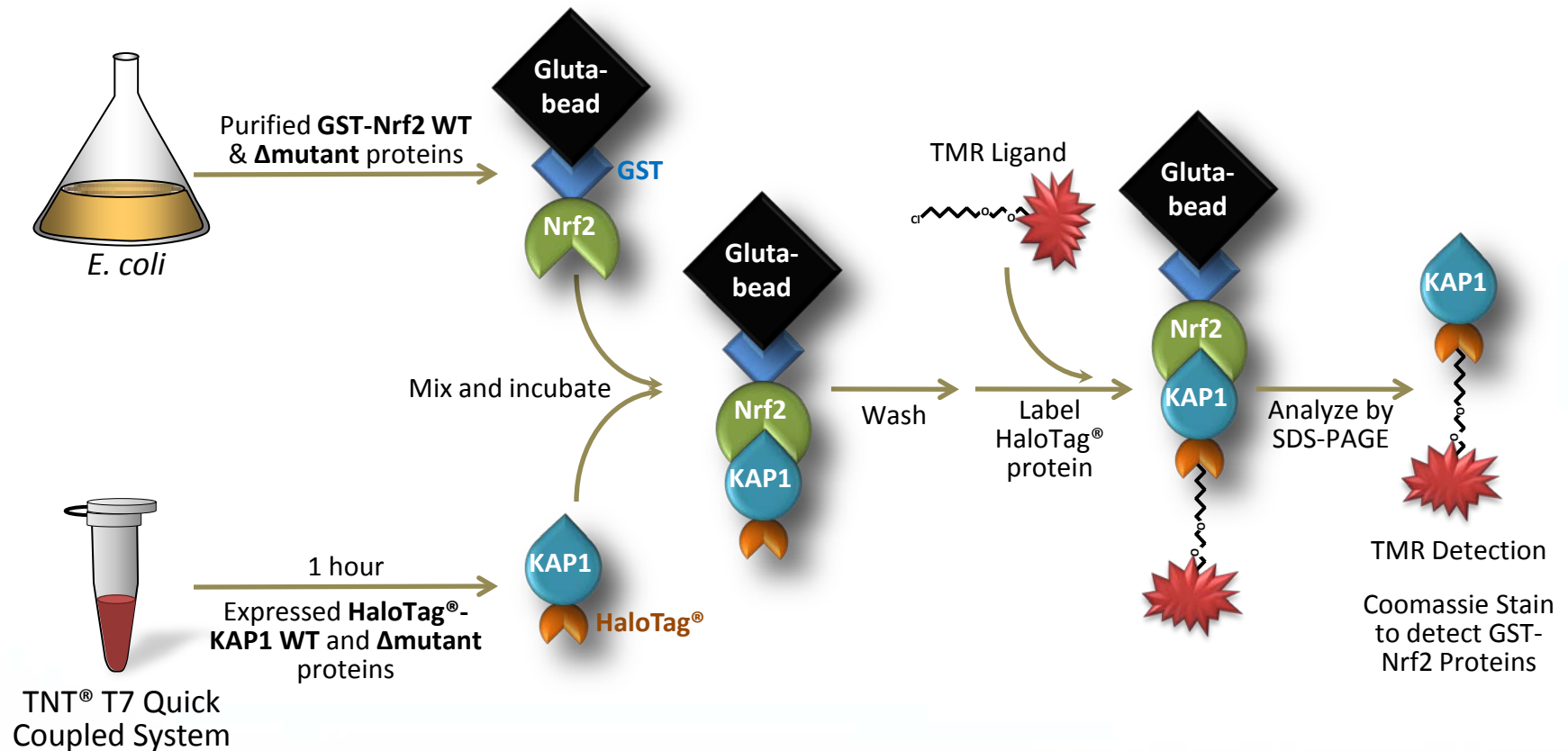
## Experimental Design - Characterizing the Interaction with GST Pulldowns



# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction



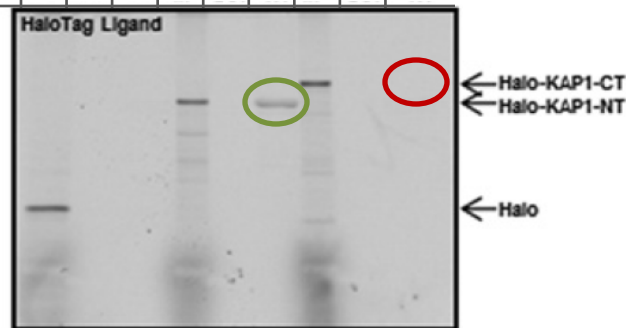
## Experimental Design - Characterizing the Interaction with GST Pulldowns



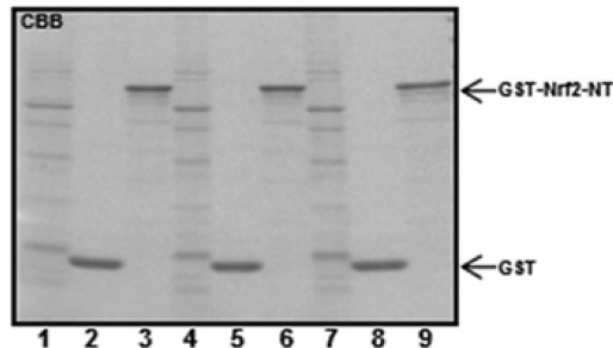
# Case Study: Verification & Characterization of a Novel Nrf2:KAP1 Interaction

HaloTag® Control	IN	+	+						
HaloTag®-KAP1 NT				IN	+	+			
HaloTag®-KAP1 CT							IN	+	+
GST		+			+			+	
GST-Nrf2 NT			+			+			+

TMR Detection  
(HaloTag®-KAP1)  
(Fluoroimager)



Coomassie Stained Gel  
(GST, GST-Nrf2)



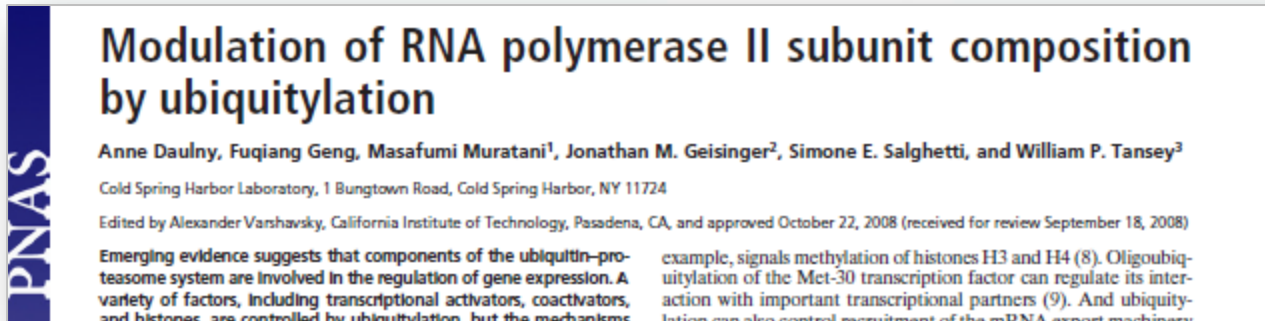
Data from Maruyama A., et al. (2011) *Biochem. J.* **436**, 387-97

## Results

1. *N-terminus (NT) of KAP1 promotes the interaction with Nrf2*

2. *Further analysis showed that activation domain within Nrf2 NT is responsible for interaction*

# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD



Daulny A., et al. (2008) *Proc. Nat. Acad. Sci.* **105**, 19649-54

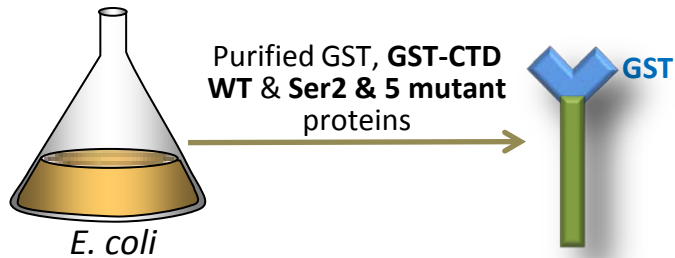
## • Asr1

- Involved in the yeast alcohol stress response
- Two-hybrid data on a mammalian protein with limited Asr1 homology suggested a possible role of Asr1 in ubiquitylation of RNAP II in yeast.
- Interested in understanding if Asr1 does indeed play a role in ubiquitylation of RNAP II and initial experiments in the paper demonstrated Asr is a ubiquitin-ligase.
- **Goal:** Demonstrate direct binding of Asr1 to the carboxy terminal domain (CTD) of RNAP II and characterize the interaction based on phosphorylation of the CTD.

# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD



## Experimental Design - Use Far Westerns to Detect Interactions

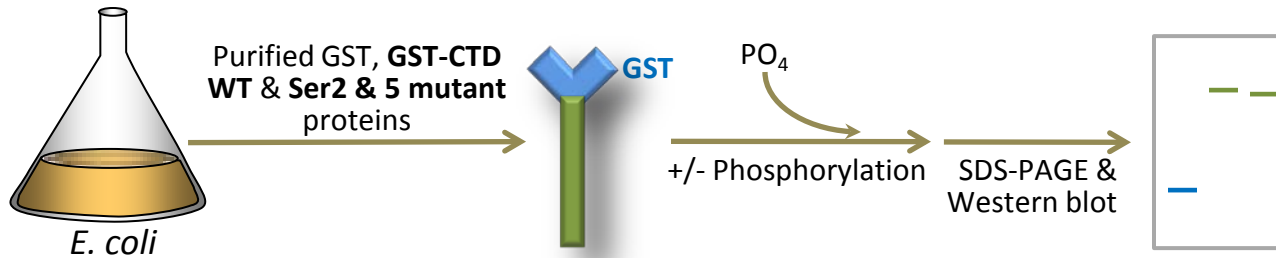




# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD

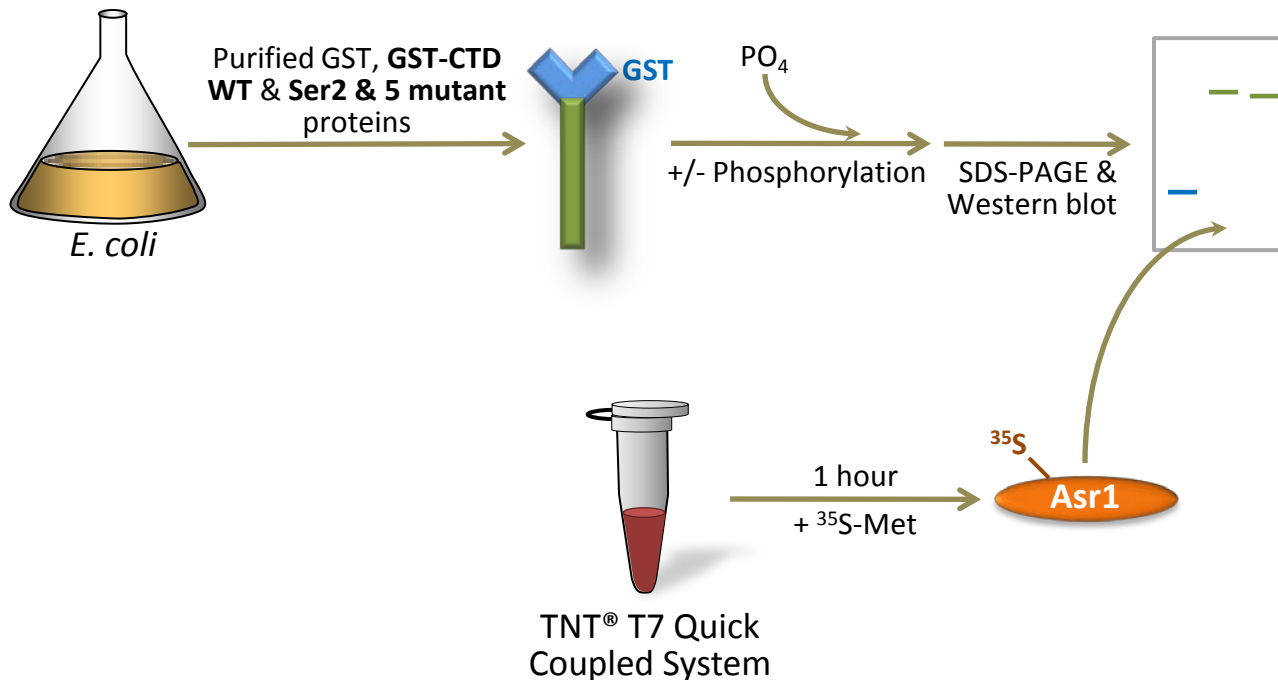


## Experimental Design - Use Far Westerns to Detect Interactions



# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD

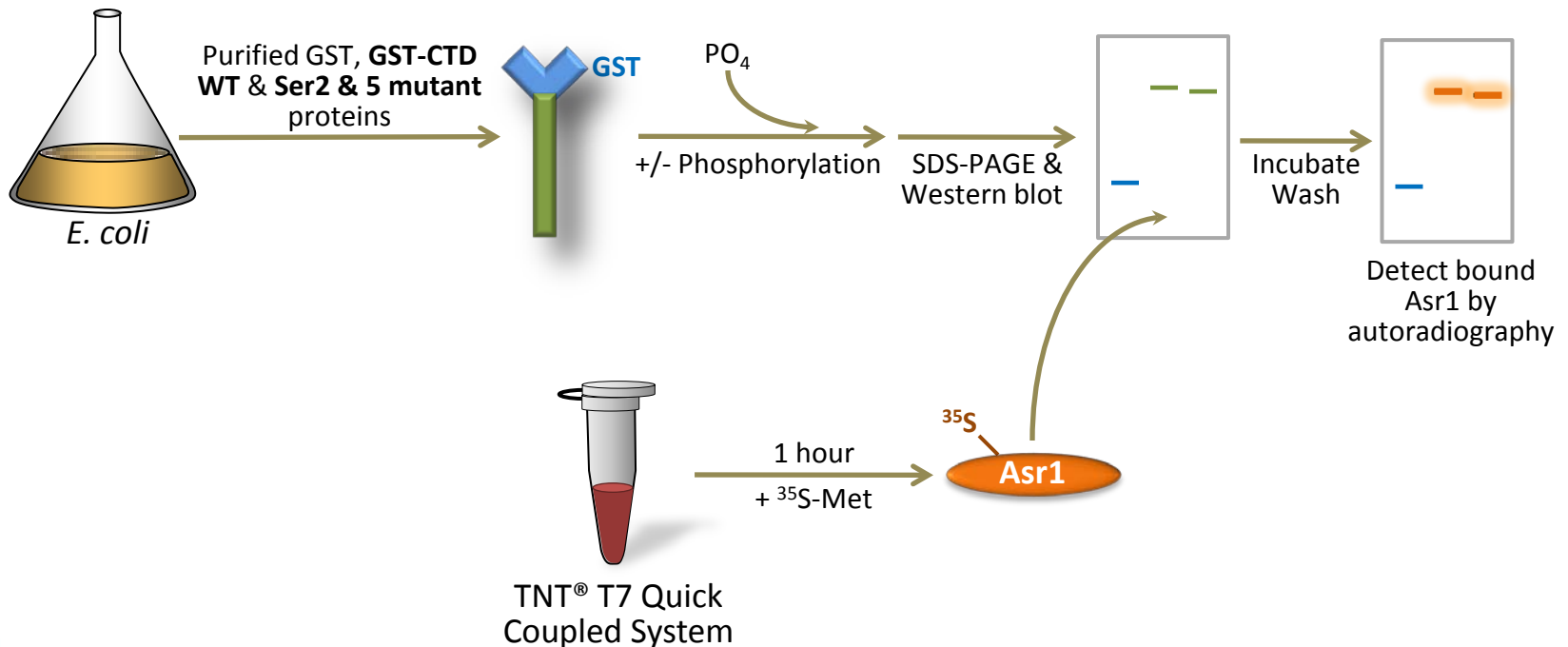
## Experimental Design - Use Far Westerns to Detect Interactions



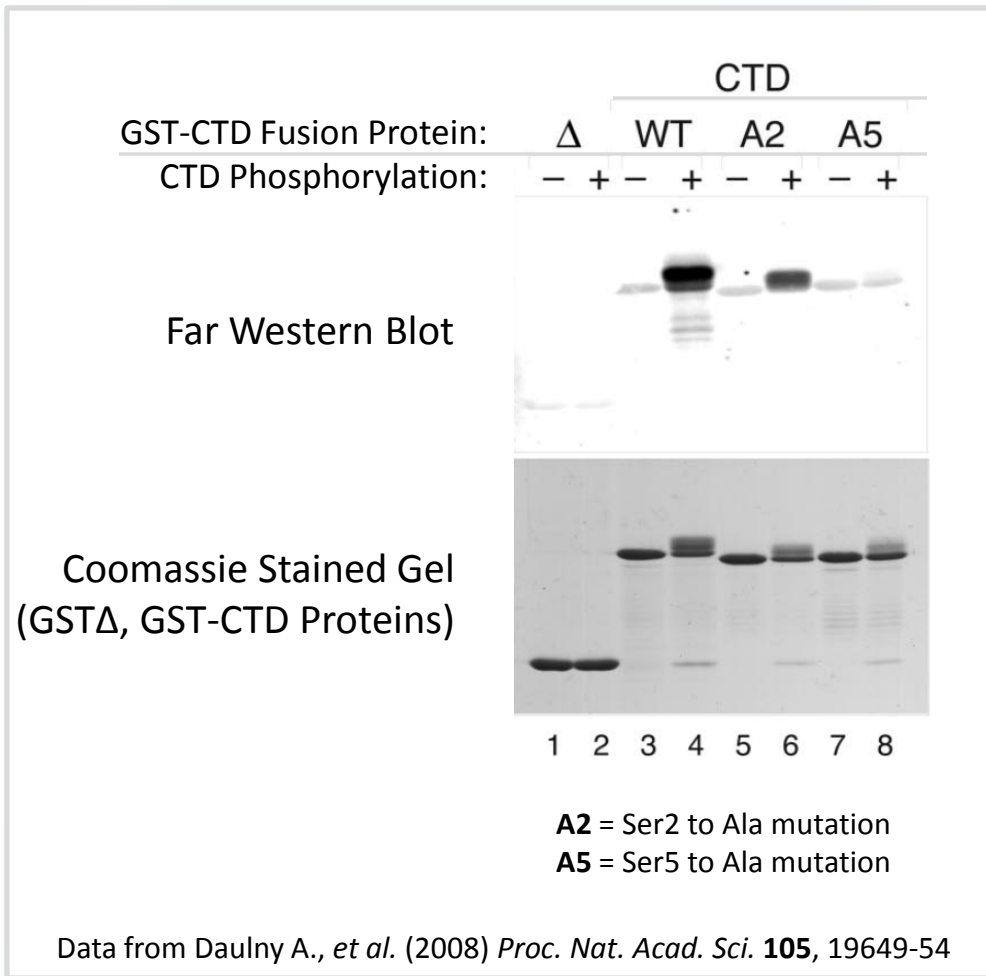
# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD



## Experimental Design - Use Far Westerns to Detect Interactions



# Case Study: Testing for Direct Interaction of Asr1 with the RNAP II CTD



## Results

1. *Asr1* binds directly to the RNAP II CTD in a phosphorylation dependent manner.
2. Mutation of Ser5 to alanine prevents phosphorylation at Ala5 & blocks *Asr1* binding.
3. Phosphorylation of Ser2 plays less of a role in promoting *Asr1* binding.

# ***Mammalian Two-Hybrid Assays***

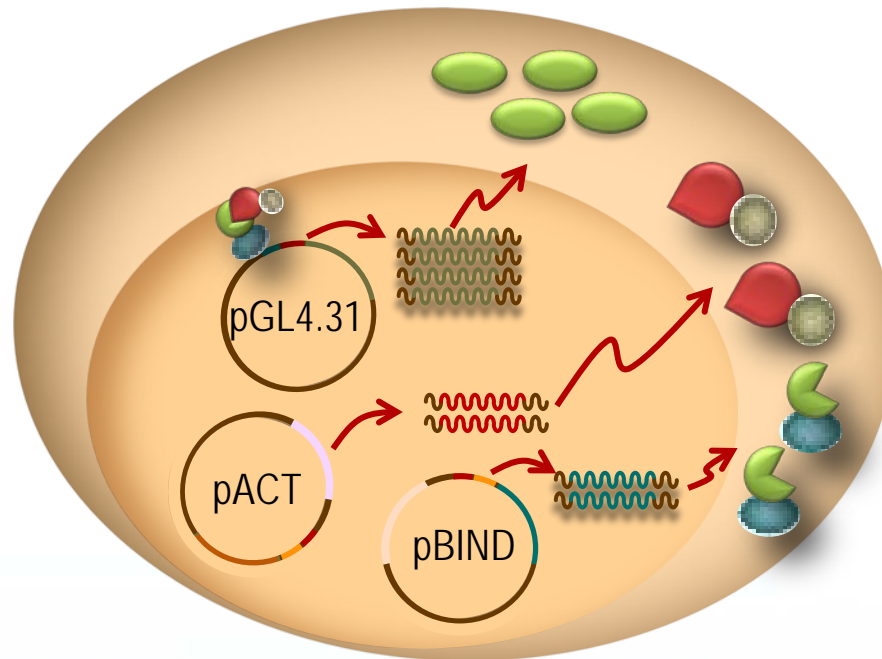
Adapting the Yeast Two-Hybrid Assay to Mammalian Cells



# Adaptation of the Yeast Two Hybrid Assay to Mammalian Cells: CheckMate™



- 1 TRANSFECT
- 2 CULTURE 2-3 DAYS
- 3 DUAL-LUCIFERASE ASSAY



15 citations  
in 2011  
Jan-July

# Case Study: Testing for Interaction of Cdk3 with ATF1 in Mammalian Cells



## Research Article

### Cyclin-Dependent Kinase 3–Mediated Activating Transcription Factor 1 Phosphorylation Enhances Cell Transformation

Duo Zheng, Yong-Yeon Cho, Andy T.Y. Lau, Jishuai Zhang, Wei-Ya Ma, Ann M. Bode, and Zigang Dong

The Hormel Institute, University of Minnesota, Austin, Minnesota

Zheng, D., *et al.* (2008) *Cancer Res.* **68**, 7650-60.

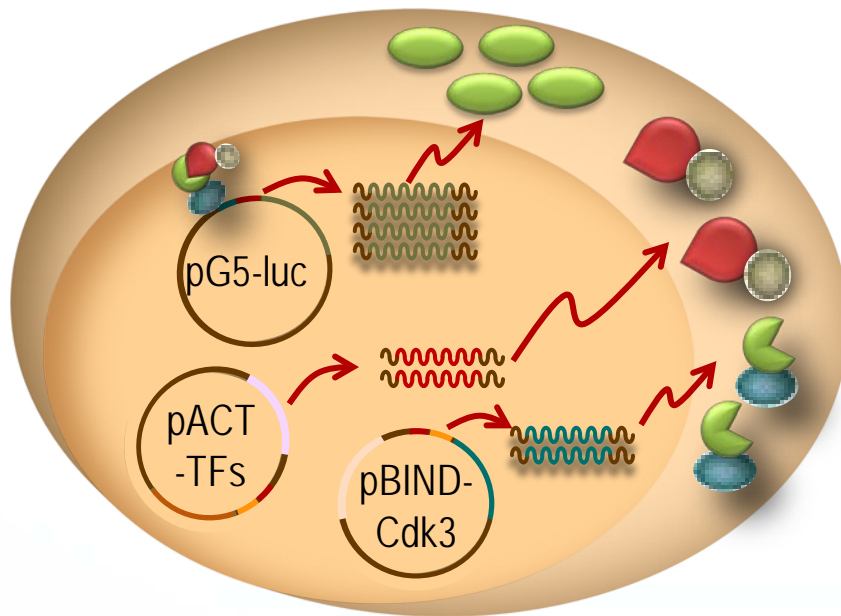
- **Cdk3** (Cyclin-dependent kinase 3)
  - Ser/Thr kinase
  - Highly expressed in glioblastoma tissues and cell lines
- Interested in determining if Cdk3 is interacting with and phosphorylating transcriptional regulatory proteins in glioblastoma cells to upregulate transcription
- **Goal:** Screen various transcription factors (TFs) using the CheckMate Mammalian Two Hybrid Assay to determine if they interact with Cdk3 in vivo.

# Case Study: Testing for Interaction of Cdk3 with ATF1 in Mammalian Cells

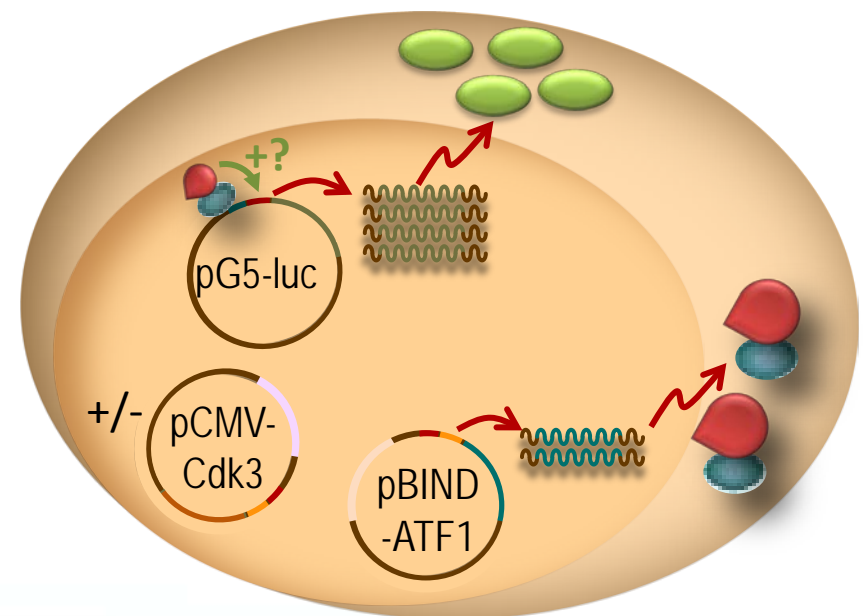


## Experimental Design

CheckMate™ Two Hybrid Assay



Modified CheckMate™ One Hybrid Assay

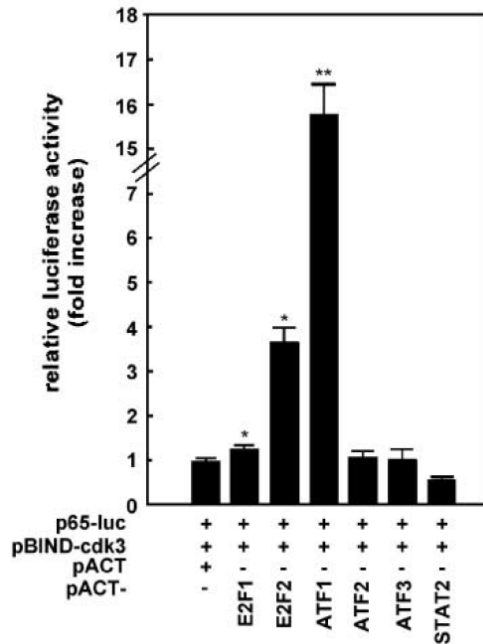




# Case Study: Testing for Interaction of Cdk3 with ATF1 in Mammalian Cells

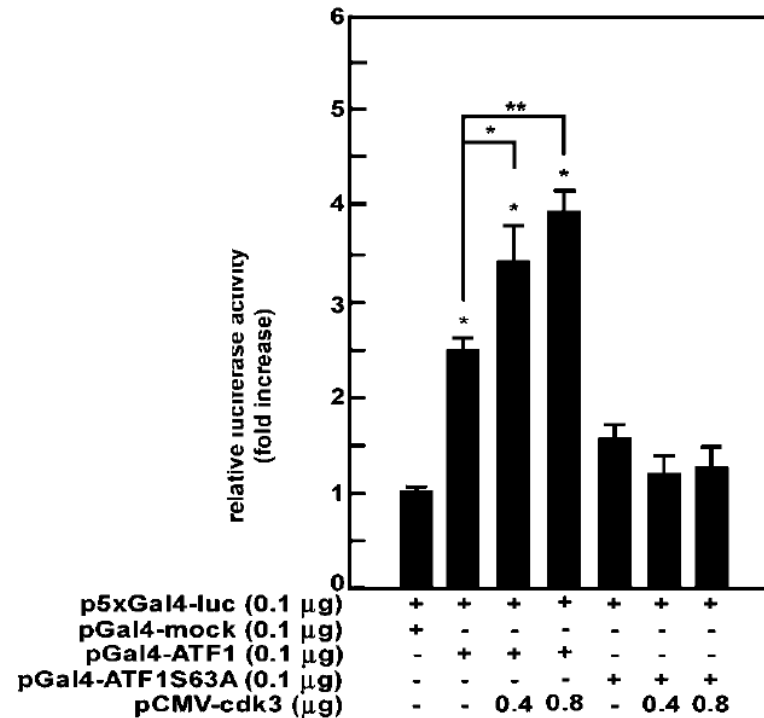


## Two-Hybrid Assay



Identified ATF1 as a new binding partner for Cdk3 (E2F2 = positive control)

## One-Hybrid Assay



Phosphorylation of ATF1 increases transcriptional activation activity

# ***Additional Applications of In Vitro Expression***



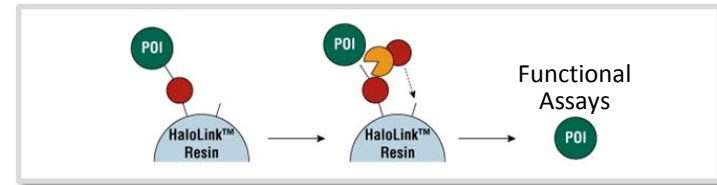
- Functional Assays
  - Enzymatic activity assays
    - Screening applications
- Post-translational modification analysis
- In Vitro Protein-DNA & Protein-RNA Interactions Studies
  - Gel-shift assays

# Additional Applications HaloTag<sup>®</sup> Technology

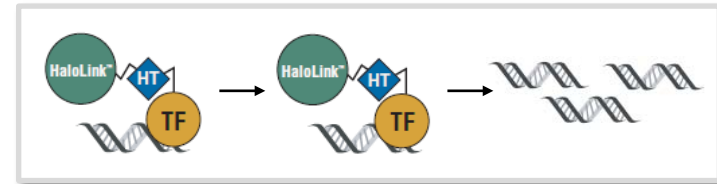


- HaloTag<sup>®</sup> Technology

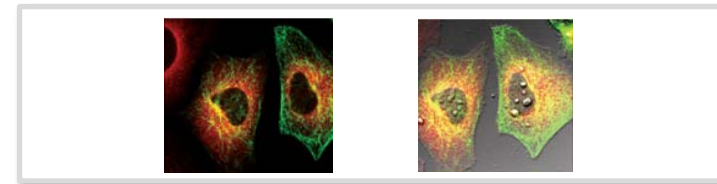
- Functional assays with protein purified from *E. coli* and mammalian cells



- Chromatin pull-down assays with HaloCHIP<sup>™</sup> System



- Protein localization, trafficking and turnover



- Target protein imaging in whole animals

## *Summary*



- In vitro (cell-free) expression provides a rapid means to produce full-length or deletions of your protein of interest for interaction studies.
- HaloTag<sup>®</sup> Fusions provide a multifunctional handle on your protein of interest to study protein interactions both in vitro and in vivo.
- The combination of in vitro (cell-free) expression and HaloTag<sup>®</sup> fusions is a powerful tool for protein interaction studies.
- The CheckMate<sup>™</sup> Mammalian Two-Hybrid System is another tool for studying protein interactions in mammalian cells.

# ***Promega Field Applications Specialists*** ***Your Local Promega Resource***



**Please Contact Us to Obtain Contact Information for Your Local FAS**

**United States:** 800-356-9526

**Canada:** 608-274-4330



# *Technical Services Scientists Ready to Help*



**By phone:** 800-356-9526

- Available 6am-5pm MST M-F

**By Online Chat @ [promega.com](https://www.promega.com)**

- Available 6am-5pm MST M-F
- Global Chat with Branch office tech serv scientists, too, after hours (language dependent)

**By email:** [techserv@promega.com](mailto:techserv@promega.com)

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



North American Branch Technical Support Team