



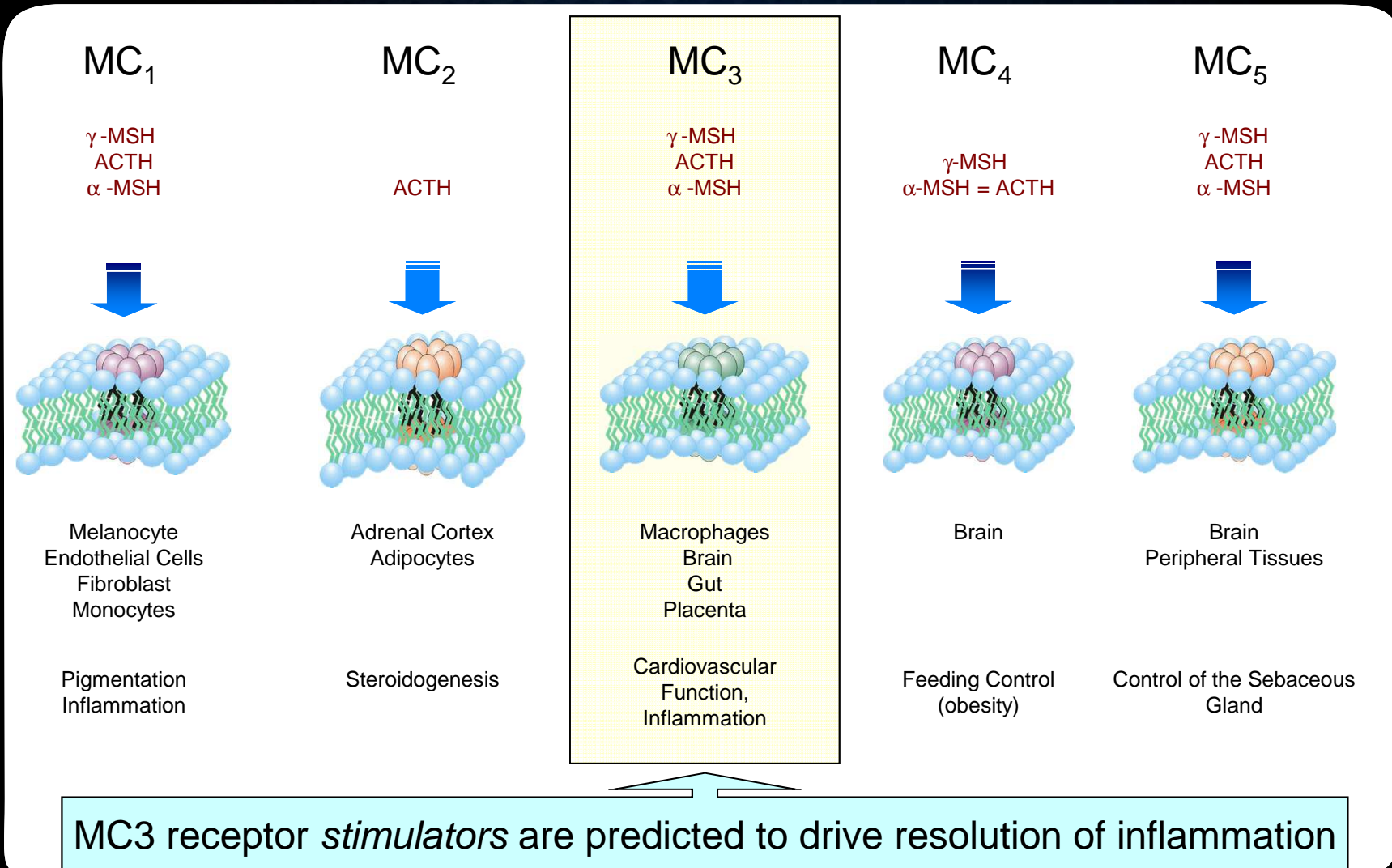
MRC Technology
Centre for Therapeutics Discovery

Identification and characterization of allosteric modulators of GPCRs: The utility of HTRF and incorporation into generalised screening strategies

Jeff Jerman
ELRIG – Cisbio Workshop
Sep 2012

- Melanocortin receptors
- Allosteric modulation of 7TM
- HTS and compound profiling considerations
 - *Major challenges*
 - *Suggested PAM screening strategies*
- Comparative Pharmacology
- Summary

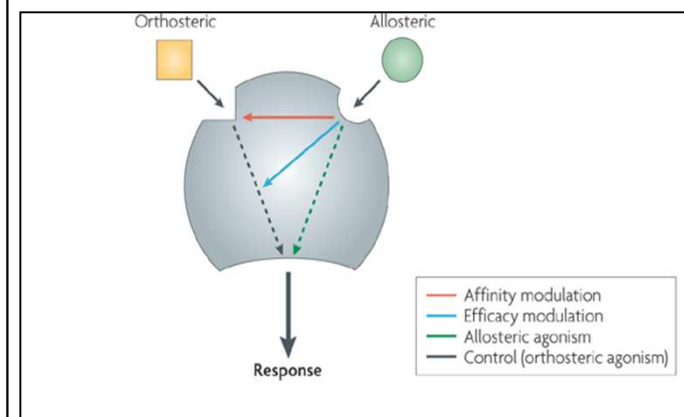
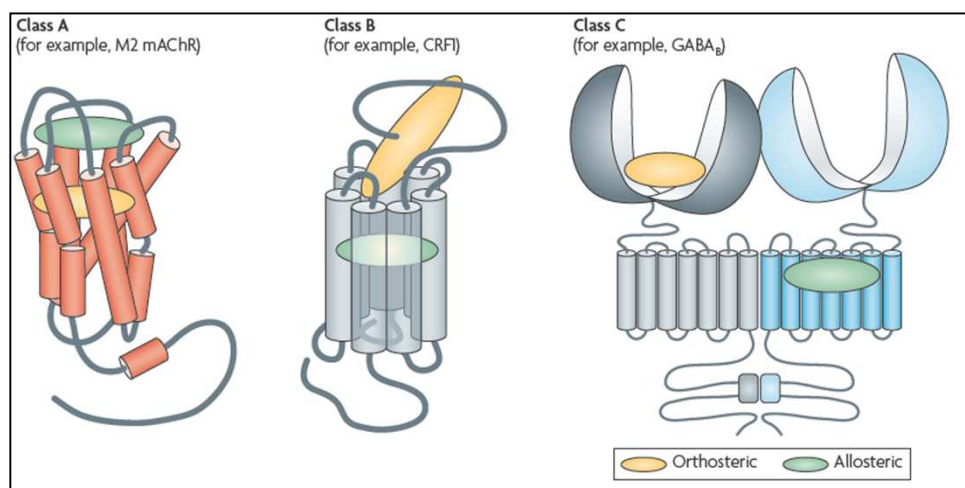
Melanocortin receptors



Targeting novel compound mechanism of action



Orthosteric vs Allosteric binding/functional modality



Advantages of Positive Allosteric Modulation (PAM)

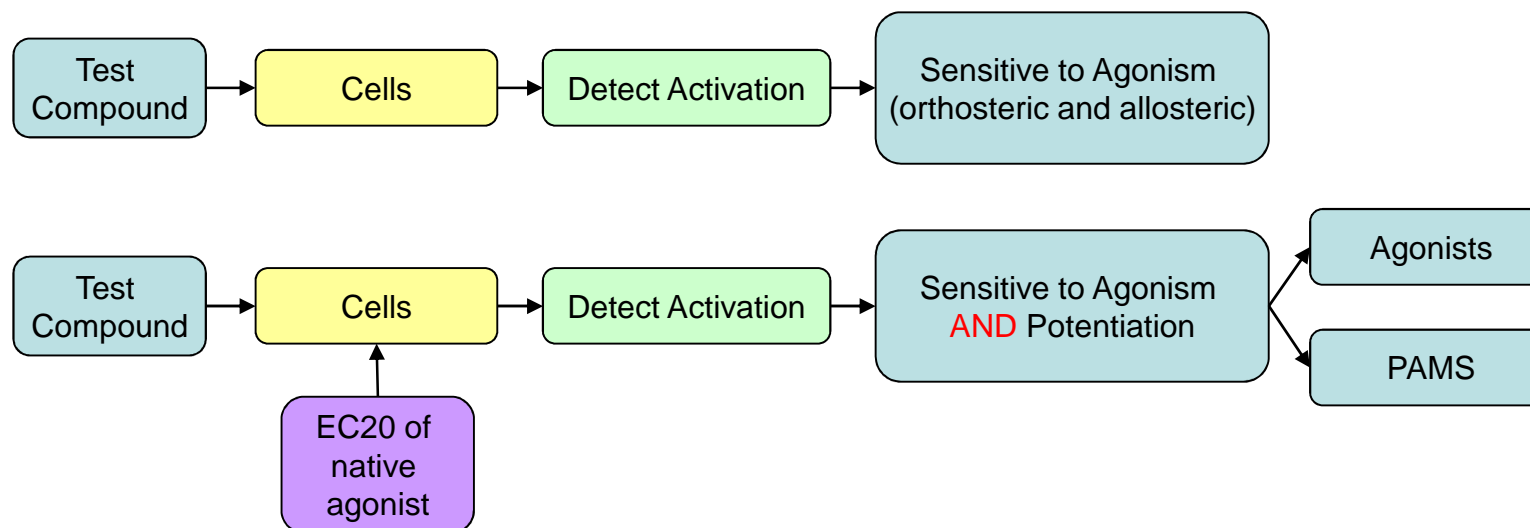
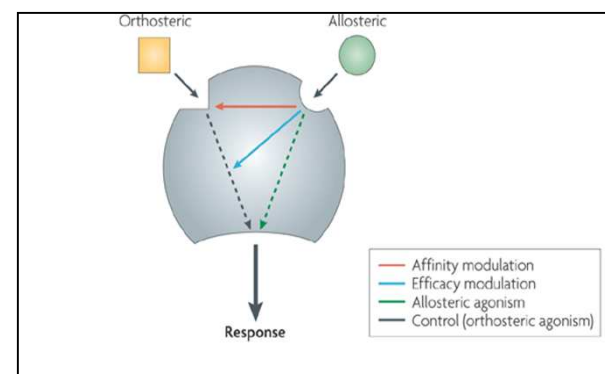
1. *Improved selectivity*
2. *Saturability (self-limiting) biological effect*
3. *Temporal and spatial resolution*

Agonist vs PAM Assay configurations



- 7TM HTS can be configured to detect both agonism and Positive Allosteric modulation simultaneously

- The 'simple' inclusion of a submaximal (EC20) of agonist facilitates this*



7TM PAM HTS/CP – Major Challenges (1)

Prediction and control of an EC20 stimulus

- *Endogenous melanocortin agonists are 'sticky' peptides (loss and/or carry over)*
- *Changes in receptor expression/coupling can dramatically affect pEC50*
- *The predictability and stability of the EC20 determines the sensitivity to PAMS*
- *HTRF affords both sensitivity and stability in response*

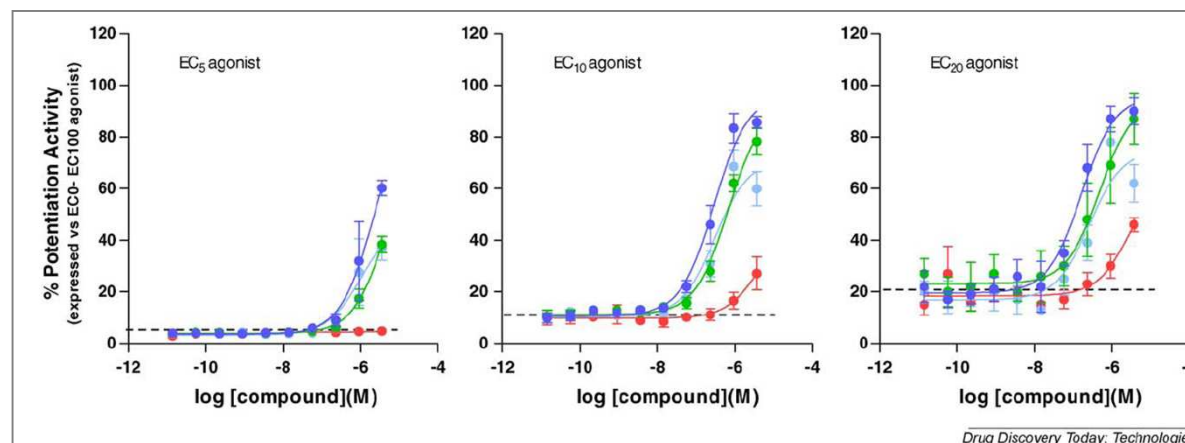
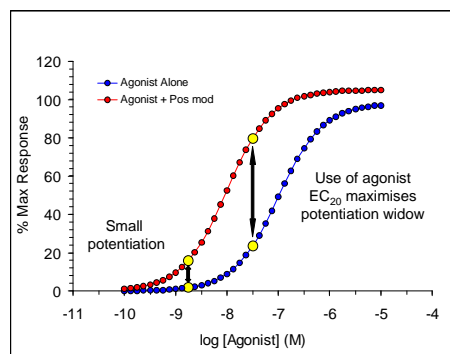


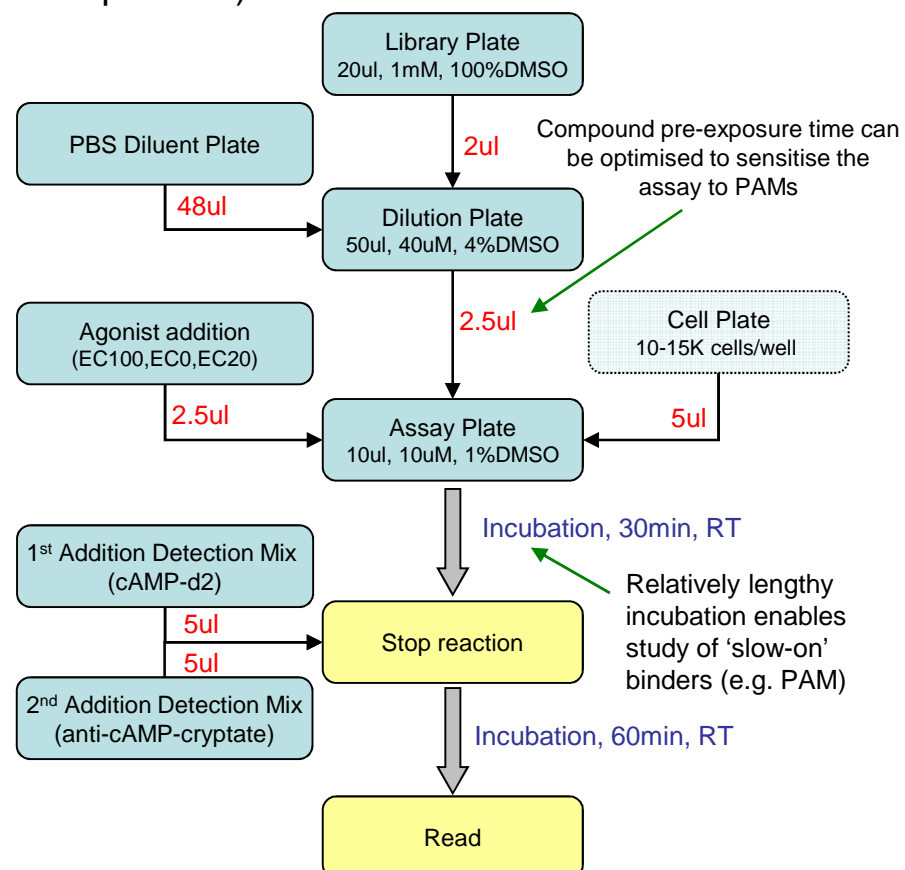
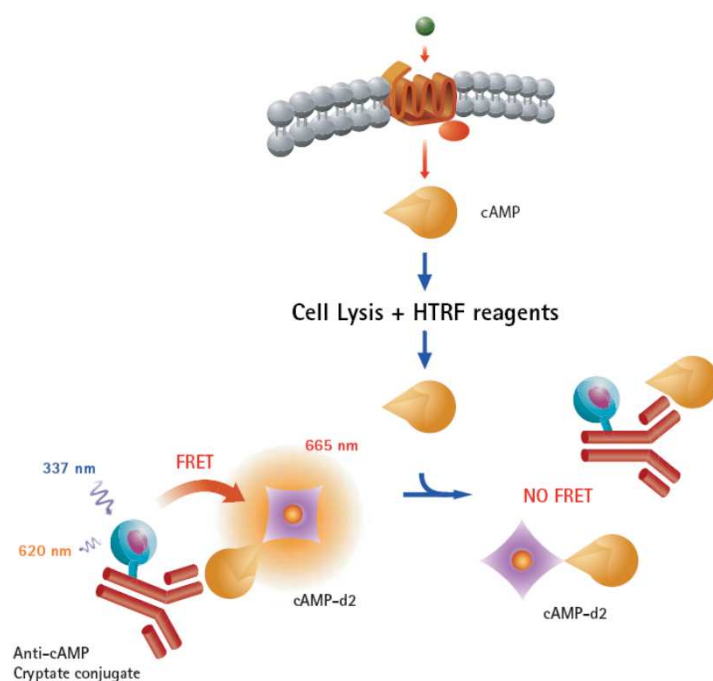
Figure 2. During screening, PAMs are typically tested as a function of a fixed agonist concentration corresponding to EC₂₀. The variability inherent to the measure of pEC₅₀ of modulation in production screening can often be under-estimated. The impact associated with using lower than anticipated agonist concentrations over plate runs and/or days on assay sensitivity is significant. The graphs below illustrate the differential modulatory profiles obtained for four compounds derived from the same chemotype series when tested at agonist EC₅, EC₁₀ and EC₂₀. Both potency and efficacy values appear to be affected to a different extent for each compound. In particular, it can be noted that one of the compounds is inactive at the lowest condition of agonist.

Drug Discovery Today: Technologies

Automated Assay Protocol



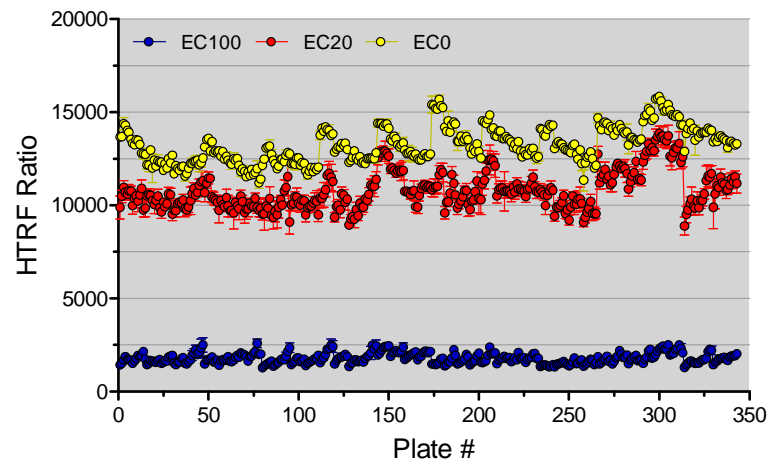
Agonist-induced cAMP accumulation (generic HTRF automation protocol)



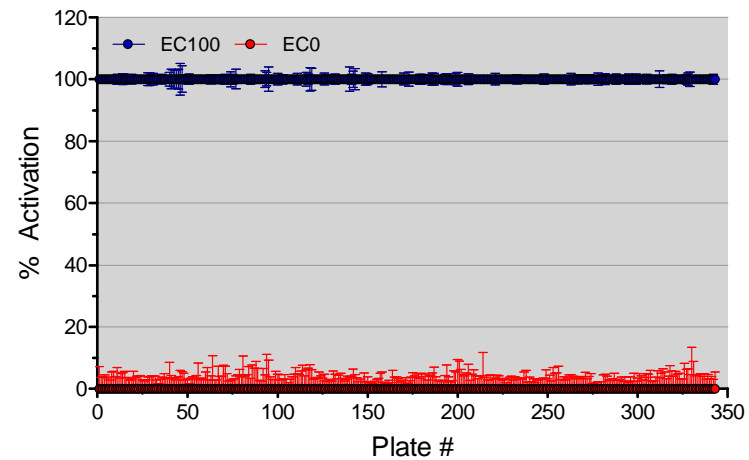
HTS Performance



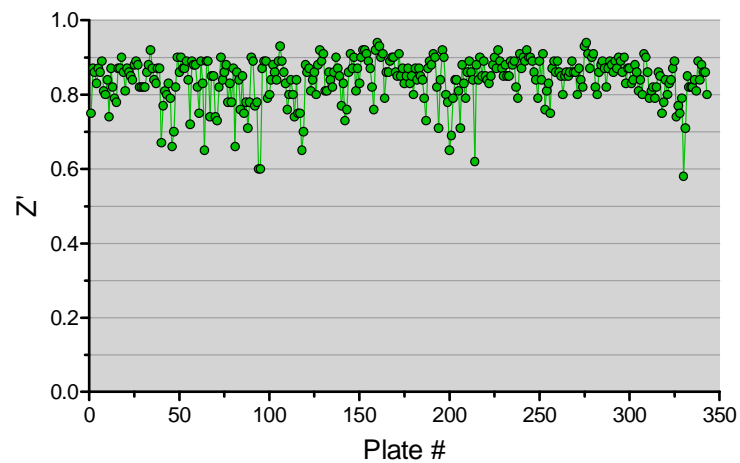
Raw Data



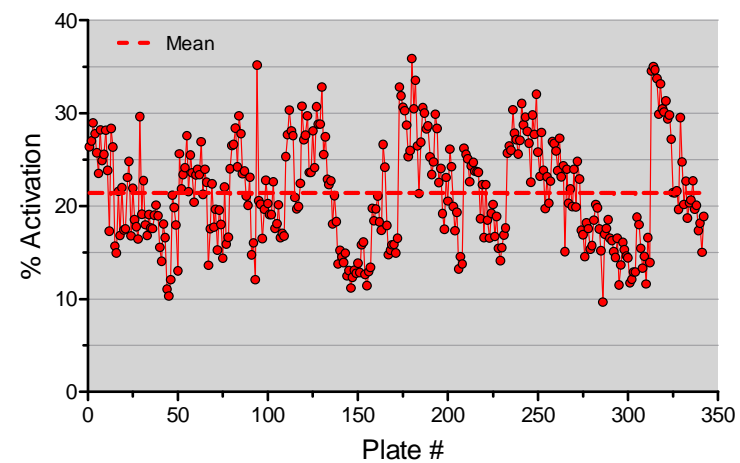
% Control (max)



Z Prime



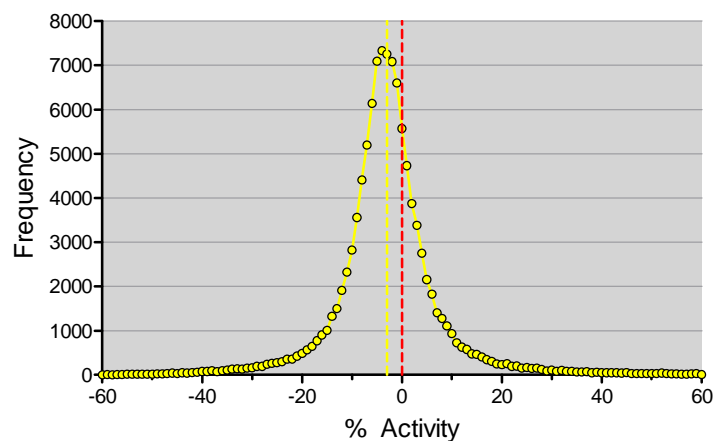
EC20 %Control (max)



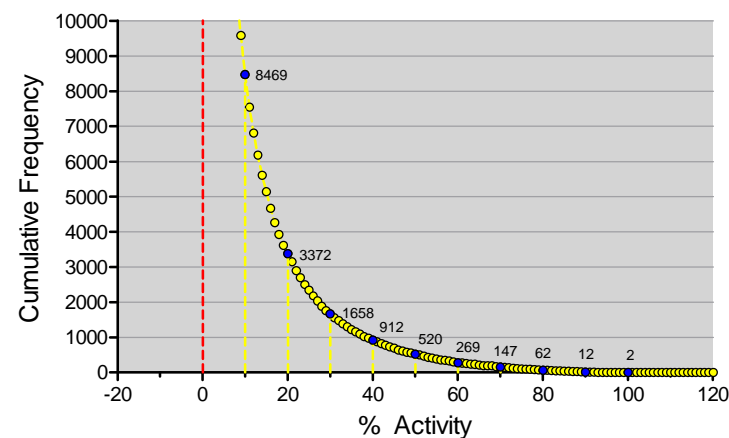
HTS Performance



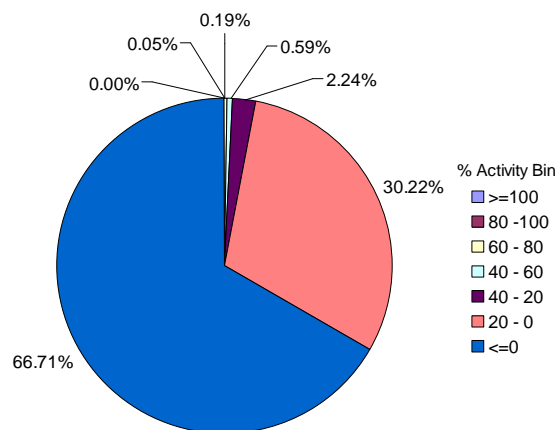
Frequency Distribution



Frequency Distribution



% Activity Distribution



109760 compounds @ 10mM (1% DMSO)

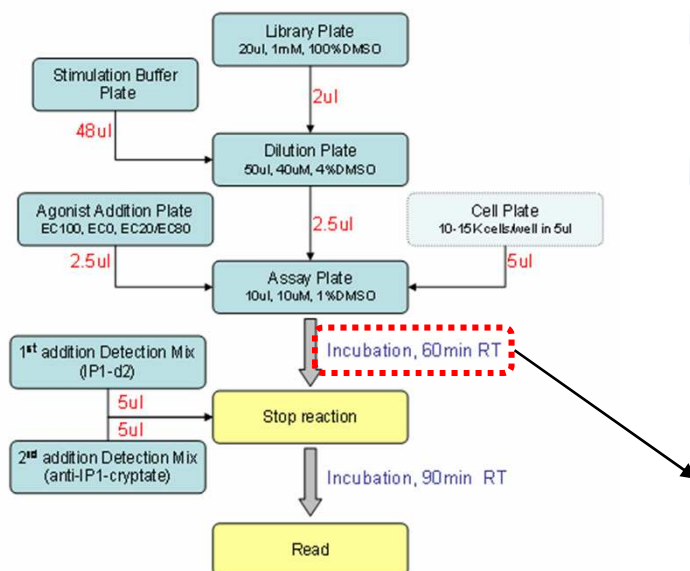
Mean Z' = 0.84 (± 0.06)

Low Control %CV = 5.3 (± 2.7)

High Control %CV = 3.6 (± 1.5)

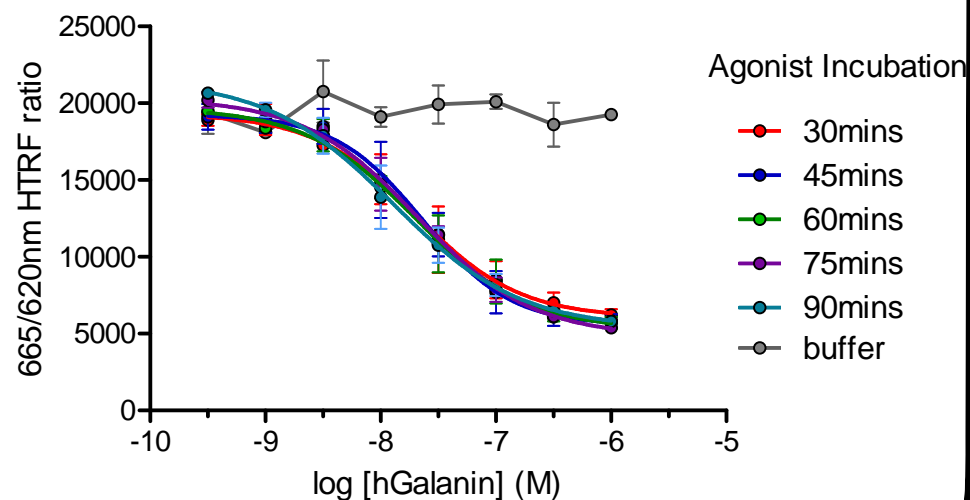
Cutoff(%)	# Hits	% HR
40	912	0.83
50	520	0.47
60	269	0.25
70	147	0.13
80	62	0.06
90	12	0.01
100	2	0.00

GalR2 IP1 HTS – Agonist incubation time

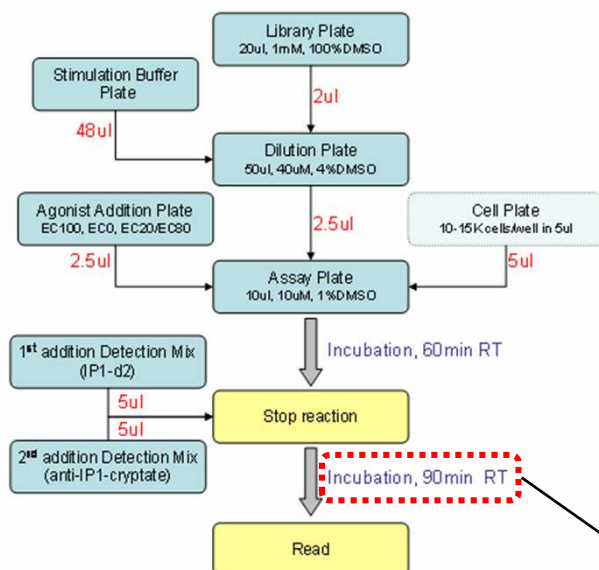


- Agonist pEC_{50} is not dependant on incubation time (within these limits)
- This suggests an EC_{20} should be stable throughout an HTS run/day

Galanin-induced IP1 accumulation in CHO (Gα16) cells stably expressing hGalR2

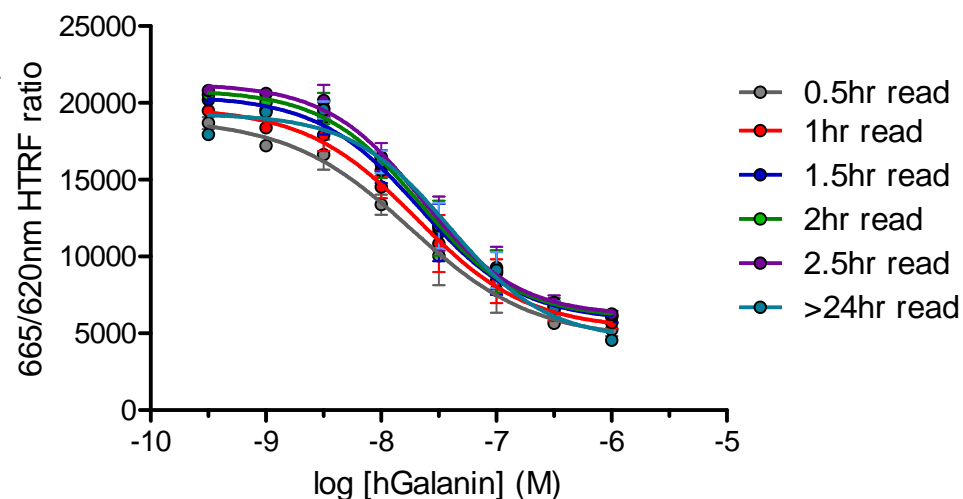


GalR2 IP1 HTS – Detection reagent stability



- Signal window is stable up to 24hrs
- Long read times (e.g. high plate numbers plates/reader delays) will not adversely affect assay performance

Galanin-induced IP1 accumulation in CHO (Gα16) cells stably expressing hGalR2

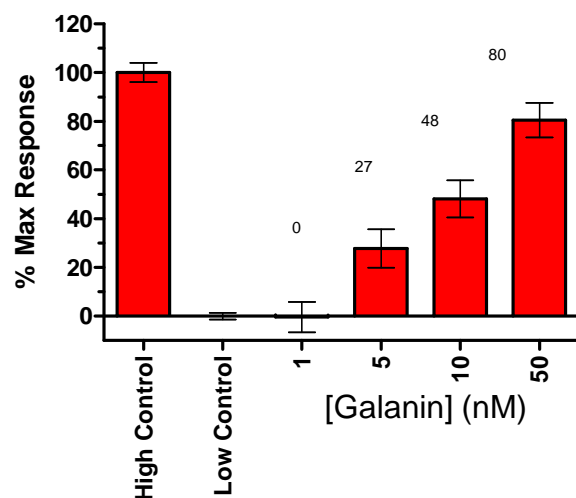


GalR2 IP1 HTS – Stability and variance of EC20



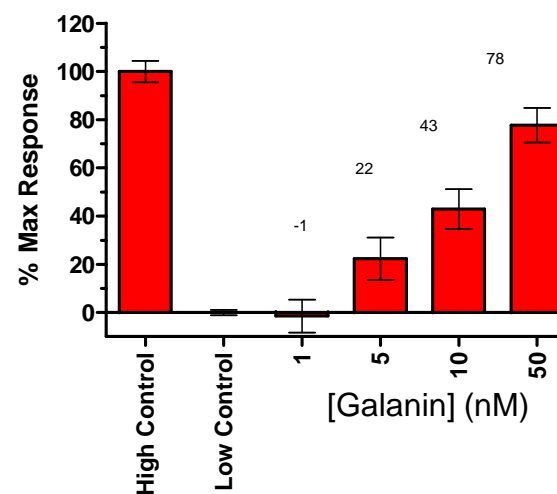
Day 1

Agonist-induced increase in [IP1] in CHO cells stably expressing GalR2 receptors



Day 2

Agonist-induced increase in [IP1] in CHO cells stably expressing GalR2 receptors



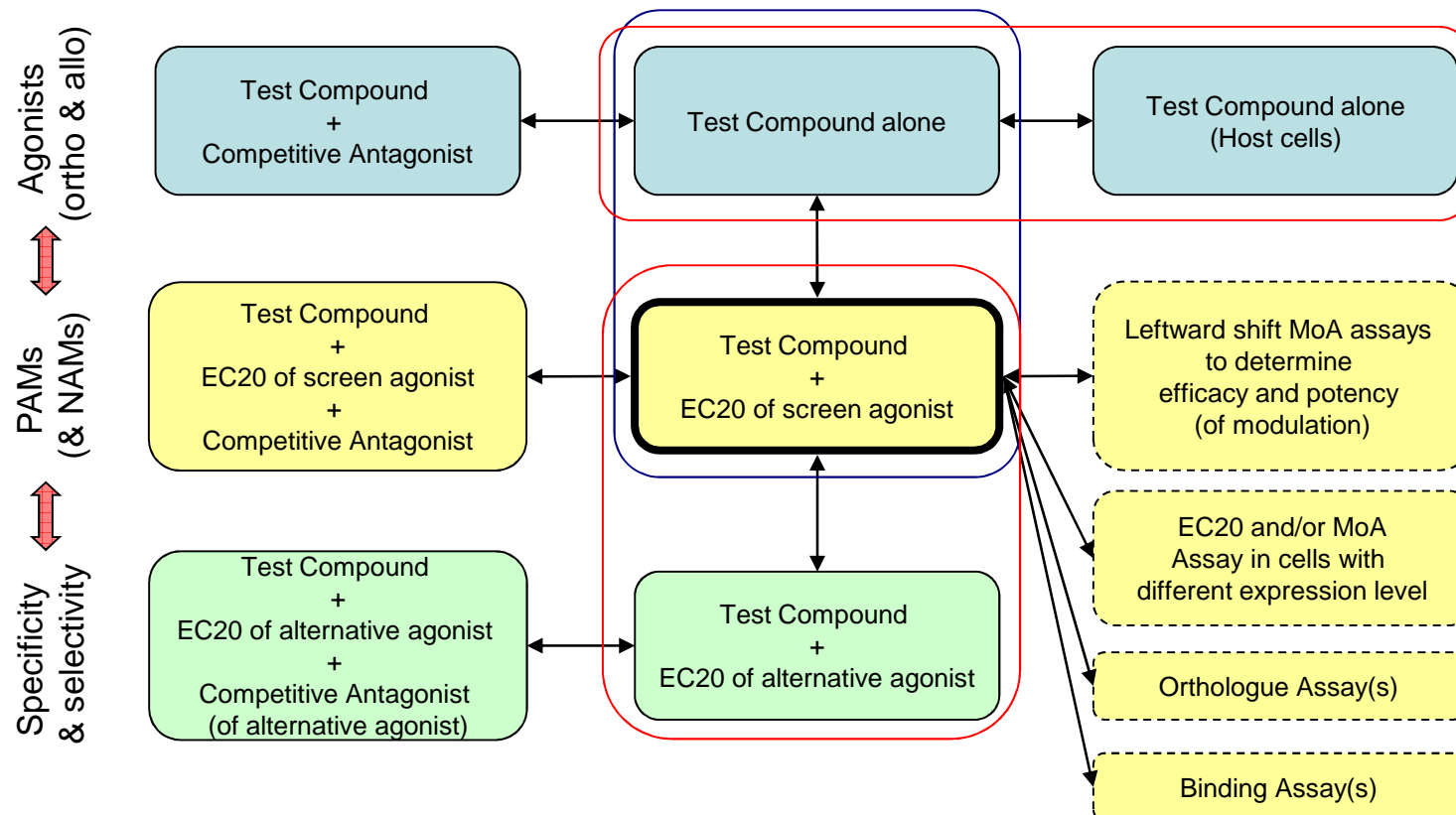
- 5nM Galanin produces a robust and stable response (~EC20)
- CVs <10%, window 2.5 fold, Z' 0.6 – 0.7

PAM HTS/CP – Major Challenges (2)

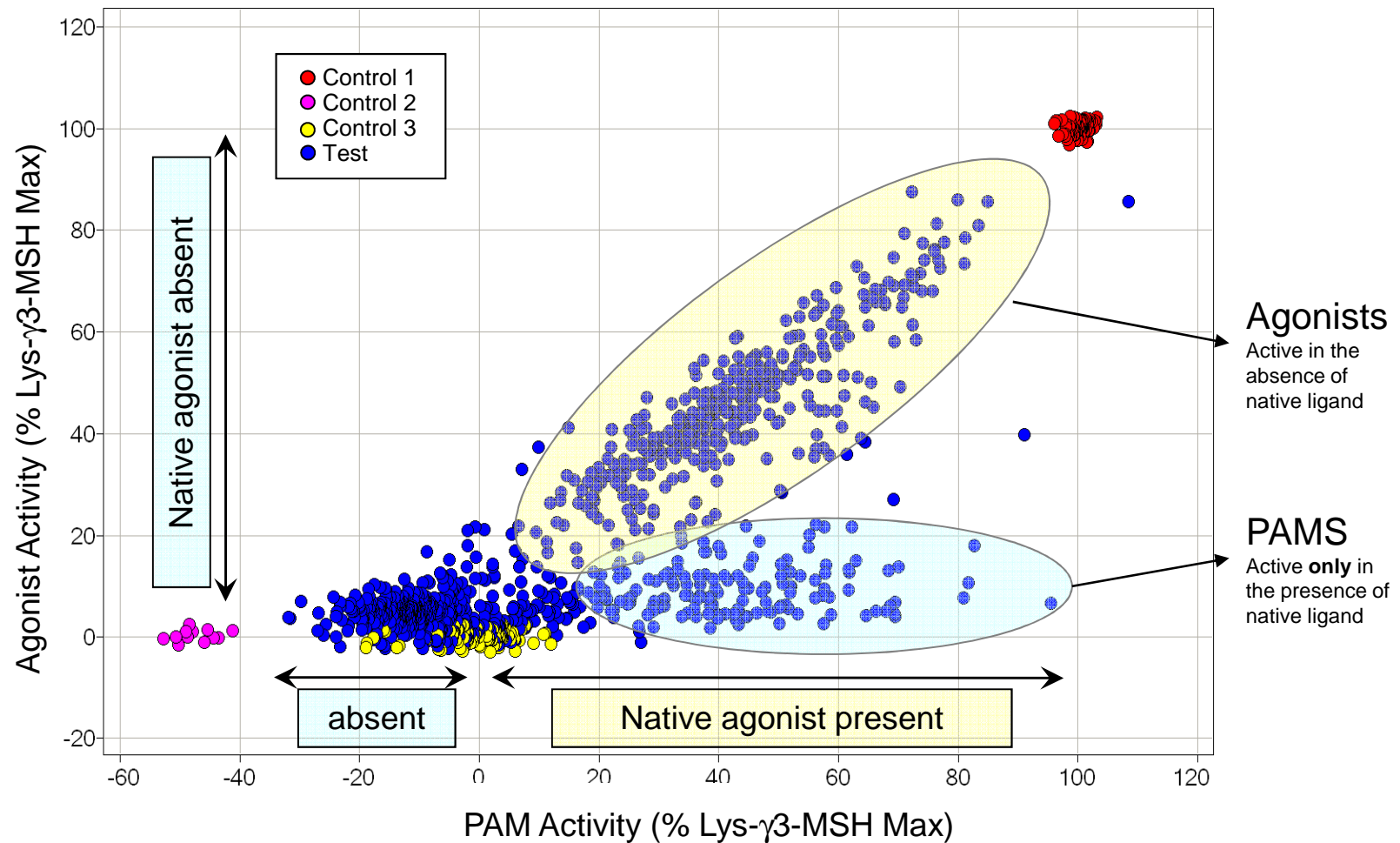


Deconvolution of PAM/Agonist hits

Removing false positives and/or non-preferred mechanism(s)



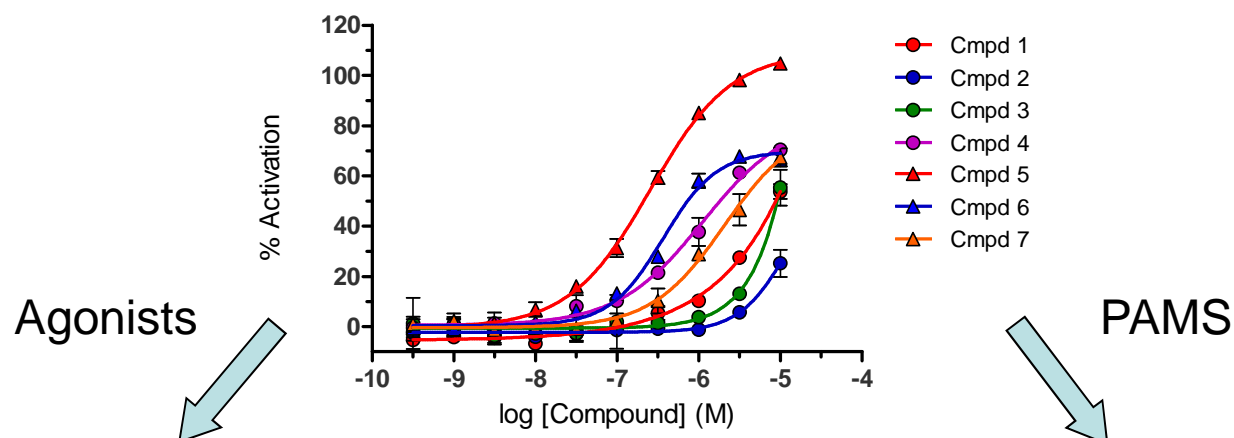
Deconvolution of PAM/Agonist modalities (Single Shot Triage)



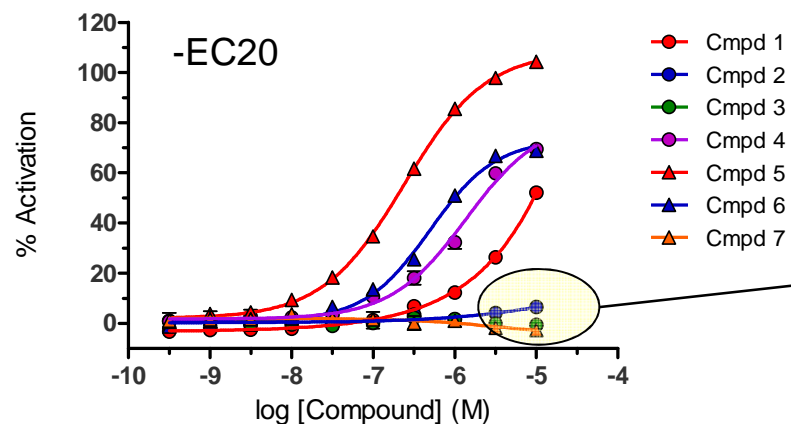
Deconvolution of PAM/Agonist modalities (Full Curve)



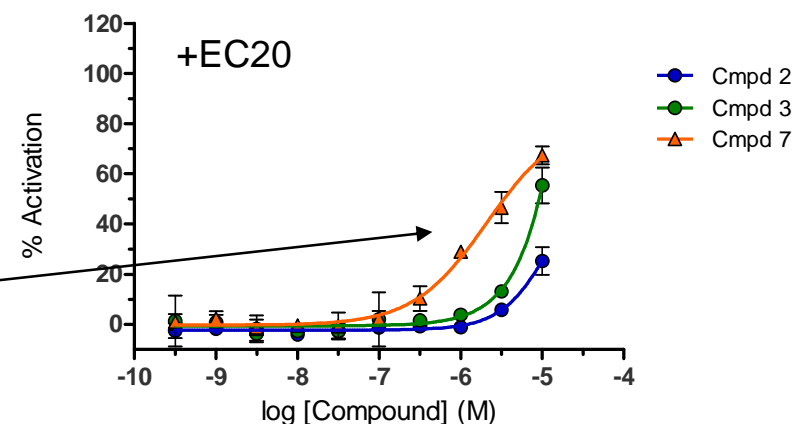
Increase in $[cAMP]_i$ in CHO cells
stably expressing MC3 receptors



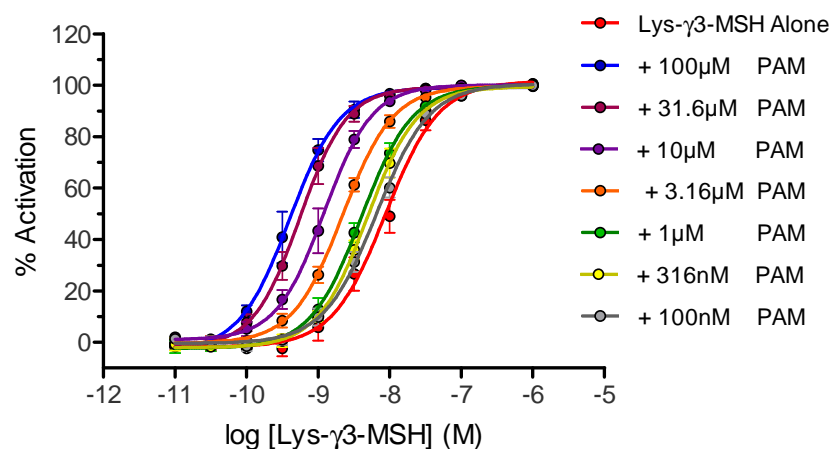
Agonist-induced increase in $[cAMP]_i$ in CHO cells
stably expressing MC3 receptors



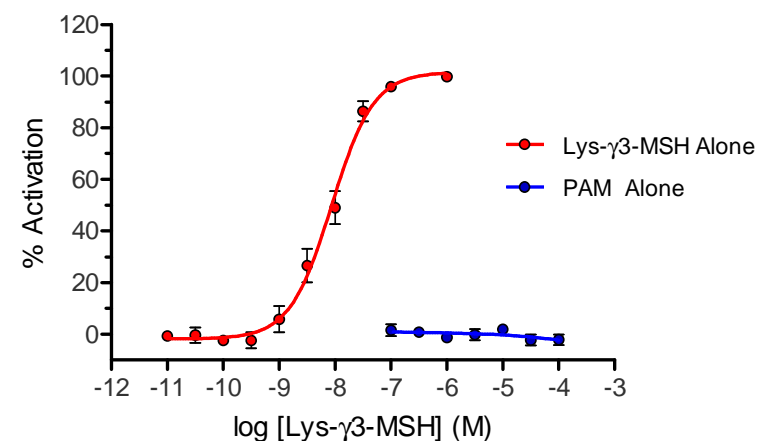
PAM-induced increase in agonist-mediated $[cAMP]_i$
in CHO cells stably expressing MC3 receptors



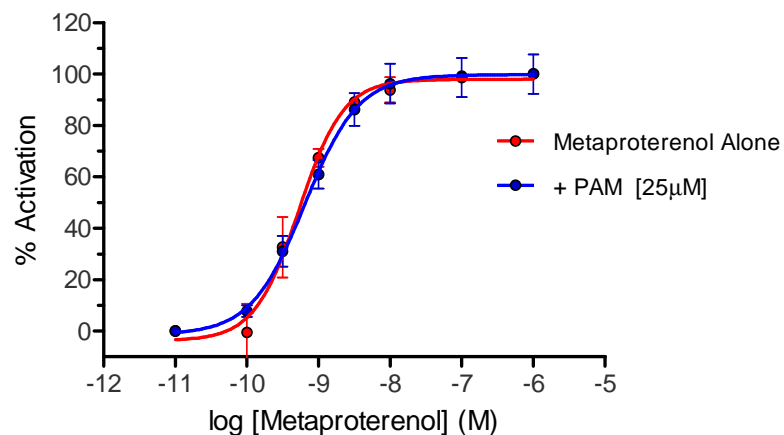
Agonist-induced increase in $[cAMP]_i$ in CHO cells
stably expressing MC3 receptors
(leftward shift)



Agonist-induced increase in $[cAMP]_i$ in CHO cells
stably expressing MC3 receptors
(leftward shift)



Agonist-induced increase in $[cAMP]_i$ in CHO cells
stably expressing β_2 -Adrenoceptors
(leftward shift)



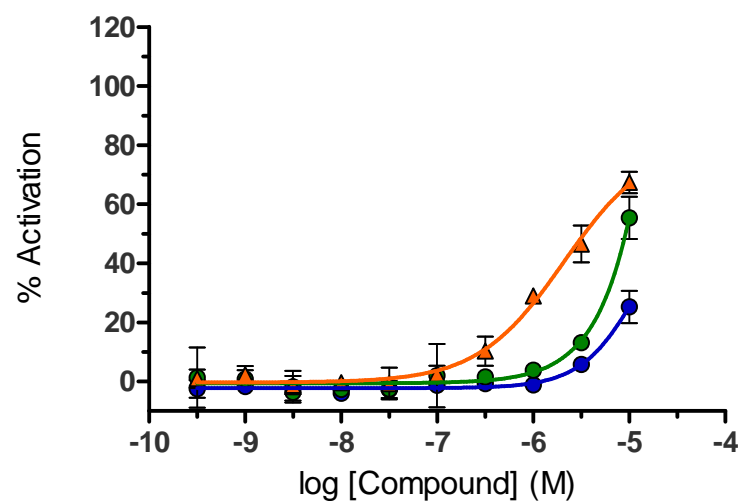
Shared robotic HTRF protocols facilitate;

- Assay in both single-point and full-curve mode (& transitions between)
- Flexibility in assay design ($pEC50_{mod}$ vs leftward shift)
- Simultaneous 'counter' assay (Ag vs PAM, off-target selectivity)

Apparent Potency in different assay formats



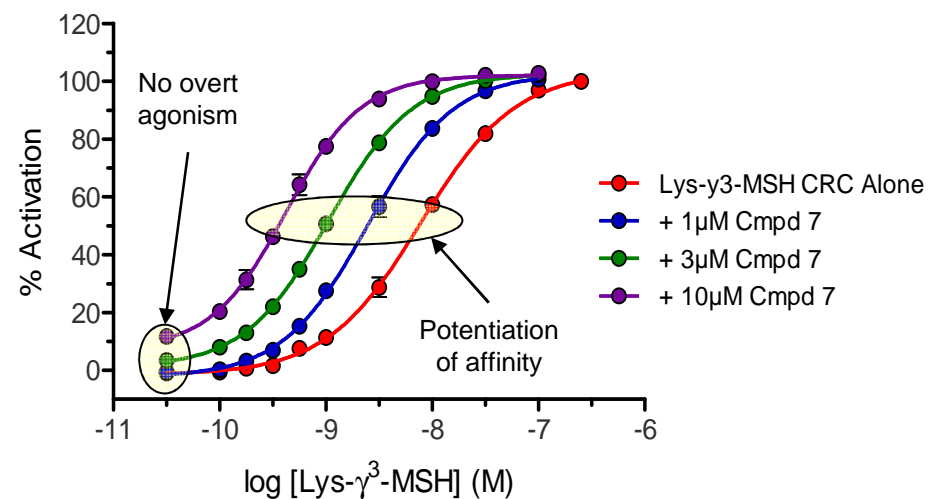
PAM-induced increase in agonist-mediated $[cAMP]_i$
in CHO cells stably expressing MC3 receptors



● Cmpd 2
● Cmpd 3
▲ Cmpd 7

Moderate pEC_{50} of modulation (EC20 mode)
translates to very effective leftward shift

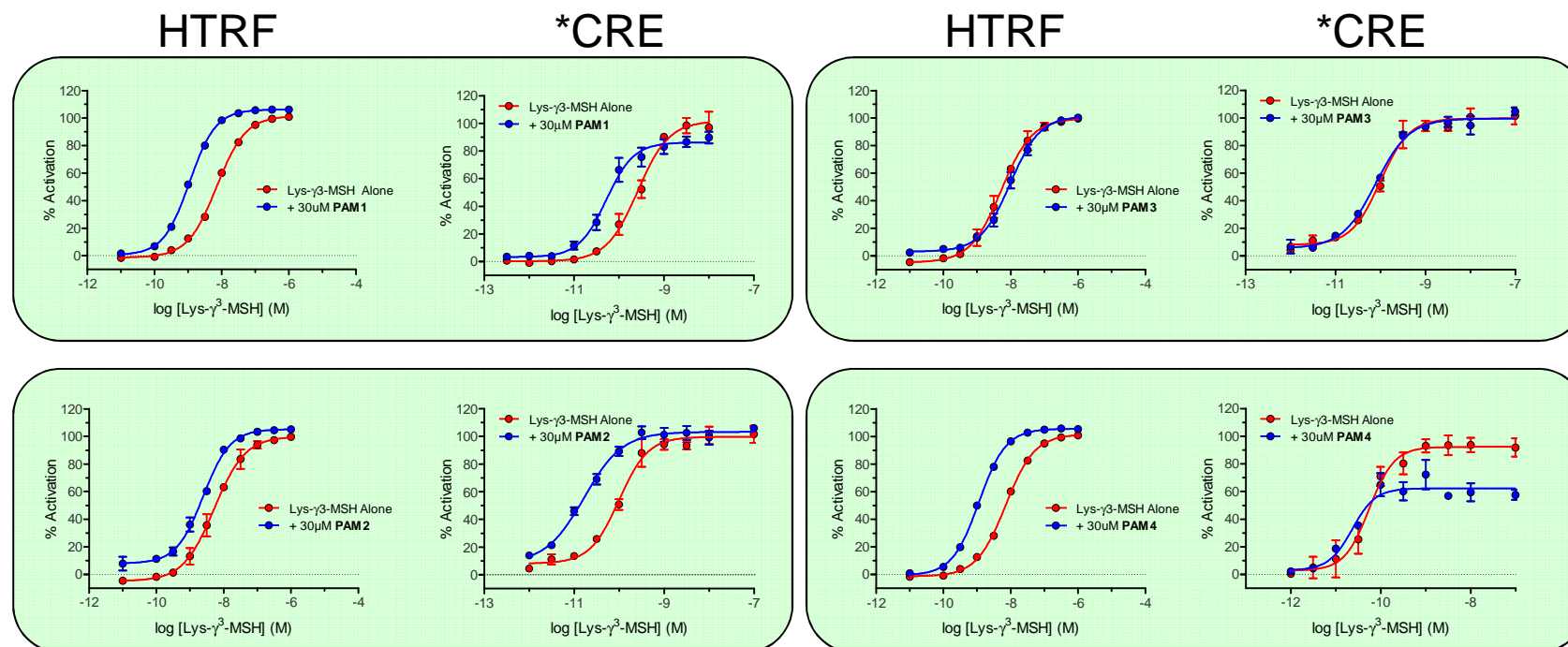
Agonist-induced increase in $[cAMP]_i$ in CHO cells
stably expressing MC3 receptors
(leftward shift)



PAM HTS/CP – Major Challenges (3)

Differences in pharmacology between cAMP detection systems

- PAM 'activity' may not necessarily align between detection formats*



*GeneBLAzer; β-lactamase coupled to a cyclic AMP response element (CRE)

- HTRF provides a sensitive and stable assay from which to configure PAM assay(s)
 - *In HTS mode the stability of EC20 is pivotal to PAM sensitivity*
- HTS and full curve hit profiling assays can be configured to share common (simple) robotised protocols
 - *pEC50 of modulation (underpinned by EC20)*
 - *Partial or full leftward shifts provide texture to PAM activity*
 - *Quantitative pharmacological analyses e.g. ETCM modelling to dissect potentiation of affinity vs efficacy*
- The technology lends itself to establishing appropriate and necessarily extensive deconvolution assays
 - *Which can share a common detection platform*

- The accumulation nature of the signal affords greater flexibility in compound pre-exposure
 - *Arguably improving the sensitivity to slow binders (PAMS)*
 - *Circumventing confounding kinetic issues with more transient detection systems (Ca^{2+})*
- The pharmacology of PAMs is complex and 'perfect' alignment with other (cAMP) detection technologies is likely to be rare
 - *A plethora of biological factors can give rise to subtle differences in apparent PAM pharmacology*

Acknowledgments



- Ahmad Kamal
- Jenny Cook
- Rachel Forfar
- Debra Taylor
- Hayley Jones
- Paul Wright
- Puneet Khurana



University of Bristol

- David Wynick



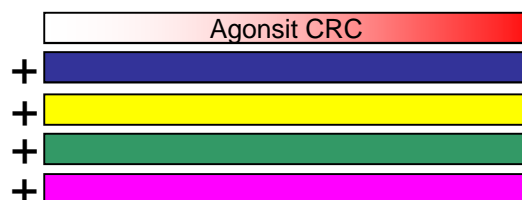
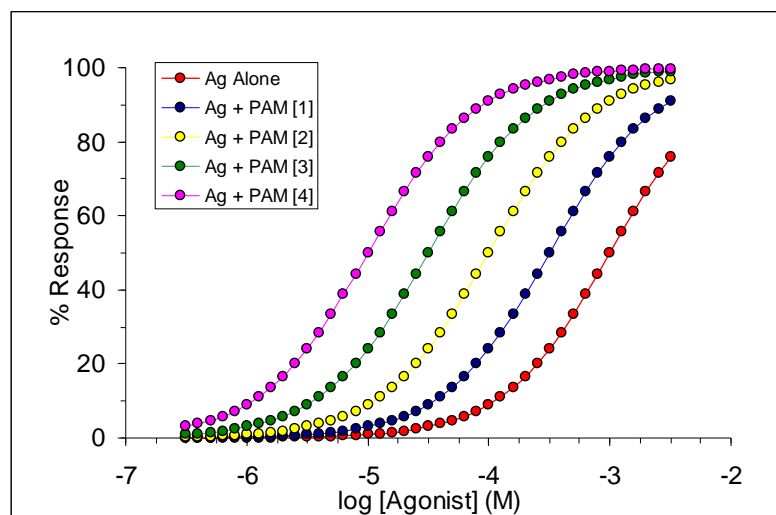
William Harvey Research Institute, Queen Mary University

- Mauro Perretti
- Trinidad Montero-Melendez

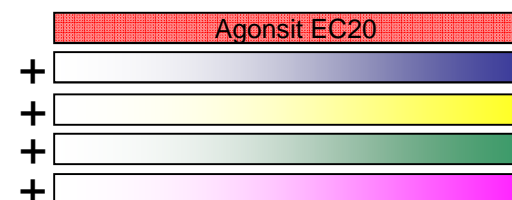
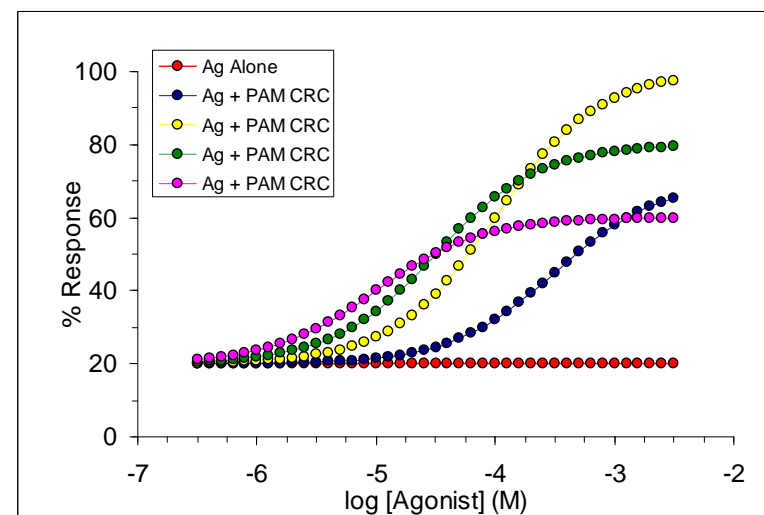
Alternative PAM assay configurations



'Leftward Shift'



'EC20'



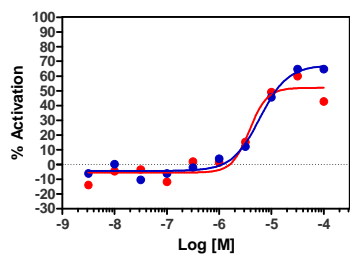
Exemplar curve signatures



PAM MC3

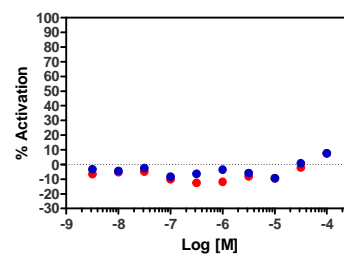
'Clean' PAM

MCR3 PAM cAMP HTRF - Plate 4, Cpd 15



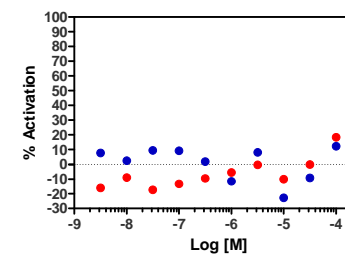
Ag MC3

MCR3 AG cAMP HTRF - Plate 4, Cpd 15



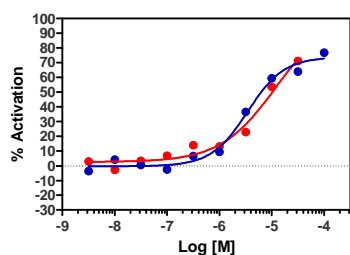
PAM β 2

ADRB2 cAMP HTRF - Plate 4, Cpd 15

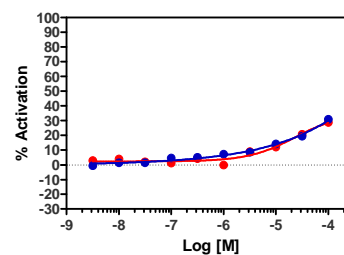


PAM with (allo) agonism

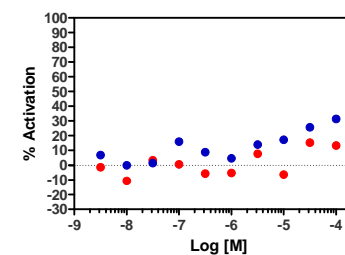
MCR3 PAM cAMP HTRF - Plate 1, Cpd 13



MCR3 AG cAMP HTRF - Plate 1, Cpd 13

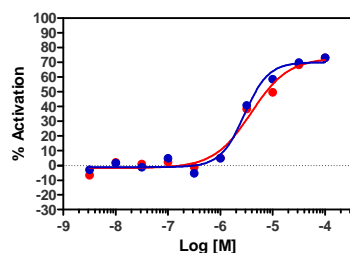


ADRB2 cAMP HTRF - Plate 1, Cpd 13

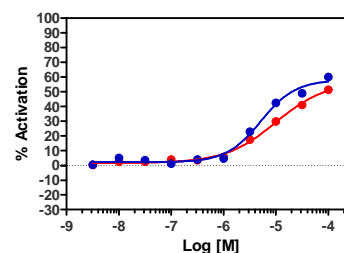


Non specific compound
(receptor and or mechanism)

MCR3 PAM cAMP HTRF - Plate 4, Cpd 11



MCR3 AG cAMP HTRF - Plate 4, Cpd 11



ADRB2 cAMP HTRF - Plate 4, Cpd 11

