

# Academia/Industry Cross-Fertilization through Translational Research



**Franck Madoux, PhD**

HTS/Lead Identification Department

The Scripps Research Institute Molecular Screening Center

Translational Research Institute

Scripps Florida, Jupiter, FL

Before we start...



Scientific thinking, Business thinking

# Cifre

Conventions

CIFREs (conventions industrielles de formation par la recherche) are funded by the French Ministry of Higher Education and Research.

Industrial contracts for training through research



## Students

Boost your future with a diploma-based first job

## Companies

Boost your performance by recruiting a PhD student

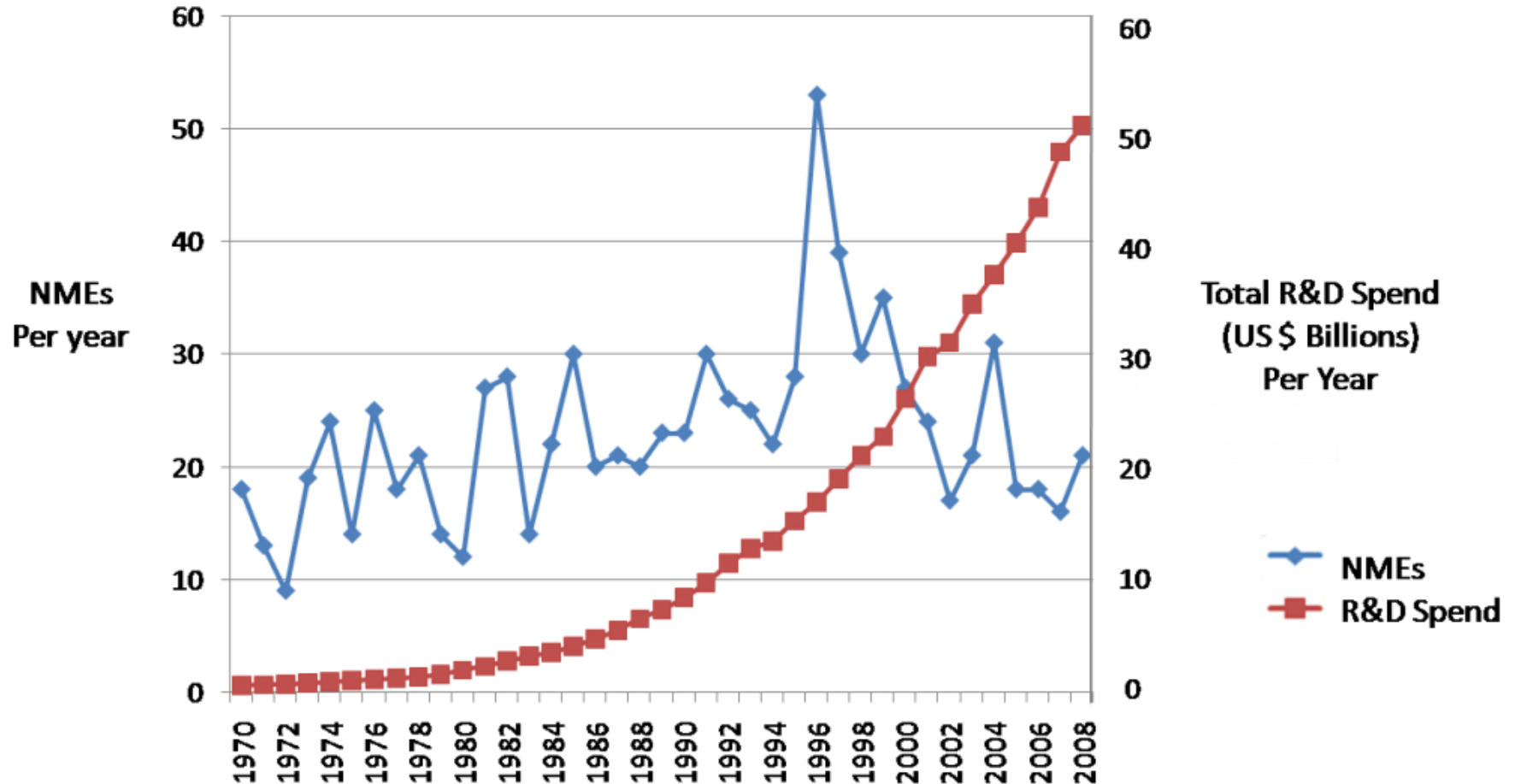
## Laboratories

Boost your research with a CIFRE

→ ***CIFRE industrial fellowship: a WIN-WIN-WIN deal!***

- **Current challenges in the Pharma industry and how to fill the innovation gap.**
- **Translational Research & its role in public-private partnerships**
- **Define the strengths of academia & industry in drug research and development**
- **The Scripps Research Institute's (TSRI) approach to Translational Research**
- **The NIH's Roadmap Initiative and its role in catalyzing translational research in the US**
- **Four examples of Industry/Academia collaborations performed at Scripps boosted through Translational Research**

# Current Challenges in Drug Discovery

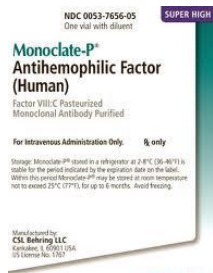


→ *How do we fill the innovation gap?*

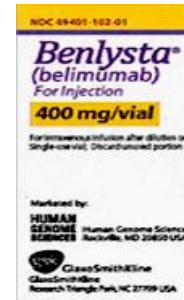
# How do we fill the Innovation Gap?



# Scripps' Industrial Partnerships



CSL Behring



## Research Collaborations



## Technology Licensing

Compound	Therapeutic Area	Company	Phase I	Phase II	Phase III	NDA
<b>Late Stage Products</b>						
Cilengitide	Oncology	Merck Serono	→	→	→	→
ch14.18	Oncology	United Therapeutics	→	→	→	→
<b>Other CAT Products</b>						
Tralokinumab	Respiratory	MedImmune/AstraZeneca	→	→	→	→
CAM-3001	Inflammation	MedImmune/AstraZeneca	→	→	→	→
Rozrolimupab	Cardiovascular	Symphogen	→	→	→	→
GC-1008	Multiple Areas	Genzyme	→	→	→	→
TB-403	Ophthalmology	BioInvent	→	→	→	→
BI-505	Oncology	BioInvent	→	→	→	→
MT-203	Inflammation	Micromet	→	→	→	→
Bertilimumab	Allergy	iCo Therapeutics	→	→	→	→
<b>CovX Products</b>						
CVX-060	Oncology	Pfizer	→	→	→	→
CVX-096	Diabetes	Pfizer	→	→	→	→
<b>Sangamo Products</b>						
SB-728	HIV	Sangamo	→	→	→	→
SB-313	Oncology	Sangamo	→	→	→	→
<b>Click Chemistry Products</b>						
Solithromycin	Antibiotic	Cempra	→	→	→	→
AZ-01	Autoimmune	Allozyne	→	→	→	→
<b>Ambrx Products</b>						
ARX-201	Growth Deficiency	Ambrx	→	→	→	→
ARX-424	Autoimmune	Ambrx	→	→	→	→
<b>Other Early Stage Products</b>						
IC-14	Respiratory	Implicit	→	→	→	→
CTL-04	Oncology	J&J	→	→	→	→
ALT-801	Oncology	Altior	→	→	→	→
RPC-1063	Autoimmune	Receptos	→	→	→	→
Shok-Pak	Organ Failure	InflammaGen	→	→	→	→
RG-2833	Friedreich's Ataxia	Repligen	→	→	→	→
HSC-835	Oncology	Novartis	→	→	→	→
3K3A-APC	Cardiovascular	ZZ Biotech	→	→	→	→

## Startups/Spinouts



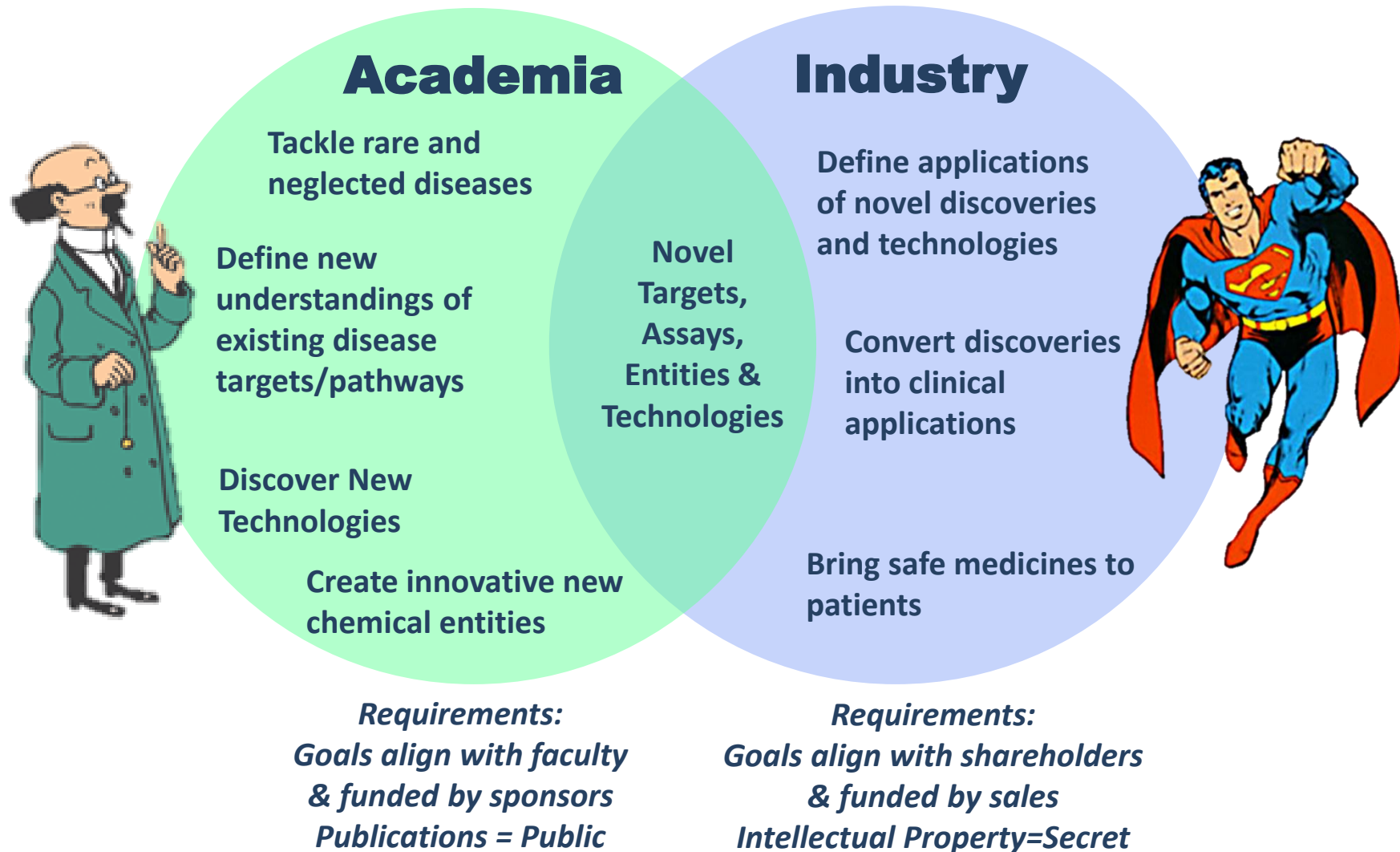
ABIDE THERAPEUTICS



ACHAOPEN



# Defining the strengths of Industry and Academia in Drug R&D



→ **Academia and Industry share common goals**

**How do we make Professor Calculus and Superman reach these common goals together?**

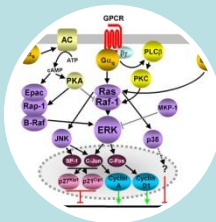


# Scripps Florida's Translational Research Institute





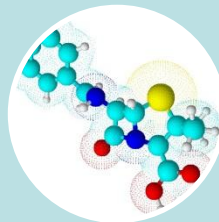
# Scripps Florida's Translational Research Institute



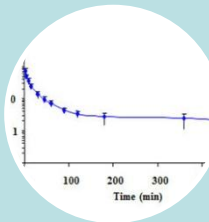
Discovery  
Biology



Lead  
Identification



Medicinal  
Chemistry

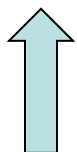


DMPK



Animal  
Studies

Translational Research Institute (TRI)

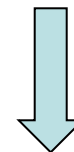


Novel or  
Neglected target



The Translational Research Institute offers:

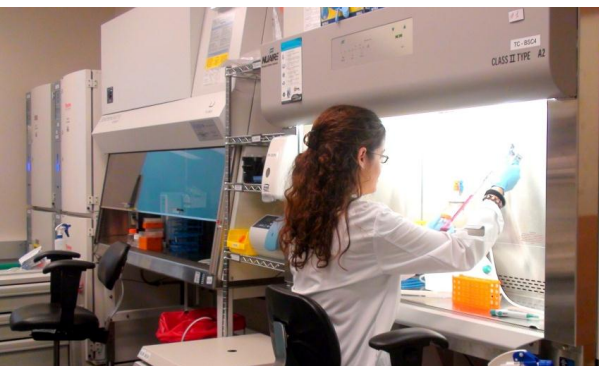
- Biotech-type organization
- A 50/50 blend of faculty from the industry and the academia
- A common language
- Drug Discovery Resources
- Pharma/Biotech expertise
- Advanced equipment and technologies



Drug  
Candidate

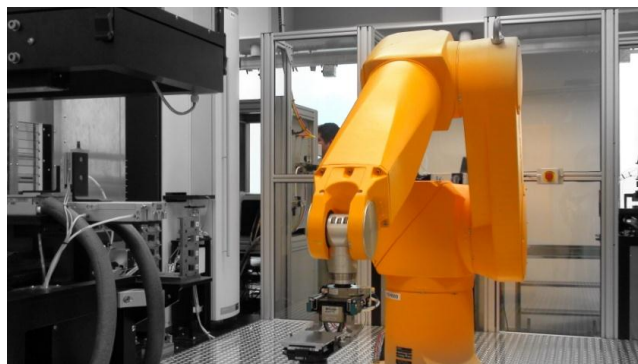


# The uHTS/Lead Identification Group @ Scripps Florida



## Assay Development Lab

“From test tube to plate”  
protein expression/purification  
>60 mammalian cell lines  
insect cells  
bacteria, yeast  
Batch transfection  
Frozen, assay-ready cells  
Small organisms



## GNF Systems uHTS Platform

>250,000 tests /day  
1,536 well format  
Homogenous & Heterogeneous Assays  
FLINT, FP, TR-FRET, Lumi, Abs., FLIPR...  
High Content Screening



## Compound Management

> 1M Screening Repository:  
~640K Proprietary to Scripps  
~370K Public Domain (NIH)  
~100K Private Collaborations  
comprising both small molecules and natural products

- The “Lead ID” facility occupies >6100 ft<sup>2</sup> of laboratory space
- 20 FTEs: cell /molecular biologists, biochemists, microbiologists, software programmers, engineers, cheminformaticists, compound managers
- Access to >50 Medicinal Chemists, DMPK, Pharmacology @ Scripps FL
- Engaged in both private & NIH-funded HTS collaborations

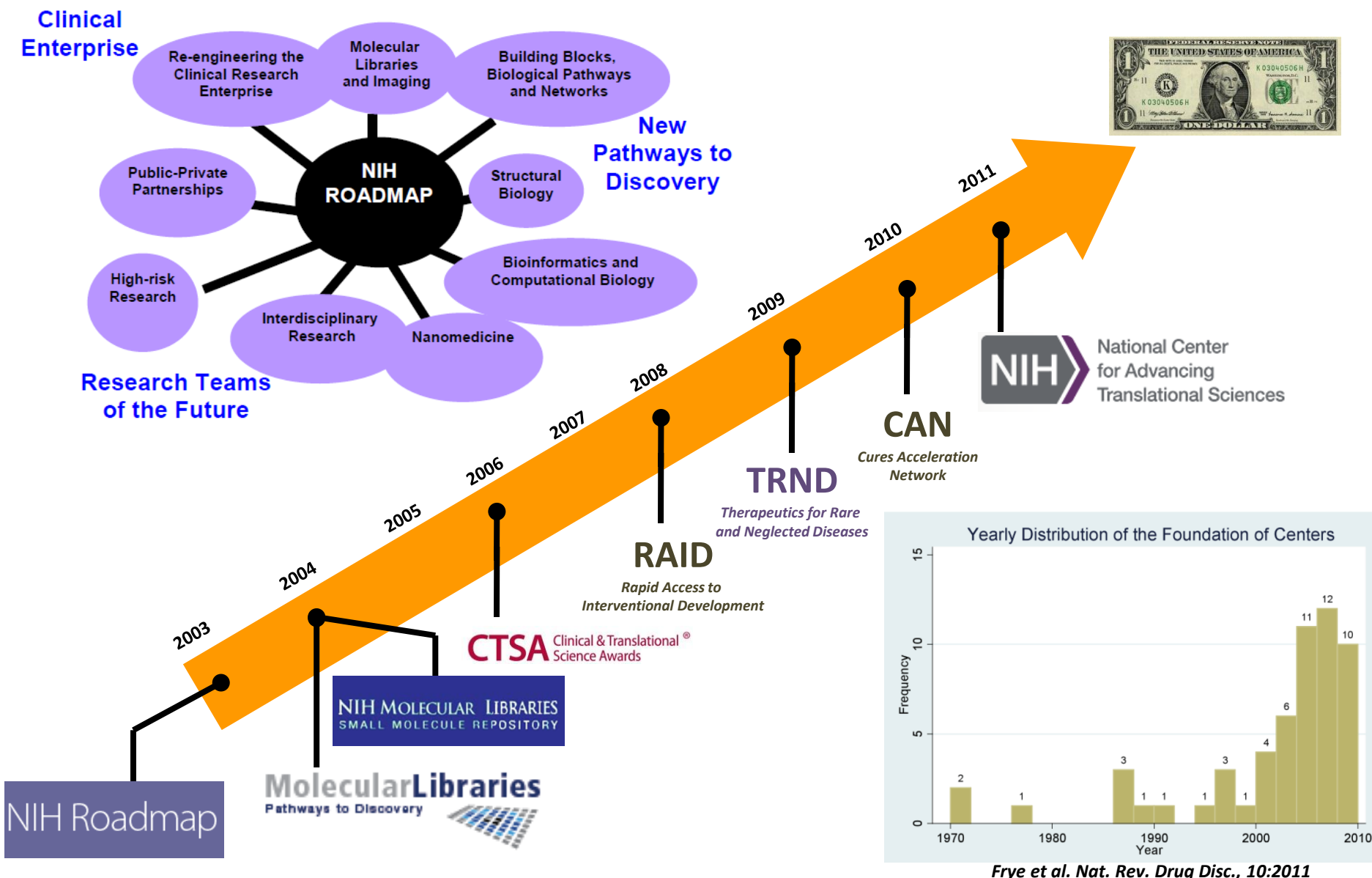
→ *“industrial-grade” expertise and equipment , as well as workflows, timelines, and QC standards makes us credible interlocutors for the pharma industry*

*“It is the responsibility of those of us involved in today’s biomedical research enterprise to translate the remarkable scientific innovations we are witnessing into health gains for the nation”*

**Elias Zerhouni, MD  
Director, NIH  
NEJM 2005**



# The Role of the NIH in catalyzing Translational Research



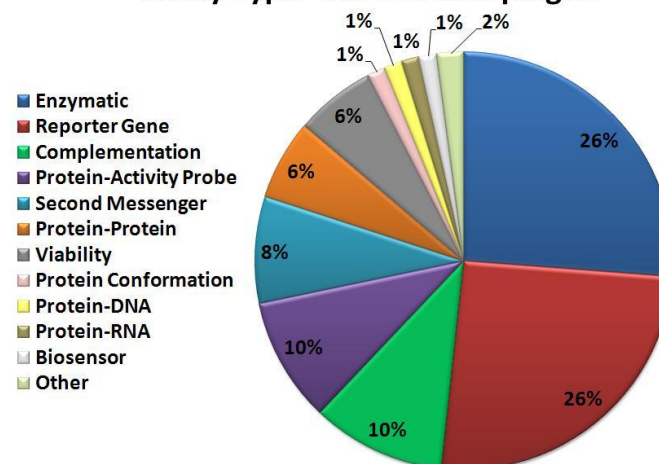
# The Molecular Libraries Probe Production Center Network (MLPCN)

- **Part of NIH's "roadmap initiative" for biomedical research in the 21<sup>st</sup> century**
- **Grants for assay development and/or HTS**
- **Peer-reviewed process**
- **Open to academia, government, non-profit and industry**
- **Screening centers provide assay development, HTS, DMPK and medicinal chemistry support A/R**
- **All results are publicly available through NCBI's PubChem website**
- **Each center has a copy of the same library (360K, still growing)**
- **Production phase:**
  - **10 specialized centers**
  - **4 comprehensive centers 25 targets/year/center**

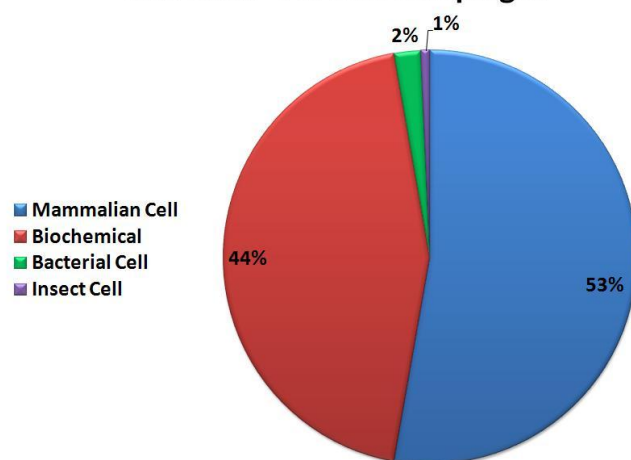
# MLPCN @ Scripps: a retrospective of the past 6 years

- \$89 million over 6 years
- >150 targets
- 71 chemical probes
- >1,300 PubChem reports
- 141 publications
- 4,155 citations

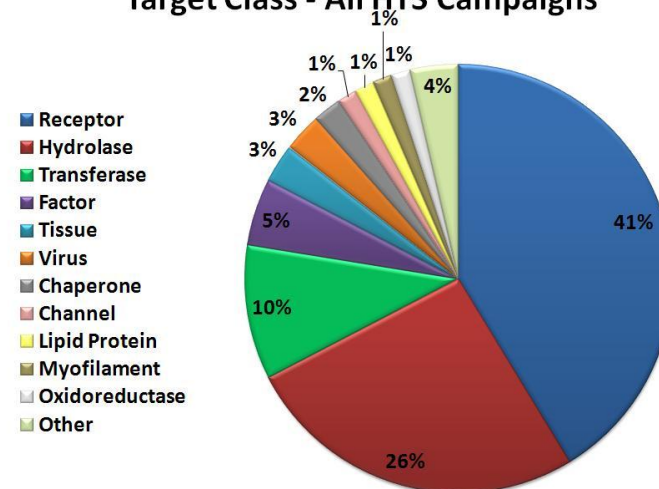
Assay Type - All HTS Campaigns



HTS Class - All HTS Campaigns



Target Class - All HTS Campaigns



**An interactive table that details all our public domain HTS collaborations can be found at:**

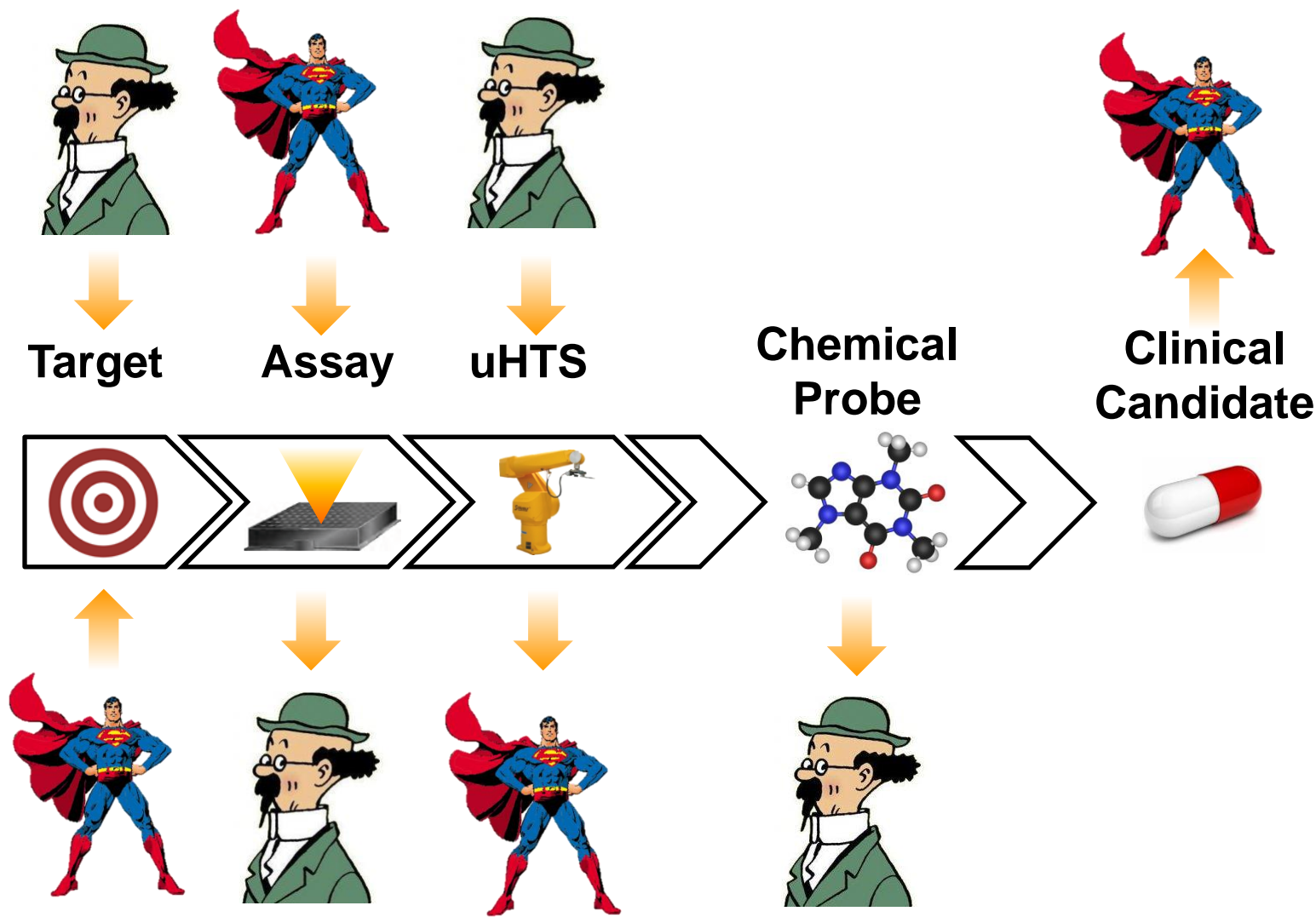
<http://hts.florida.scripps.edu/index.php/our-capabilities/hts-campaigns/pipeline.html>



# 4 examples of Academia/Industry partnerships

Therapeutic area	Target	Target Class	Assay Technology	Academic Partner	Industrial Partner
Infectious Diseases	HCV-CORE	Viral protein	HTRF	Scripps FL Boston University	Pfizer
Cancer	SF-1	NHR	Luciferase Reporter	Scripps FL - CNRS	Orphagen + Biotech 'X'
Autoimmune Disorders	S1P1	GPCR	BLA	Scripps LJ	Receptos
CNS	x	x	x	Scripps FL	Envoy Therapeutics

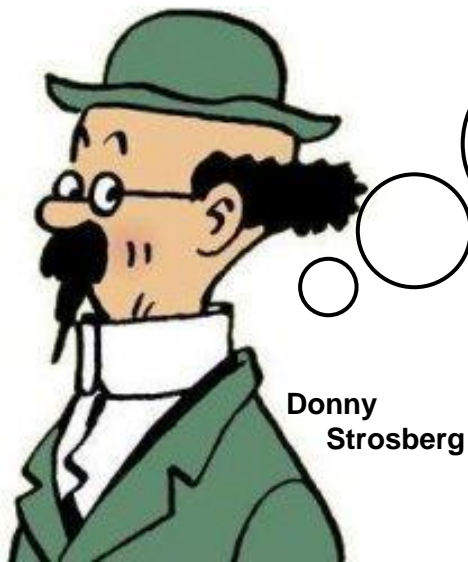
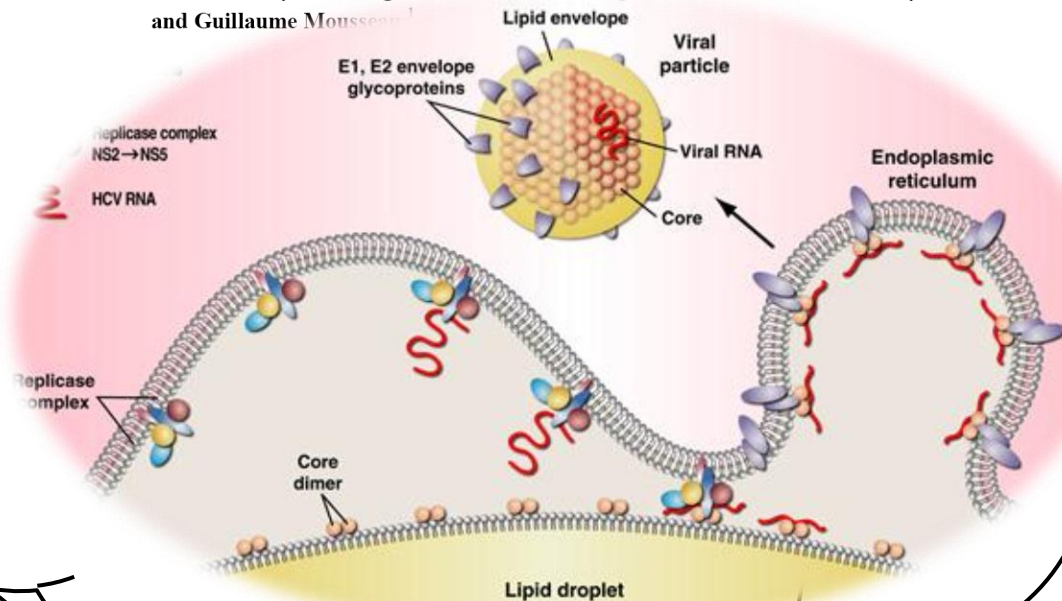
# Academia/Industry Cross-Fertilization



Review

## Core as a Novel Viral Target for Hepatitis C Drugs

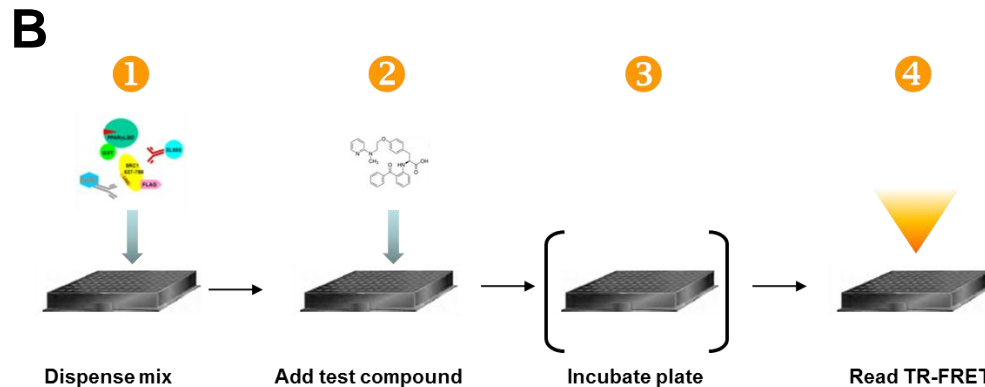
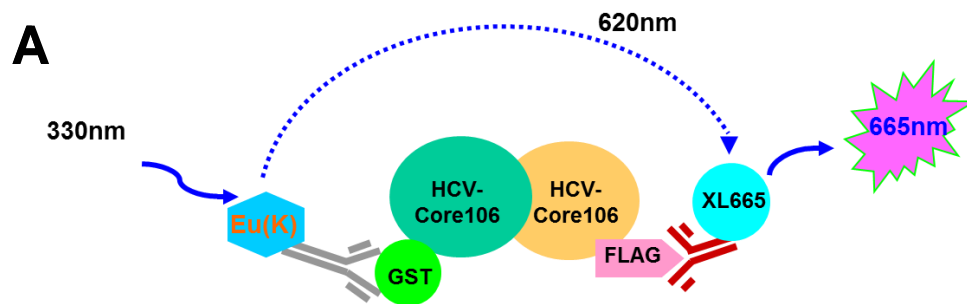
Arthur Donny Strosberg<sup>1,\*</sup>, Smitha Kota<sup>1</sup>, Virginia Takahashi<sup>1</sup>, John K. Snyder<sup>2</sup>  
and Guillaume Mousseau<sup>1</sup>





# A Time-Resolved Fluorescence-Resonance Energy Transfer Assay for Identifying Inhibitors of Hepatitis C Virus Core Dimerization

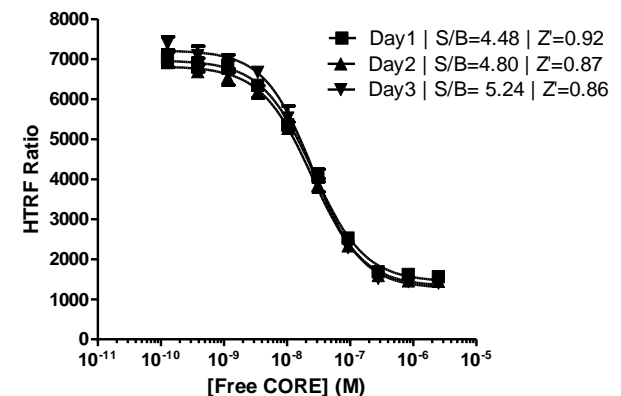
Smitha Kota,<sup>1</sup> Louis Scampavia,<sup>2</sup> Timothy Spicer,<sup>2</sup> Aaron B. Beeler,<sup>3</sup>  
Virginia Takahashi,<sup>1</sup> John K. Snyder,<sup>3</sup> John A. Porco Jr.,<sup>3</sup>  
Peter Hodder,<sup>2</sup> and Arthur Donny Strosberg<sup>1</sup>

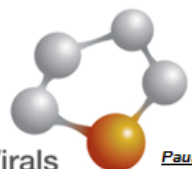


**C**

		EuD01				EuD02				EuD03				EuD04			
		XL665-1	XL665-2	XL665-3	XL665-4	XL665-1	XL665-2	XL665-3	XL665-4	XL665-1	XL665-2	XL665-3	XL665-4	XL665-1	XL665-2	XL665-3	XL665-4
FLAG-1	OST-1	1.93	1.76	1.63	1.39	1.92	1.69	1.62	1.38	1.59	1.40	1.45	1.36	1.33	1.34	1.33	1.35
	OST-2	2.05	1.73	1.63	1.45	1.84	1.73	1.65	1.47	1.44	1.60	1.55	1.43	1.34	1.33	1.33	1.26
	OST-3	1.89	1.70	1.61	1.41	1.81	1.69	1.60	1.46	1.58	1.51	1.56	1.44	1.31	1.33	1.34	1.33
	OST-4	1.70	1.61	1.50	1.40	1.63	1.62	1.60	1.51	1.42	1.43	1.46	1.41	1.29	1.24	1.31	1.29
FLAG-2	OST-1	2.20	1.97	1.78	1.56	2.33	1.95	1.85	1.64	1.82	1.72	1.68	1.55	1.40	1.46	1.42	1.38
	OST-2	2.36	2.08	1.89	1.65	2.36	2.05	2.06	1.72	1.86	1.68	1.82	1.61	1.56	1.52	1.48	1.47
	OST-3	2.33	2.04	1.89	1.65	2.39	2.28	2.00	1.81	1.80	1.82	1.72	1.64	1.61	1.55	1.55	1.50
	OST-4	2.05	1.96	1.83	1.62	1.99	2.05	1.95	1.79	1.74	1.78	2.02	1.66	1.50	1.55	1.53	1.51
FLAG-3	OST-1	2.17	2.03	2.03	1.71	2.23	2.07	1.97	1.75	1.91	1.87	1.77	1.66	1.62	1.56	1.48	1.43
	OST-2	2.51	2.23	2.08	1.83	2.40	2.41	2.22	1.94	2.07	1.99	1.93	1.80	1.75	1.74	1.65	1.53
	OST-3	2.50	2.15	2.16	1.88	2.43	2.40	2.30	2.06	2.04	1.98	2.00	1.93	1.77	1.73	1.74	1.65
	OST-4	2.41	2.24	2.15	1.89	2.42	2.40	2.33	2.32	2.05	2.07	2.00	1.94	1.78	1.78	1.78	1.71
FLAG-4	OST-1	2.03	1.99	1.95	1.74	2.08	2.06	2.05	1.73	1.91	1.94	1.87	1.73	1.61	1.59	1.55	1.49
	OST-2	2.33	2.33	2.20	1.86	2.46	2.44	2.32	2.30	2.38	2.19	2.16	2.02	1.86	1.86	1.81	1.65
	OST-3	2.46	2.41	2.35	2.08	2.62	2.63	2.56	2.27	2.31	2.29	2.28	2.14	1.95	2.06	1.96	1.89
	OST-4	2.50	2.50	2.34	2.09	2.77	2.74	2.69	2.35	2.20	2.27	2.27	2.15	2.07	2.16	2.06	1.97

**D**





## The identification of novel small molecule inhibitors of the hepatitis C virus core protein.

Paul Tarquett-Adams<sup>1</sup>, Jared Milbank<sup>1</sup>, Helen Waller<sup>1</sup>, Francois Bertelli<sup>2</sup>, Smitha Kota<sup>3</sup>, Donny Strosberg<sup>2</sup>, Chris Pickford<sup>1</sup>, Tanya Parkinson<sup>1</sup>, Manos Perros<sup>1</sup>.

**SandwichResearch**

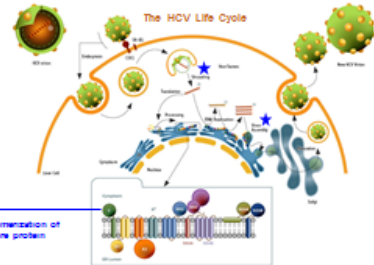


<sup>1</sup> Pfizer Antivirals Research Unit, Pfizer PGRD Laboratories, Sandwich, UK, CT13 9NU; <sup>2</sup>Pfizer HTS Centre of Emphasis, Pfizer PGRD Laboratories, Sandwich, UK, CT13 9NU; <sup>3</sup>The Scripps Research Institute, 130 Scripps Way, Jupiter, Florida 33458, USA.

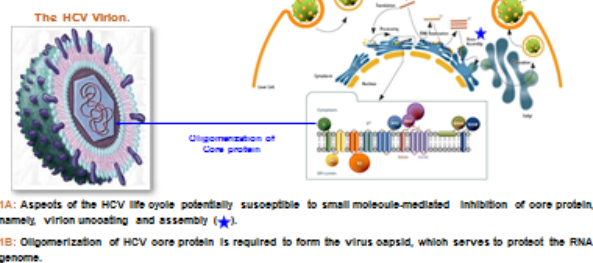
**Summary:** Hepatitis C virus (HCV) is a global health problem. Treatment with interferon and ribavirin is the current standard of care; however, it is not universally effective and is associated with considerable side effects. Consequently, there's an urgent unmet medical need for directed, well-tolerated antiviral medicines to combat the growing burden of HCV infection. The first generation of specific anti-HCV medicines are likely to be small-molecules targeting the virus-encoded enzymatic gene products. However, preliminary *in vitro* and *in vivo* data indicate the emergence of compound-resistant HCV quasi-species are likely to be a considerable complication. To improve our chances of successfully developing new medicines to treat HCV-infected patients, the Antivirals Research Unit is seeking to identify small-molecule inhibitors of core protein; a non-enzymatic HCV gene product. Core protein is the major structural component of the virus and is essential for productive infection; no virus is produced in its absence. Core oligomerizes around the RNA genome of the virus to form a protein shell (capsid), which protects the virus genome. The HCV core protein programme is a result of collaboration with The Scripps Research Institute, where a novel *in vitro* assay for core protein dimerization was developed. The assay was transferred to Pfizer and screened against a compound library composed of >2 million small molecules. Currently, we are progressing through the screening cascade and have identified a panel of compounds, which inhibit dimerization of the core protein *in vitro* and display no adverse effects in cell culture. Preliminary findings demonstrate a subset of this compound panel exhibits antiviral activity in the HCV virus-infectivity assay with low micromolar IC<sub>50</sub>s. These data suggest inhibition of the core-core interaction *in vitro* can translate into antiviral activity in HCV-infected tissue culture cells and provides confidence in targeting core protein oligomerization as a mechanism for potential therapeutic intervention.

### 1. Introduction:

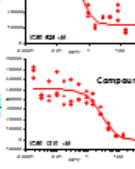
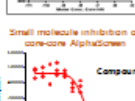
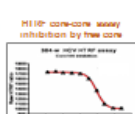
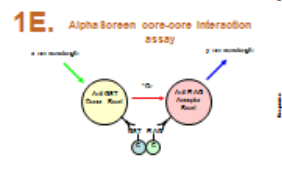
#### 1A.



#### 1B.



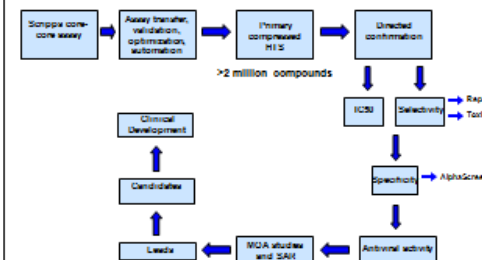
1A: Aspects of the HCV life cycle potentially susceptible to small molecule-mediated inhibition of core protein, namely: virion uncoating and assembly (★).  
1B: Oligomerization of HCV core protein is required to form the virus capsid, which serves to protect the RNA genome.



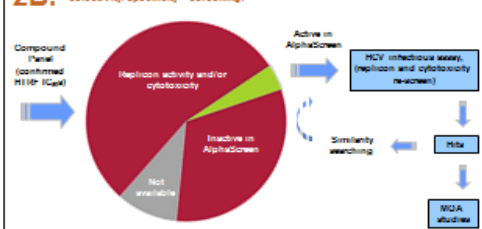
1C: HTRF energy transfer assay designed to recreate the capacity for HCV core protein to self assemble *in vitro*. Assay measures the interaction between GST-core108 and FLAG-core108.  
1D: Inhibition of the core-core HTRF assay by free (untagged) core108 protein (IC<sub>50</sub> 80nM).  
1E: AlphaScreen core-core interaction assay used as a selectivity counter-screen and for follow-up screening of HTS hits.  
1F: Example of inhibition of core-core interaction using the AlphaScreen assay and 2 compounds identified as core-core inhibitors in the HTRF HTS.

### 2. Screening:

#### 2A. Screening Cascade:



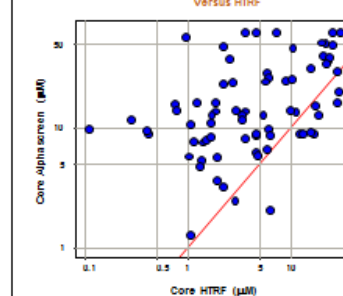
#### 2B. Selectivity/Specificity Screening:



2B: 60% of the compound panel, which was composed of confirmed hits from the *In vitro* core-core HTRF interaction assay, exhibited cytotoxicity or off-target effects as judged by activities in cytotoxicity and/or HCV subgenomic replicon assays. 6% of the compound panel were active in the AlphaScreen selectivity assay and displayed no activity in the cell-based assays; these compounds were progressed to the HCVoo assay to identify molecules with antiviral activity. Similarity searching was used to identify series of compounds with antiviral activity.

### 3. Follow-up:

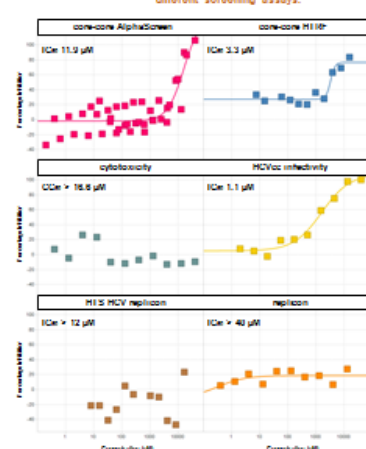
#### 3A. Core-Core Interaction assays: AlphaScreen versus HTRF



3A: Approximately 86% of compounds, confirmed as hits in the HTS HTRF core-core interaction assay, were also active in the AlphaScreen core-core selectivity screen. The scatter plot above describes the activity of a selection of compounds. In both assays, generally compounds appeared more active in the HTRF assay compared to the AlphaScreen assay; a feature presumably related to the different assay technologies.

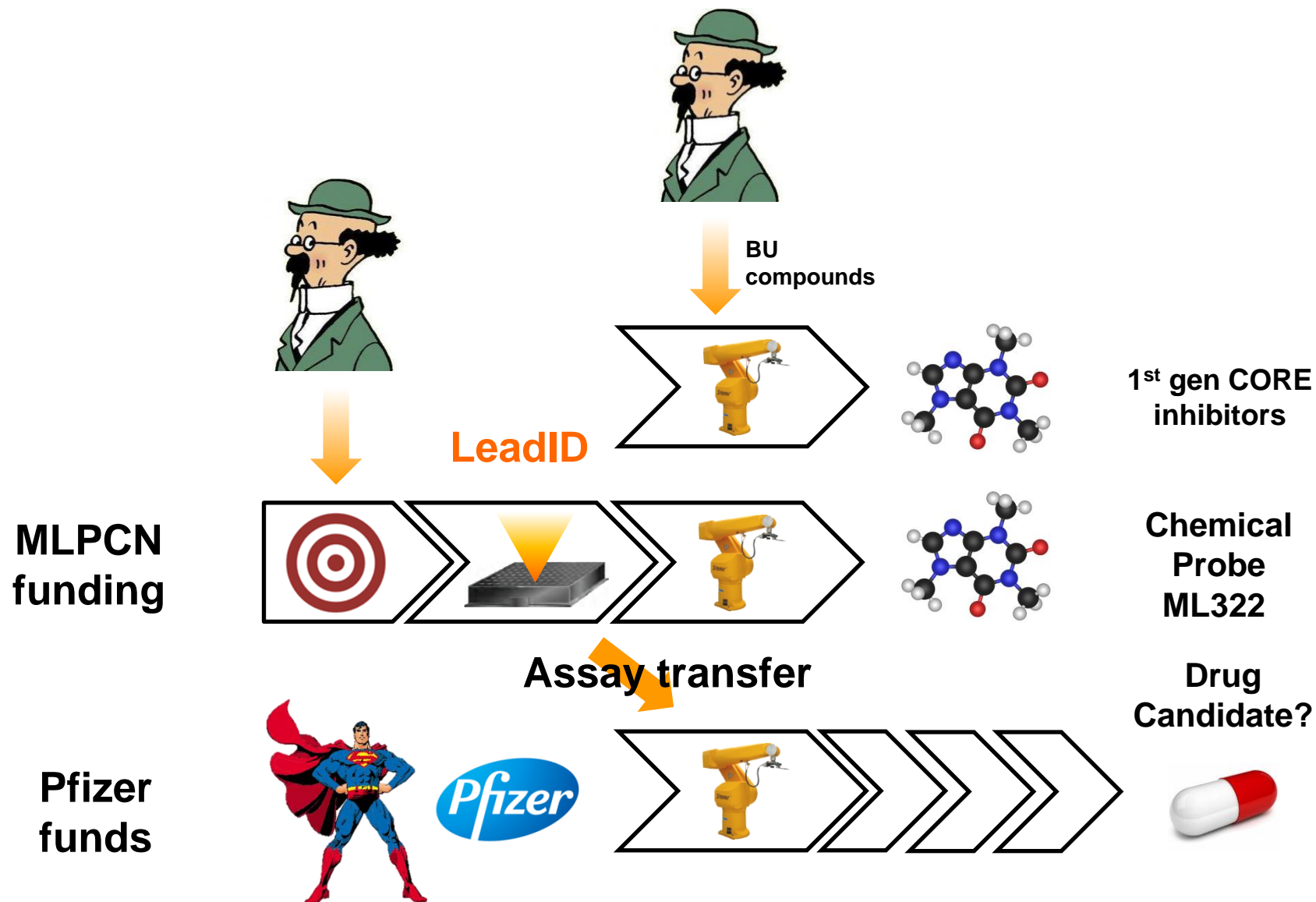
3B: Assay profiles of a representative anti-HCV core compound: Active in both *In vitro* core-core interaction assays (AlphaScreen and HTRF; IC<sub>50</sub>s of 11.8 and 3.3 µM respectively), no apparent cellular toxicity (CC<sub>50</sub> of over 18.8 µM), no off-target effects on HCV RNA replication (inactive in HCV subgenomic replicon assays), and active in the JFH-1 HCVoo infectivity assay (IC<sub>50</sub> of 1.1 µM).

#### 3B. Activity of a representative HTS hit in the different screening assays.



### 4. Conclusions:

Oligomerization of the HCV core protein has potential as a new target for the development of future therapies directed against HCV. Consequently, we have screened the Pfizer compound library for small molecule inhibitors of the core-core interaction and have identified compounds that inhibit the interaction *in vitro* as evidenced by activity in the primary HTRF assay, and the AlphaScreen selectivity assay. To rule out any off-target effects or cellular toxicity, the core actives were screened in replicon and WT-1 toxicity assays. 'Core-targeting' compounds, inactive in the cell-based selectivity screens were tested for antiviral activity in the HCVoo assay. A subset of the core actives inhibited HCV infectivity with low micromolar IC<sub>50</sub>s. Thus, inhibitors of the *In vitro* core-core interaction assay have the capacity to inhibit HCV infectivity in tissue culture cells, providing confidence that oligomerization of core protein represents a mechanism amenable to drug targeting.

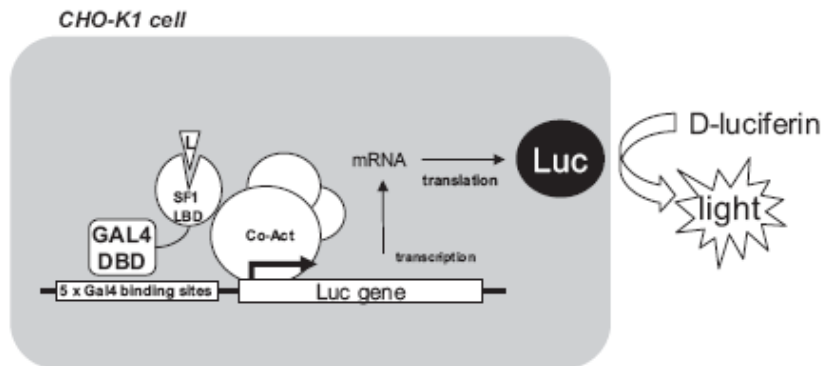


- **RO1 screening proposal submitted by Orphagen Pharmaceuticals**
- **SF-1 belongs to the nuclear hormone receptor (NHR) superfamily (druggable)**
- **Functional ligands are still unknown = orphan receptor**
- **SF-1 expression levels are higher in patients with Childhood Adrenocortical Tumors (ACTs)**
- **ACTs constitute a rare, yet very aggressive and poorly understood type of pediatric cancer**

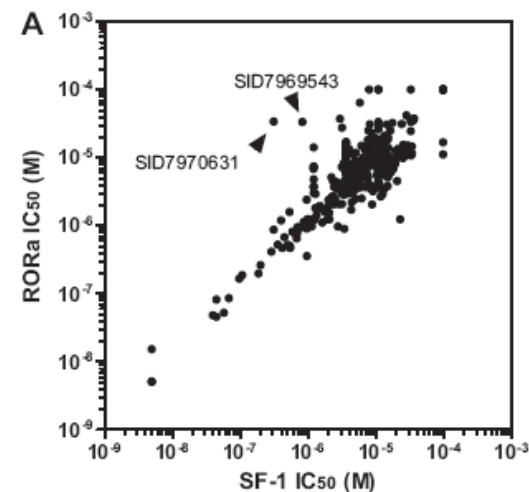


# Example #2: SF-1 | Lead Identification

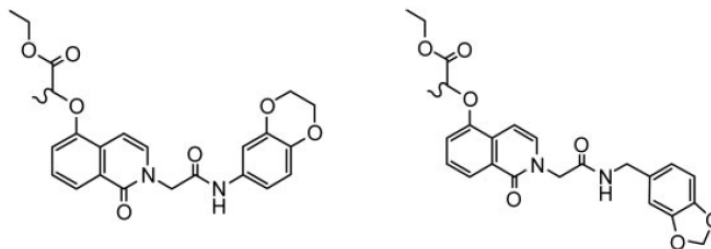
**A**



**B**



**C**



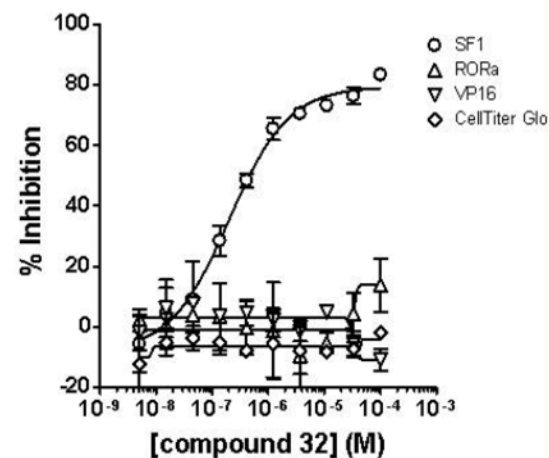
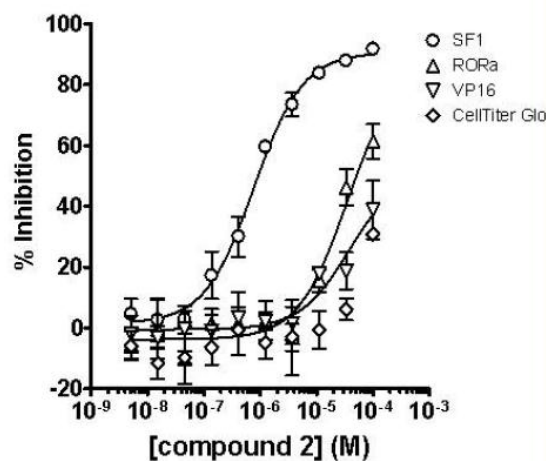
	SID7969543		SID7970631	
	Inhibition (Cytotoxicity)	IC <sub>50</sub> (CC <sub>50</sub> )	Inhibition (Cytotoxicity)	IC <sub>50</sub> (CC <sub>50</sub> )
	%	nM	%	nM
Gal4-fusion assays				
SF-1	84 ± 2	760 ± 102	80 ± 2	255 ± 63
ROR-α	16 ± 4	>33,333	20 ± 13	>33,333
VP16	18 ± 2	>33,333	-48 ± 21	>33,333
SFRE promoter assays with full-length proteins				
SF-1	136 ± 4	30 ± 15	126 ± 3	16 ± 7
LRH-1	0 ± 3	N.A.	-9 ± 9	N.A.
Cytotoxicity assay				
Cytotoxicity	-1 ± 6	>99,000	19 ± 1	>33,333

N.A., not applicable, because compound did not reach 50% inhibition; N.T., not tested.

**→ a transient transfection cell-based screen led to the identification of 2 selective compounds**

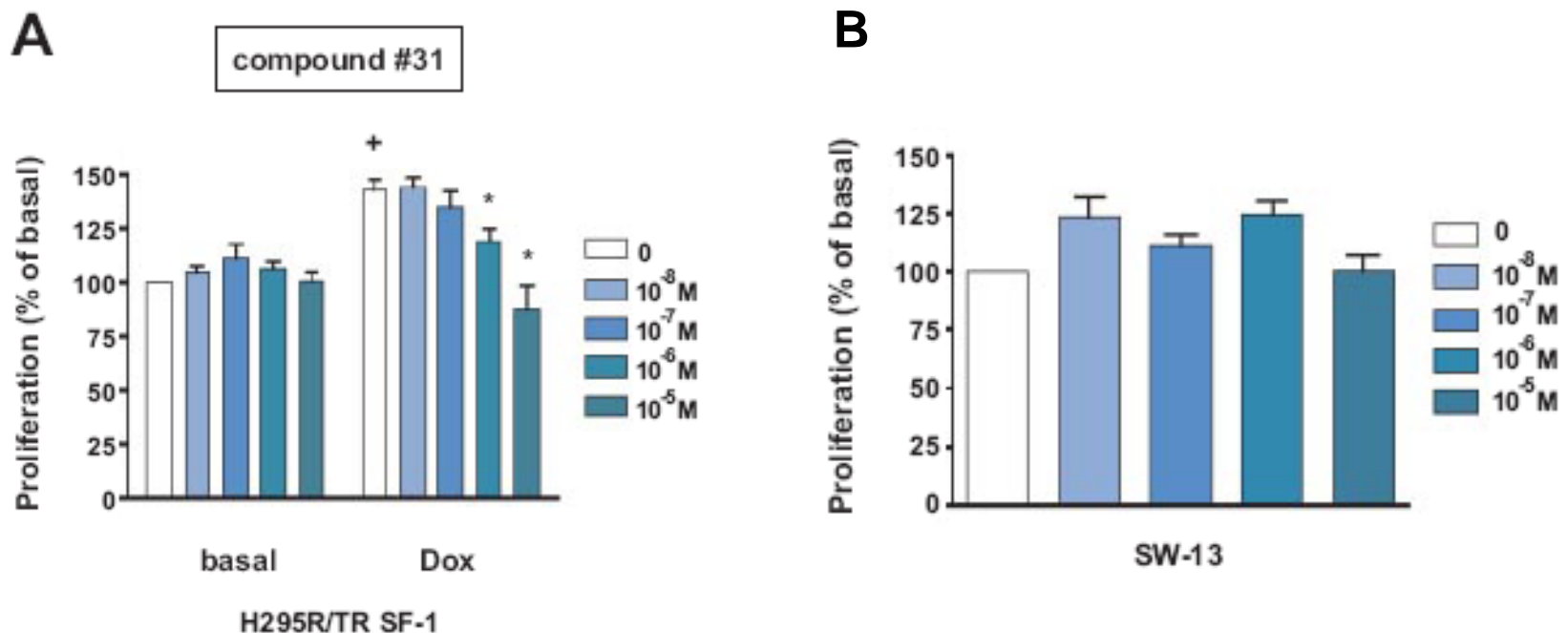
**Madoux et al., Mol Pharm 2008**

## Example #2: SF-1 | Lead Optimization



Roth et al., Bioorg. Med. Chem. Lett. 2008

**→ a SAR-based chemistry optimization effort based on  $\approx 50$  analogs led to the generation of compounds exhibiting improved potency and a cleaner activity profile in the counterscreen assays**



Doghman et al. 2009

→ Compounds of the Isoquinolinone family inhibit adrenocortical cell proliferation in a SF-1 dependent



- Target
- Assays



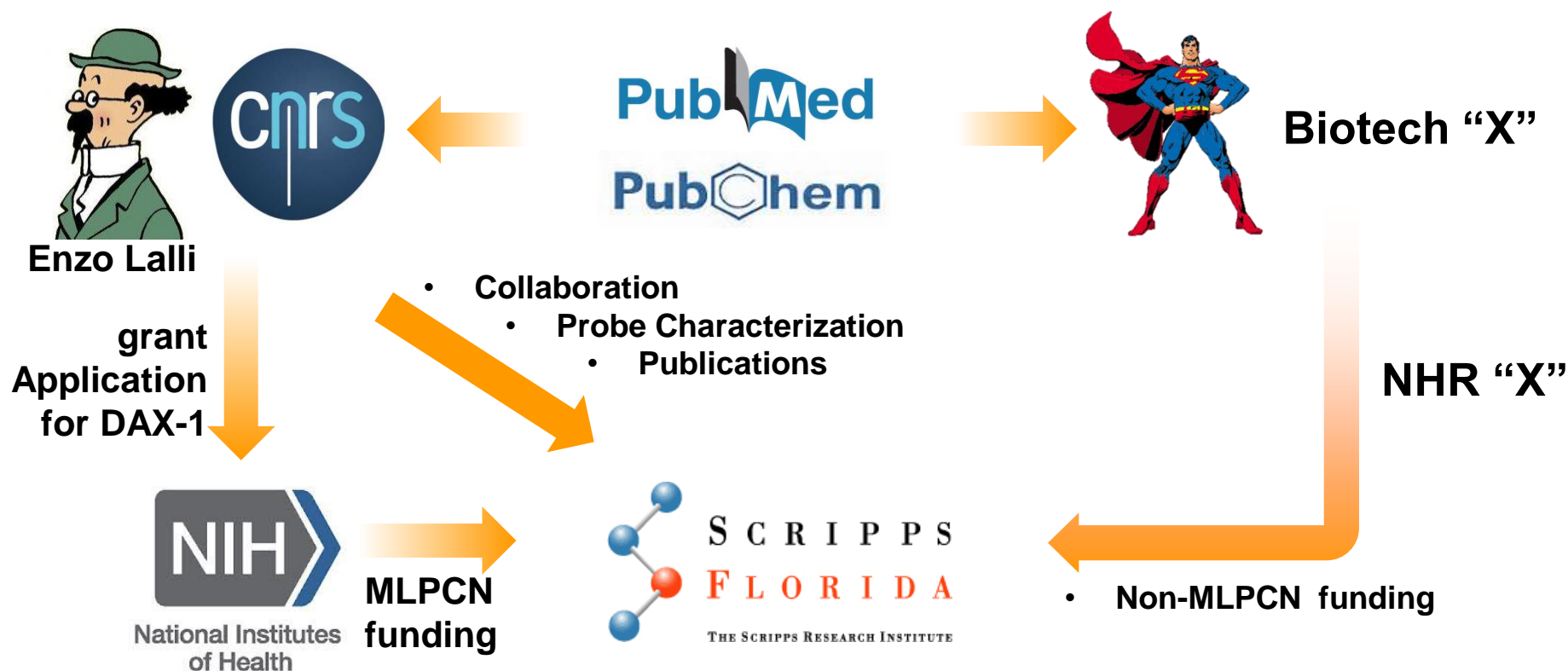
- Proof of concept
- Titration Results (1-year embargo)
- Chemical Probe
- Publications

- Screening Results
- Probe Report

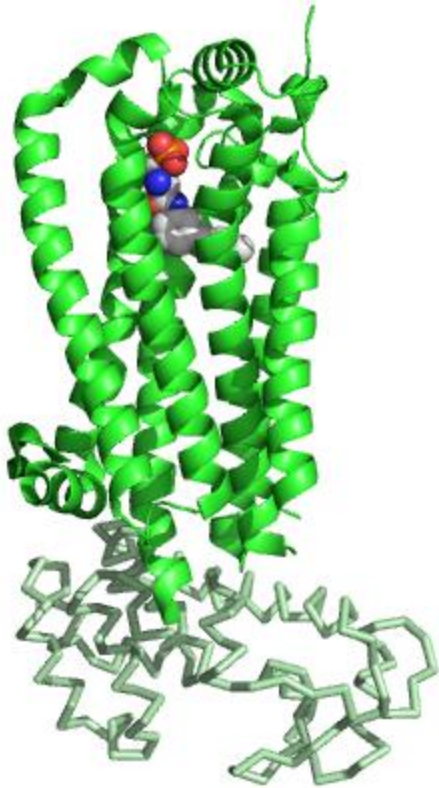




## Example #2: initiation of a virtuous cycle

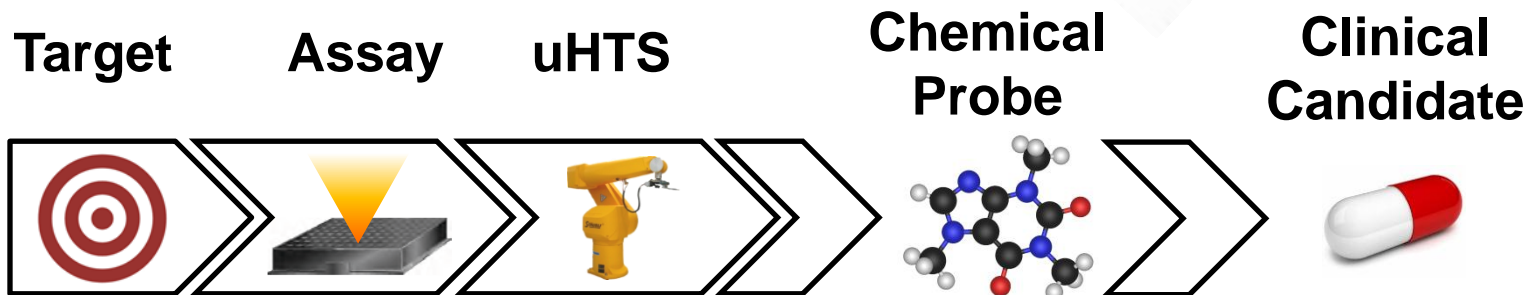
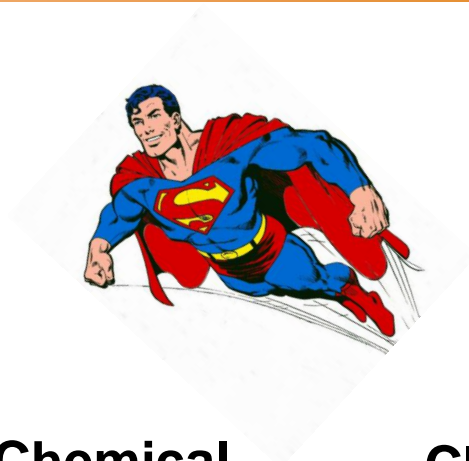


## Example #3: S1P1 modulator



- The sphingosine-1-phosphate (S1P)-driven signaling regulates fundamental biological functions, including cell proliferation, angiogenesis, endothelial cell chemotaxis, immune cell trafficking and mitogenesis.
- S1P1 receptor (S1P1-R) agonists can be used as immunosuppressants.
- S1P1R agonists can be used to treat autoimmune disorders such as relapsing–remitting multiple sclerosis

# Example #3: S1P1 modulator



**Hugh Rosen**



receptos

## Receptos Doses First Patient in Phase 2/3 Trial for RPC1063 in Multiple Sclerosis

SAN DIEGO, CA, October 22, 2012 – Receptos, Inc., announced today that its selective sphingosine-1-phosphate receptor 1 (S1P1) modulator, RPC1063, has been administered to the first patient in a Phase 2/3 study. RPC01-201, a Phase 2/3 placebo-controlled (Phase 2) and active comparator-controlled (Phase 3) trial, is the first of two planned pivotal studies for RPC1063 in the indication of relapsing multiple sclerosis (RMS).

## Example #4: Collaboration with Envoy



**Created a portfolio of CNS targets using their proprietary bacTRAP technology**

### **Envoy and Scripps Add Three Drug Screening Programs**

---

Envoy Therapeutics, Inc., a recently-formed drug discovery company, today announced that it has expanded its research collaboration with The Scripps Research Institute aimed at identifying new drugs for neurological and psychiatric diseases. Augmenting the collaboration Envoy and Scripps commenced in July focused on improved treatments for Parkinson's disease, scientists at the two organizations will carry out three additional drug discovery programs over the coming year. These programs will employ Scripps-Florida's high-throughput screening capabilities to discover compounds that modulate target proteins identified by Envoy.

# Example #4: Collaboration with Envoy



## Takeda buys Envoy for \$140m

7 November 2012

Andrew Turley



8



5



0

Japanese drug maker [Takeda](#) has struck a \$140 million (£88 million) deal to buy privately owned US drug discovery company Envoy.

The move gives Takeda access to Envoy's 'BacTrap' technology for labeling and extracting the protein-making components of specific types of cells. The technique is particularly useful in the brain, so its acquisition will strengthen Takeda's potential in the central nervous system drugs market.

Takeda plans to continue operating Envoy in Jupiter, Florida, until March 2013 and then transfer the majority of the Envoy scientific staff and management team to Takeda in San Diego, California.

Envoy was established in 2009 and now employs 21 people. Takeda invested in the company in October 2009 through its corporate investment wing, Takeda Ventures.



# Example #4: Collaboration with Envoy

Mar 13, 2013

## Takeda Expands Envoy-Scripps Collaboration

The [Scripps](#) Research Institute (TSRI) and the [Takeda](#) Pharmaceutical Company are expanding an initial collaboration launched in 2010 between scientists on the Florida campus of TSRI and [Envoy](#) Therapeutics, one that reportedly led to several breakthroughs in identifying potential new compounds for neurological and psychiatric diseases, to search for new drug targets for a variety of diseases. [Envoy was acquired by Takeda Pharmaceuticals last November](#) to obtain access to Envoy's CNS drug pipeline.

"We originally came to Jupiter because of Scripps Florida and are thrilled that the potential of our original collaboration has been realized," Stephen Hitchcock, svp, drug discovery at Envoy said. "Now we're moving into new therapeutic areas with different biological targets. The first step is to find small molecules that can validate those targets—and Scripps Florida is amongst the very best places to do that."

Takeda will be utilizing Scripps Florida's [high-throughput screening](#) facility, which is part of its larger translational research infrastructure. The facility reportedly has expertise in transforming slow, labor-intensive biological and biochemical bench-top experiments into high-throughput [screening](#) experiments ("screens"). Fully automated robotic screening platforms can then test more than 650,000 drug-like compounds for pharmacologic activity. After completion of the screens, the facility uses other technologies to support the development of clinically relevant compounds.

Takeda isn't just working with Scripps to develop new drugs: in February, they announced a partnership with Resolve Therapeutics to develop compounds for the treatment of lupus and other autoimmune diseases. Takeda will help fund continued development of Resolve's lead compound RSLV-132 through an initial payment of \$8 million to Resolve and will pay Resolve an option exercise fee, plus the potential for additional development milestones totaling \$247 million. Resolve is also eligible to receive royalties on product sales.



- **Academia and Industry have complementary talents in Drug Discovery**
- **Ensuring parties speak the same language, align their goals and understand their differences are key to a successful collaboration**
- **A dedicated infrastructure and personnel (Translational Research Institute) helps interfacing between industry and academia**
- **A vision/ignition spark is needed: the US federal government (NIH) has played an integral role in the creation and implementation of translational research in academia in general and at Scripps in particular**
- **Successful collaborations exist in a variety of different forms. Be creative!**
- **Innovative targets, risky strategies, as well as rare and neglected diseases are excellent starting points for a successful collaboration**
- **Discoveries from either the Industry or Academia can fertilize the other, and vice-versa (virtuous cycle)**



**Founder/Co-Founder of several biotechnology companies including:**

- **HepCCo LLC.**
- **BioRelix (2005)**
- **Eventus DX (Lab Discoveries) (2005)**
- **Ocean Ridge BioSciences (2005)**
- **Symansis (2002)**
- **Axxima (1998)**
- **Hybrigenics (1998)**
- **Small Molecule Therapeutics (1997)**
- **Peptide ImmunoLigands (1996)**
- **Neurotech SA (1995)**
- **Pharmaceutical Peptides Inc./Praecis (1994)**
- **Vetigen (1992)**
- **Ideon/Incyte Inc. (1989/1991)**
- **AES-Chemunex (Chemunex SA) (1984)**



# Acknowledgments



## Translational Research Institute

Pat Griffin  
Mike Cameron  
Bill Roush  
Tom Bannister  
Patsy McDonalds  
Derek Ducket  
and many more!



National Institutes  
of Health

1 X01 MH079861-01  
U54-MH074404  
U54-MH084512

## Lead ID Team

Joe Datsko  
Christina Eberhart  
Pierre Baillargeon  
Peter Chase  
Lina DeLuca  
Laura Pedro-Rosa  
Katharine Emery  
Virneliz Fernandez-Vega  
Emery Smith  
Ross Ensteinder  
Louis Scampavia  
Tim Spicer  
**Peter Hodder**

## SF-1 project

Xiaolin Li  
Scott Thacher  
Enzo Lalli  
Mabrouka Doghman  
Julie Cazareth  
Dominique Douguet  
Joshua Roth

## HCV-CORE project

Smitha Kota  
Virginie Takahashi  
Guillaume Mousseau  
Donny Strosberg

## S1P1 project

Stephan Shurer  
Hugh Rosen  
Steve Brown  
Jackie Chapman  
MT Schaeffer



Frédérique Hudon  
François Degorce  
Suzanne Graham  
Ryan Boyle  
Estelle N'Garwate  
Patrice Robinson  
Glenn Knapik



→ Visit our website @ <http://hts.florida.scripps.edu/>