

Allosteric modulator assay with IP-One.

Definition

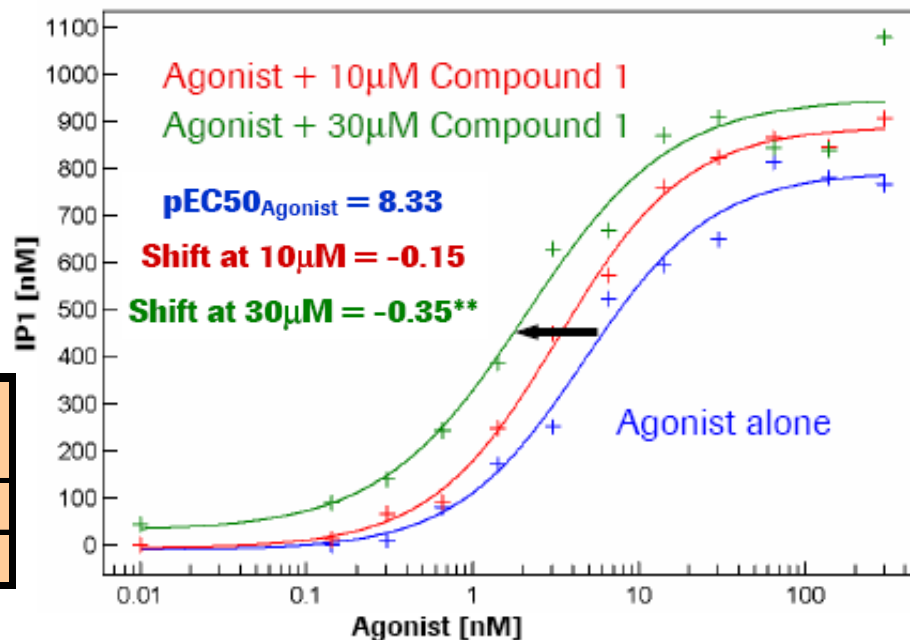
An **allosteric modulator** binds to its own binding site on the receptor and produces an effect on the receptor response to agonist stimulation by inducing a change in receptor conformation. Allosteric modulators can effect the efficacy of the receptor's affinity for the agonist.



Assay in action

Comparison:
Agonist Shifts FLIPR vs. IP-One

@ 10 μ M	Shift FLIPR	Shift IP-One
Cmpd 1	-0.36	-0.15
Cmpd 2	-0.40	-0.32



Conclusion

IP-One HTRF is an alternative method to intracellular Ca²⁺ measurement for assessing allosteric modulator activity upon agonist stimulation.