

Loison, S.<sup>1</sup>; Cottet, M.<sup>2</sup>; Orcel, H.<sup>2</sup>; Lamarque, L.<sup>3</sup>; Trinquet, E.<sup>3</sup>; Dupuis, E.; Mouillac, B.<sup>2</sup>; Hibert, M.<sup>1</sup>; Durroux, T.<sup>2</sup>; **Bonnet, D.<sup>1</sup>**

<sup>1</sup> Laboratoire d'Innovation Thérapeutique, UMR 7200, Faculté de Pharmacie, 74 route du Rhin, 67401 Illkirch, France;

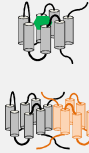
<sup>2</sup> Institut de Génomique Fonctionnelle, CNRS UMR 5203, INSERM U661, Université Montpellier I et II, 141 rue de la Cardonille, 34094 Montpellier

<sup>3</sup> Cisbio Bioassays, 30200 Codolet.

## Objectives

### G Protein-Coupled Receptor

- Site of action of about 30% of all drugs on the market.
- Drugs target only 60 GPCR (over 390 non-olfactory GPCR): potential to exploit the remaining family members.<sup>1</sup>



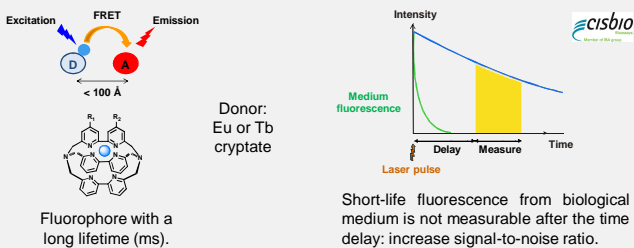
- 1) Accelerate GPCR ligands discovery: new receptor-selective HTS assays.
- 2) Gain a better understanding of their functioning and molecular structure.<sup>2</sup>

(1) Congreve et al. *Trends Pharmacol. Sci.* **2012**, *33*, 249-260; (2) Durroux, T. et al. *Nat. Chem. Biol.* **2010**, *6*, 587-594.

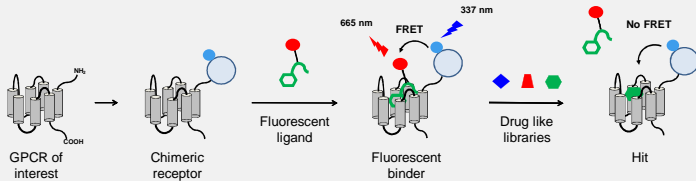
## Strategies

### Time Resolved-FRET based assay (principle)

TR-FRET: combines standard FRET with the time resolved measurement of fluorescence



### Application to GPCR for hit discovery

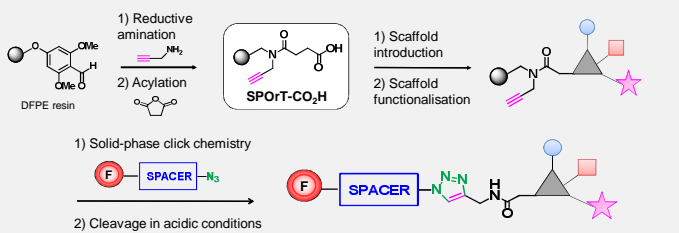


TR-FRET: high sensitivity and reduced environmental safety risk/radioactivity

Challenge: design and synthesis of high affinity fluorescent GPCR ligands

### Novel Solid-phase Organic Tagging (SPOT) resins

Facilitate the labelling of peptides and small organic compounds with fluorescent dyes



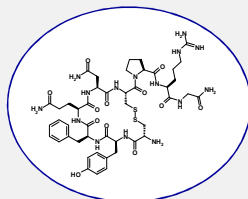
"SPOT" resin: rapid and high yielding access to fluorescent probes, readily amenable to parallel synthesis.

Bonnet, D. et al. *Chem. Eur. J.* **2008**, *14*, 6247-6254.

## Application to Vasopressin V<sub>2</sub> GPCR

### Arginine-vasopressin (AVP) hormone

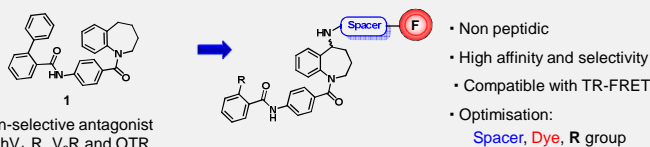
Cardiovascular control, osmoregulation and water homeostasis



### Human AVP V<sub>2</sub> receptor (hV<sub>2</sub>R)

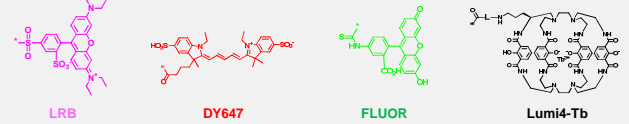
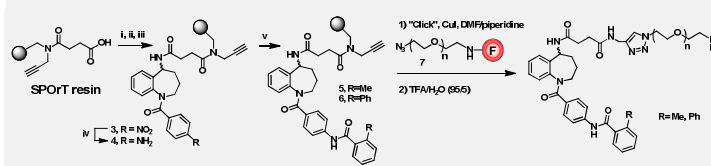
Potential target for the treatment of congestive heart failure, hypertension, or congenital nephrogenic diabetes insipidus

### Design of non peptidic fluorescent probes for hV<sub>2</sub>R



Non-selective antagonist for hV<sub>1a</sub>R, V<sub>2</sub>R and OTR (*Chem. Pharm. Bull.* **1997**, *45*, 1870-1874)

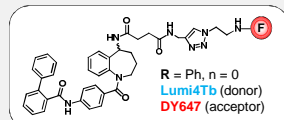
## V<sub>2</sub>R fluorescent probes synthesis



## Biological evaluation

Cpd	R	n	Binding (K <sub>i</sub> , nM)			
			V <sub>2</sub>	V <sub>1a</sub>	V <sub>1b</sub>	OT
AVP	-	-	1.48 ± 0.08	0.7 ± 0.17	0.49 ± 0.06	1.65 ± 0.49
1	Ph	-	0.57 ± 0.18	2.74 ± 0.19	nd	109 ± 34
8	Me	3	54.3 ± 6.6	>10000	>10000	495 ± 101
9	Ph	3	4.0 ± 0.7	>10000	>10000	>10000
40	Ph	3	26.1 ± 5.2	nd	nd	nd
44	Ph	0	5.69 ± 1.05	621 ± 65	>10000	184 ± 31
45	Ph	3	9.02 ± 0.24	1753 ± 338	>10000	808 ± 61
46	Ph	6	14.22 ± 0.19	nd	nd	nd
47	Ph	0	5.86 ± 0.83	180 ± 34	>10000	155 ± 8
48	Ph	3	16.1 ± 2.5	750 ± 89	>10000	304 ± 70
49	Ph	6	30.65 ± 5.03	nd	nd	nd

V <sub>2</sub> R affinity	Selectivity
R	Ph>Me
F	LRB>DY647~
	Lumi4Tb>Fluor
n	0>3>6



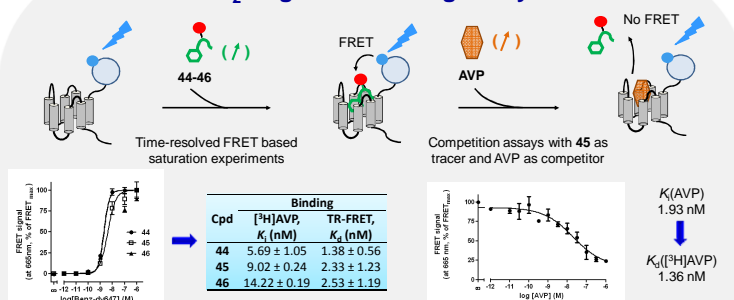
LRB: not compatible with TR-FRET technics

Lumi<sub>4</sub>Tb (donor)  
K<sub>i</sub> = 5.86 nM (hV<sub>2</sub>R)  
V<sub>1a</sub>R/V<sub>2</sub>R selectivity = 31 antagonist

DY647 (acceptor)  
K<sub>i</sub> = 5.69 nM (hV<sub>2</sub>R)  
V<sub>1a</sub>R/V<sub>2</sub>R selectivity = 109 antagonist

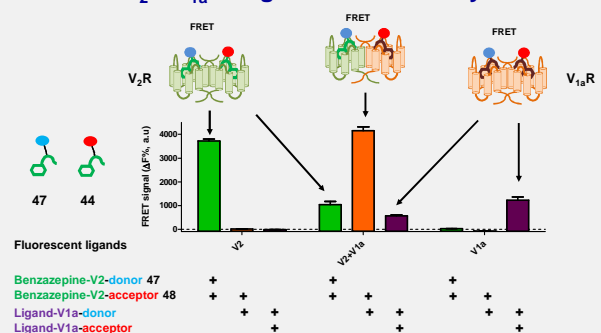
First selective fluorescent non peptidic antagonists for hV<sub>2</sub>R

## V<sub>2</sub>R ligand screening assay



Validation: K<sub>i</sub> and K<sub>d</sub> similar to those determined by radioactive assay  
New V<sub>2</sub>R TR-FRET based assay readily amenable to HTS

## V<sub>2</sub>R-V<sub>1a</sub>R oligomerization assay



V<sub>2</sub>R-V<sub>1a</sub>R heterodimerization at the cell surface

Perspective: studies in native tissues