



# **Evaluation of a Novel Biogenic Amine Transporter Assay.**

**Philip G. Jones  
Discovery Pharmacology,  
Sepracor.**

# A Novel Transporter Assay

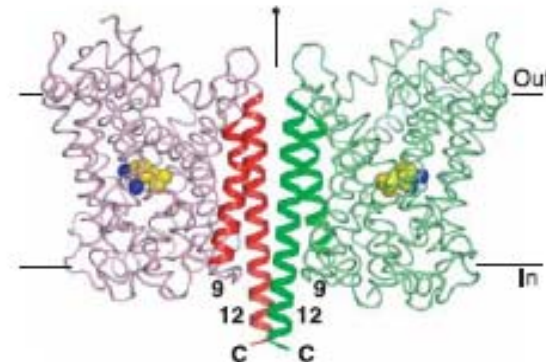
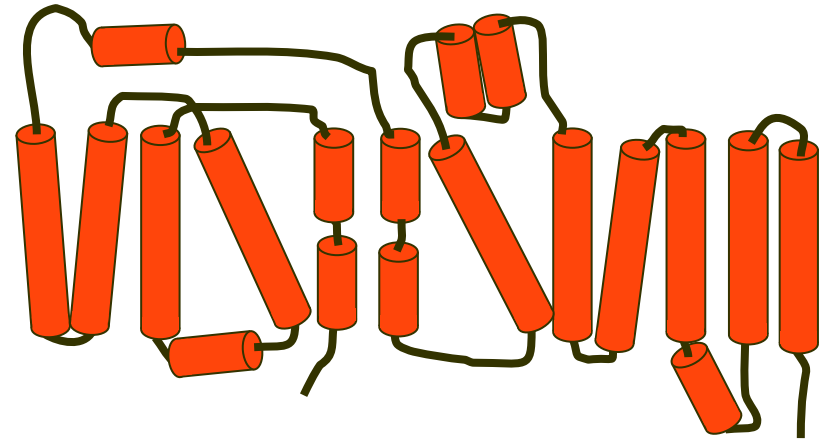


- **Introduction**
  - Neurotransmitter transporters
  - Assays
  
- **Molecular Devices' Fluorescent Assay**
  - Overview
  - Validation
  - Summary

# Neurotransmitter Transporters



- **Transporters**
  - Vesicular
  - Plasma Membrane
- **SLC6 family**
  - Na<sup>+</sup> Cl<sup>-</sup> coupled transporters
  - Dopamine, 5HT, norepinephrine, glycine, GABA
  - Reduce synaptic concentrations of neurotransmitters
  - Terminate Signaling
  - Important targets for pharmaceuticals and drugs of abuse



Taken from Yamashita et al (2005) Nature 437 215-233

# Pharmacology of Biogenic Amine Transporters



<b>Transporter</b>	<b>Endogenous Substrate(s)</b>	<b>Synthetic Substrates</b>	<b>Potent Inhibitors</b>	<b>Therapeutic Use/Potential</b>
<b>DAT</b>	<b>DA, NE,EPI</b>	<b>Amphetamines MPP+</b>	<b>Win35428, mazindol, nomifensine, cocaine</b>	<b>ADHD, Depression, PD, Tourettes, Schizophrenia</b>
<b>NET</b>	<b>DA, NE,EPI</b>	<b>Amphetamines MPP+</b>	<b>nisoxetine, nortripyline, desipramine, duloxetine</b>	<b>Depression, Pain</b>
<b>SERT</b>	<b>5HT</b>	<b>MDMA</b>	<b>sertraline, fluoxetine, duloxetine citalopram, imipramine,</b>	<b>Depression, Anxiety, OCD, Panic, Pain</b>

# Assay Requirements

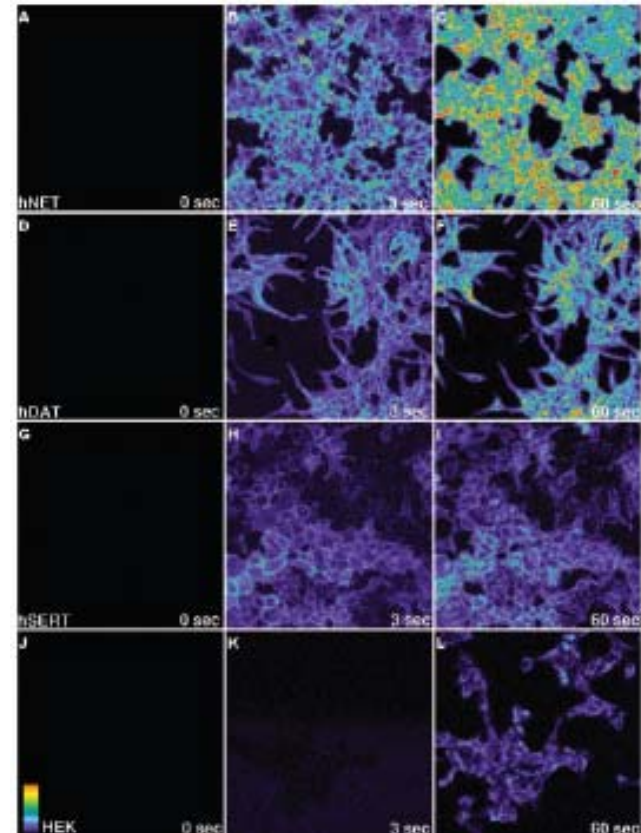


- **High-throughput**
- **Sensitive**
- **Reproducible**
- **Homogeneous**
- **Non-radioactive**
- **Solvent compatible**
- **Compatible with multiple targets**

# ASP+ is a Fluorescent substrate



- **ASP+, fluorescent substrate for DAT and NET (1-3 $\mu$ M  $K_M$  values)** Schwartz et al. 2003, Mason et al 2005
  - $K_M$  for SERT 10-20 $\mu$ M Fowler et al. 2006
  - Trypan blue added to quench external fluorescence
- **Valuable for transporters studies**
  - Transfected cell lines
  - Synaptosomes
  - Primary cultures
- **Compatible with different platforms**
  - Fluorescent screening assays
  - High content assays
  - Confocal microscopy

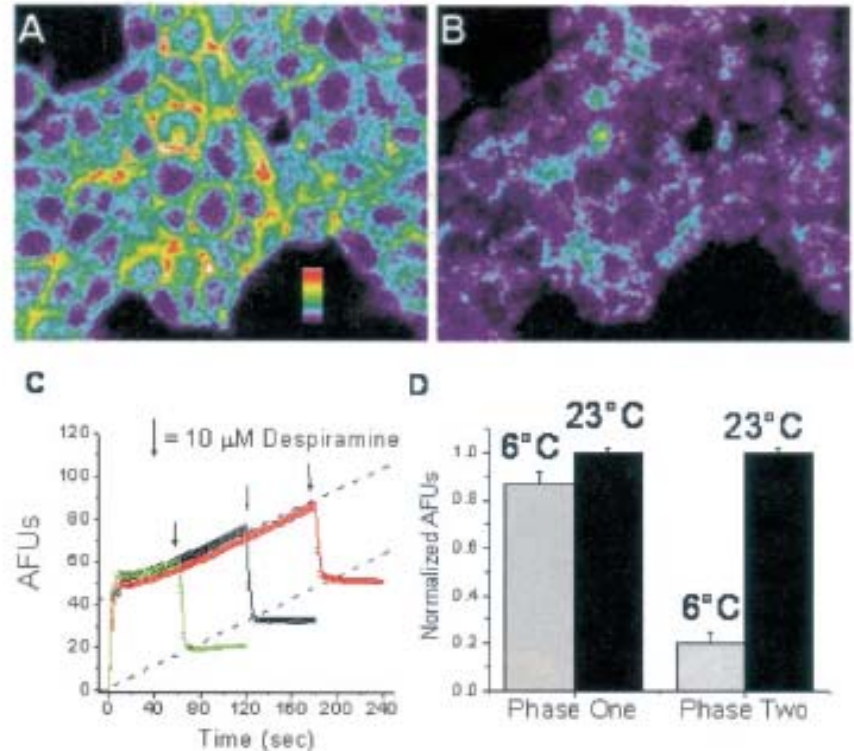


Taken from Schwartz et al. 2003 JBC 278, 9768

# ASP+ Can Be Used For Mechanistic Studies

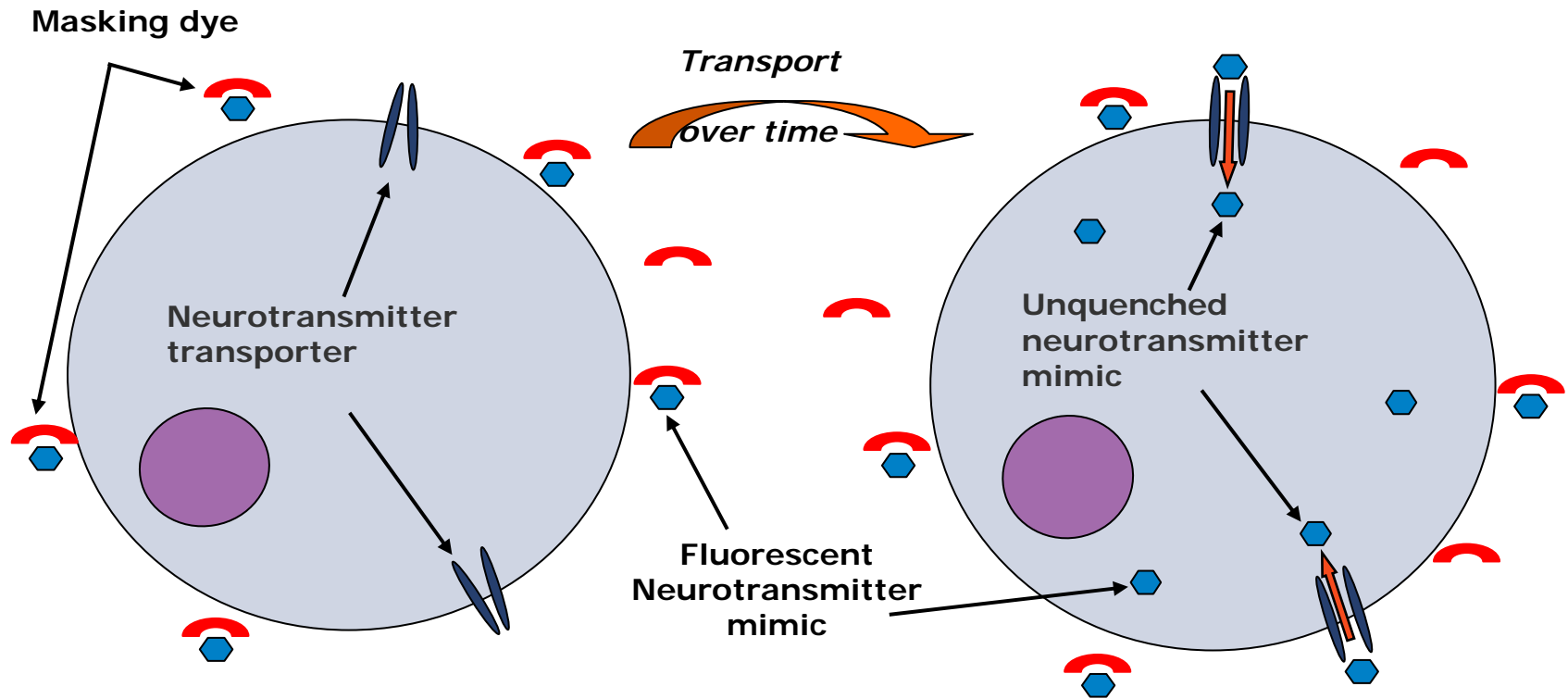


- **ASP+ rapid on rate <1sec**
  - Suitable for binding
- **ASP+ Binding is Na<sup>+</sup> independent**
  - Cell surface labeling
  - Temperature independent
  - Binding rate 100> transport rates



Taken from Schwartz et al. 2003 JBC 278, 9768

# Neurotransmitter Transporter Uptake Assay Principle



# Fluorescent Assay is Simple



Plate cells 2 days before assay

↓ Replace growth media with assay buffer

+/- Test compound

↓ Pre-incubate

Add radioactive substrate

↓ Allow uptake reaction

↓ Quench w/inhibitor

Remove radiolabel & wash

↓ Dry plates

Add scintillation fluid

↓

Read

Plate cells 1 day before assay

↓ Replace growth media with assay buffer

+/- Test compound

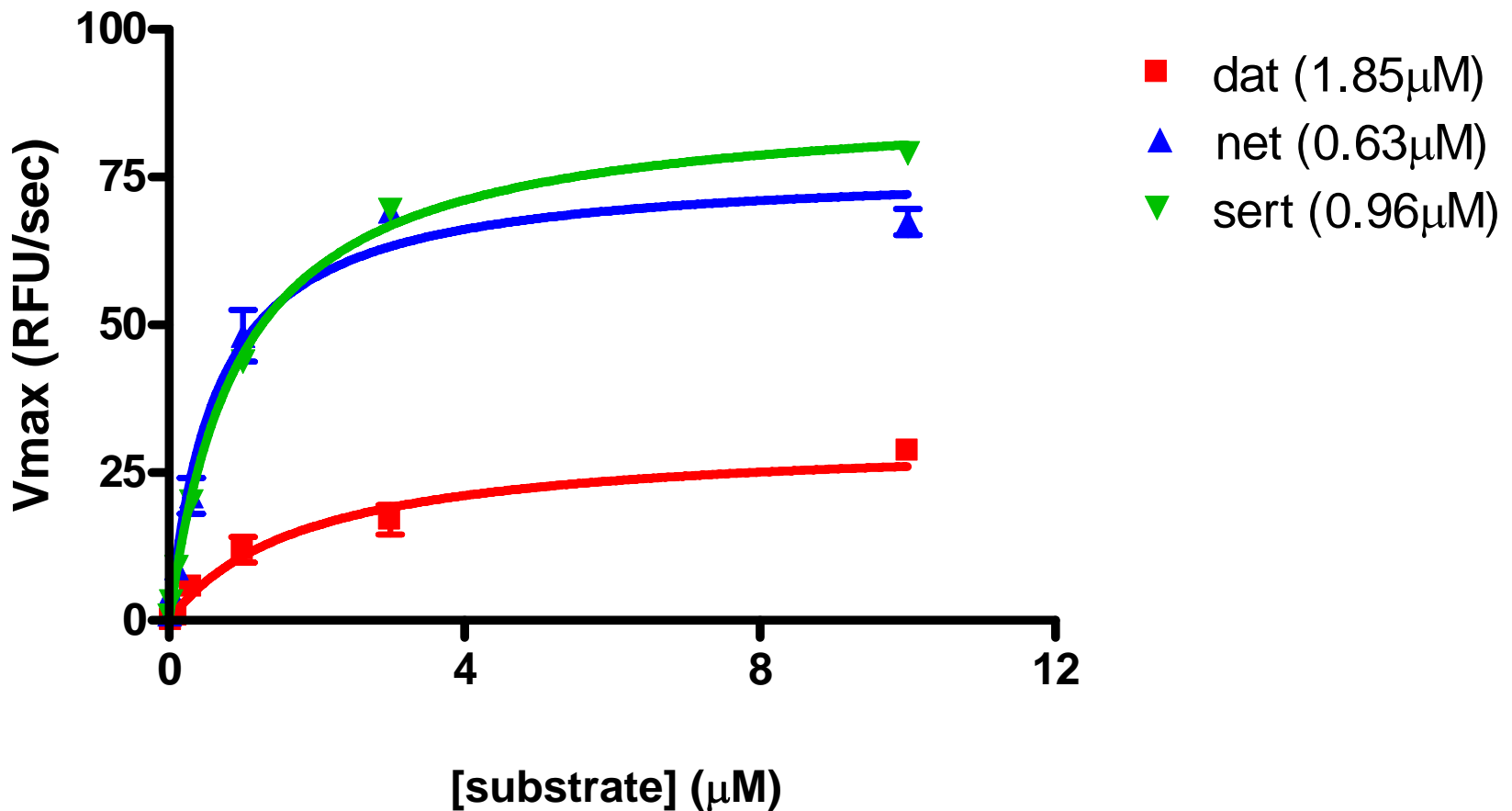
↓ Pre-incubate

+ Fluorescent substrate

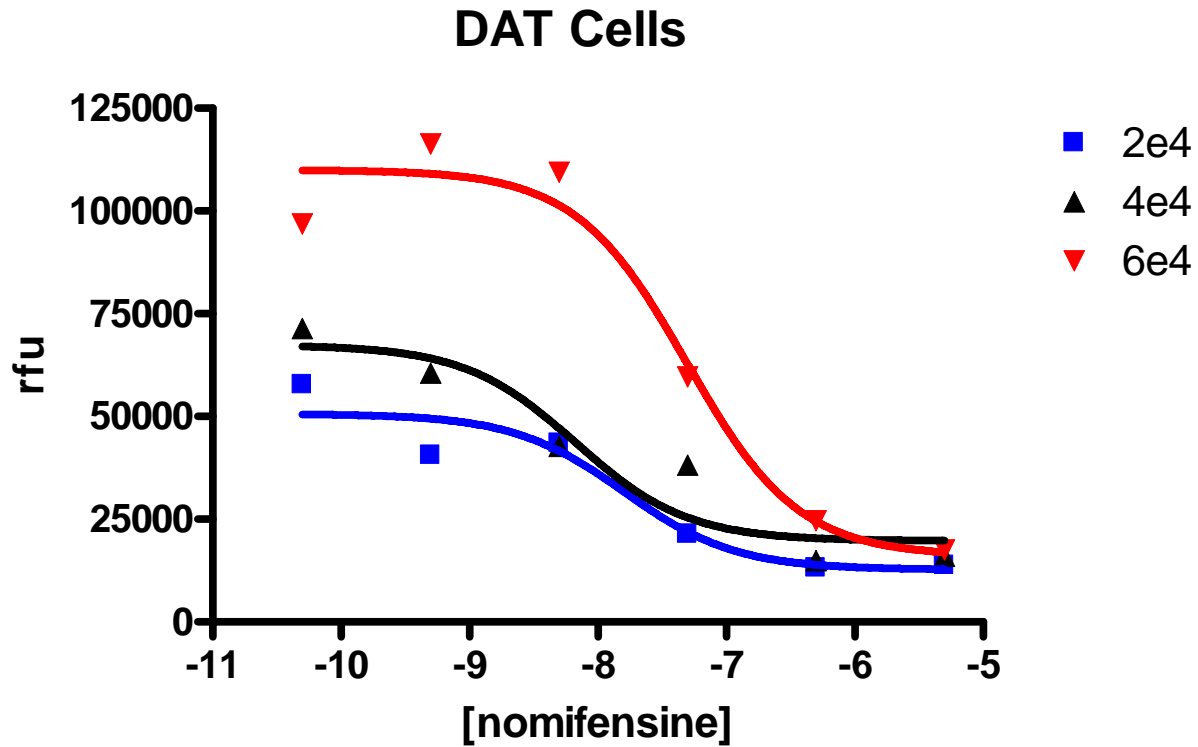
↓

Read

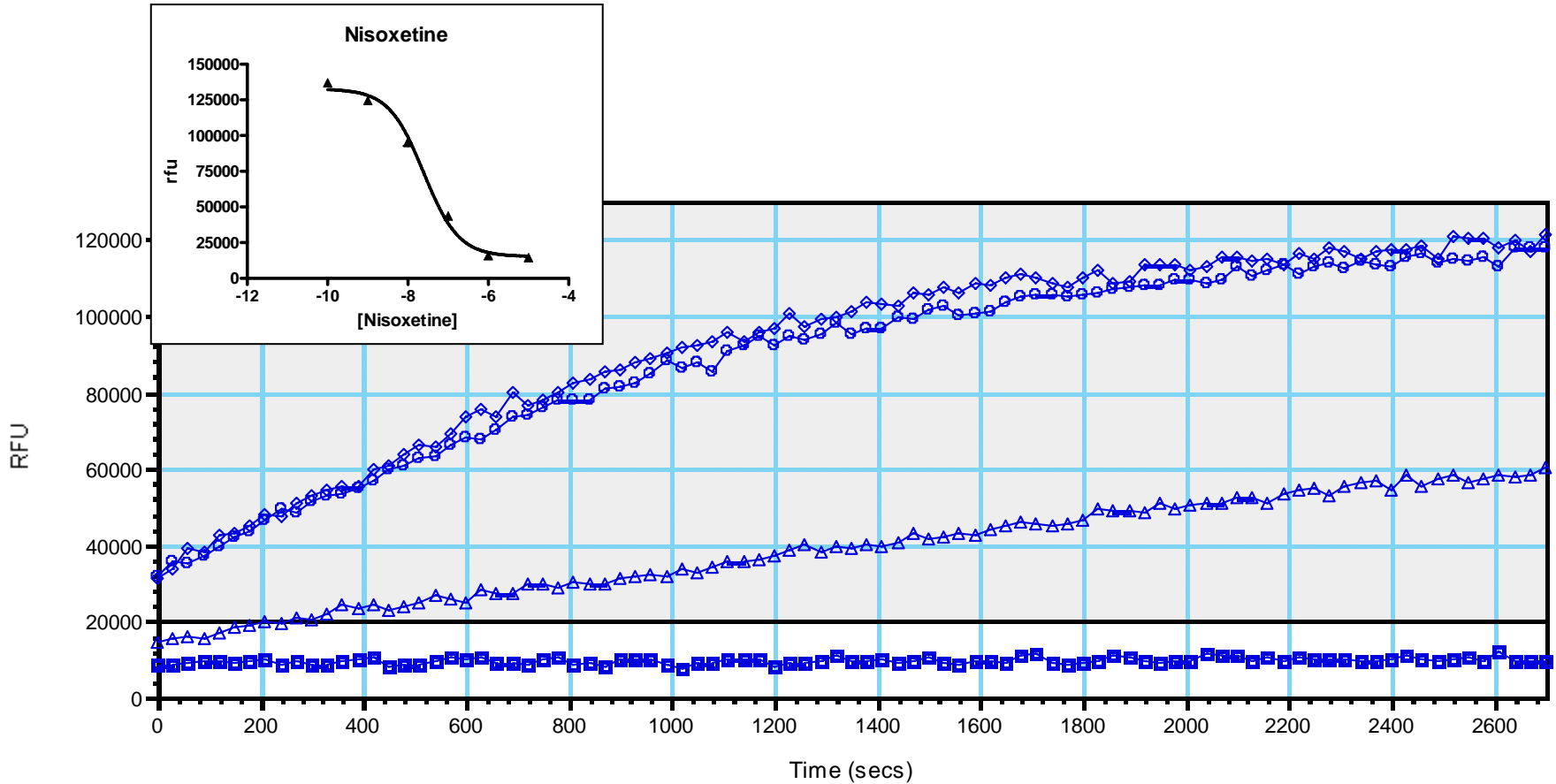
# $K_M$ is Similar for DAT, NET & SERT



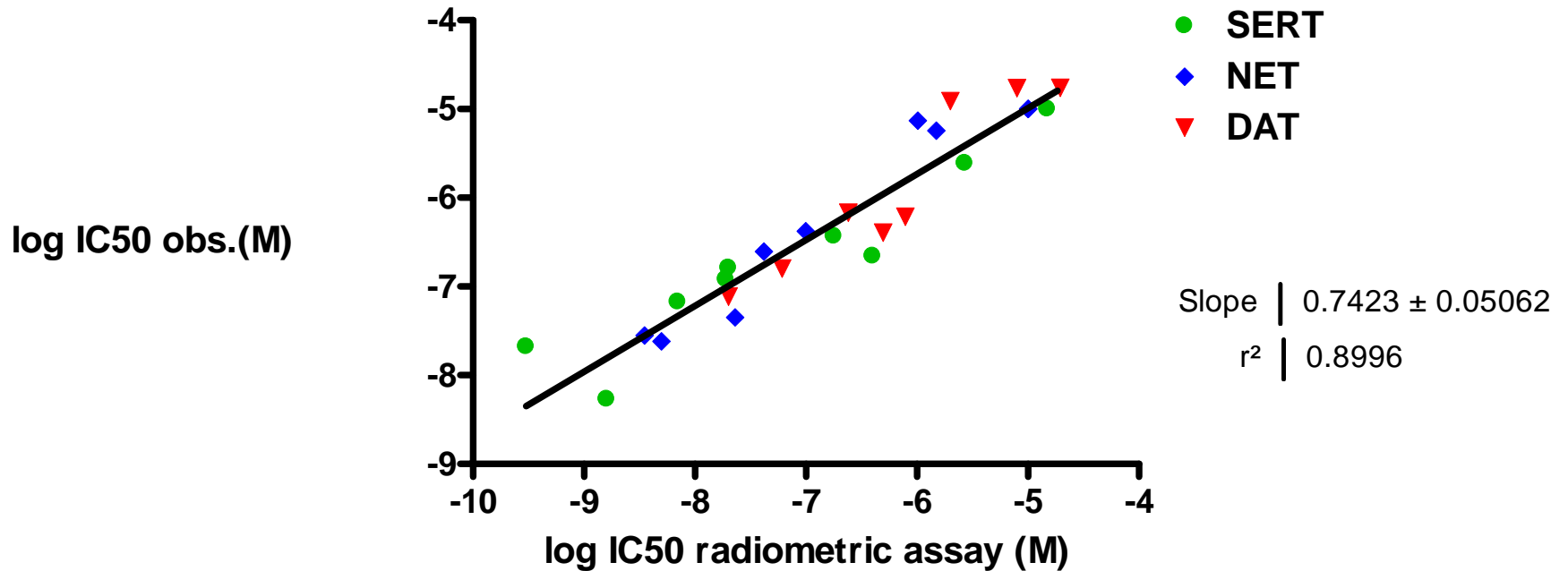
# 60,000 Cells/Well is Optimal (96 well)



# Representative Data



# Assay Yields Expected IC50s





# Comparison of Assays



## Radioactive Assay

- Uses natural substrate
- High-throughput
- Sensitive
- Reproducible
- Homogeneous
- Non-radioactive
- Solvent compatible
- Real-time analysis
- Compatible w/ multiple targets

## Fluorescent Assay

- Uses natural substrate
- High-throughput
- Sensitive
- Reproducible
- Homogeneous
- Non-radioactive
- Solvent compatible
- Real-time analysis
- Compatible w/ multiple targets

# Summary



- **Simple assay for all 3 transporters**
- **Homogeneous**
- **Sensitive (>8 fold window)**
- **Reproducible (Z' factors 0.64-0.72)**
- **IC<sub>50</sub> values comparable to radioactive assay**
- **Amenable to high throughput**
- **Commercially Available!**

# Acknowledgements



## **Sepracor**

**Nancy Spicer**

**Larry Hardy**

## **Molecular Devices**

**Rich Sportsman**

**Peter Fung**

**Annegret Boge**

**Susan Williams-Clark**