

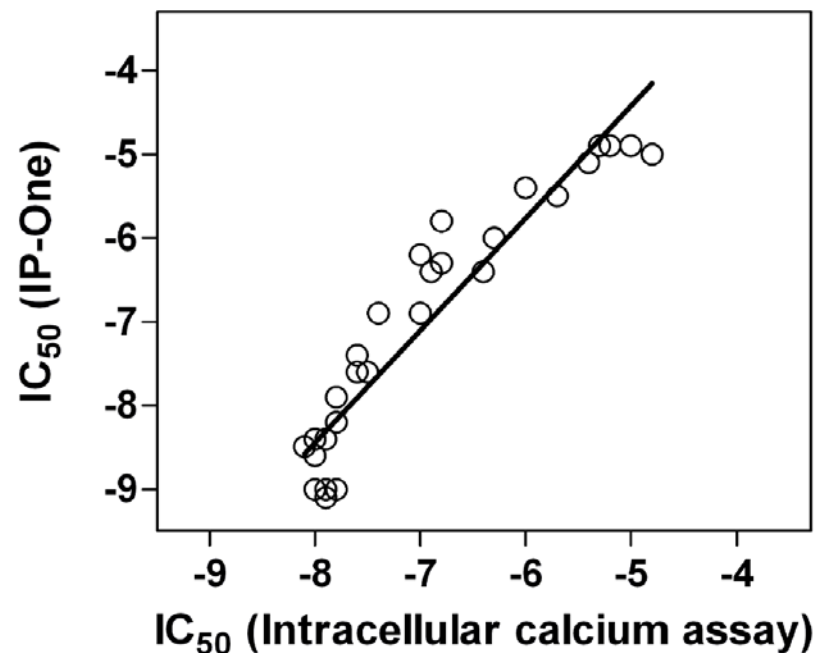
A comparison of IP-One and FLIPR (calcium) assays. Correlation of 30 known cholinergic antagonists activities between IP-One HTRF and FLIPR assays.

Assay description

Cellular model: Frozen CHO-M1 resuspended and plated in 1536 well plates 2,000 cell / well for the calcium and the IP-one assays and were incubated at 37° C overnight.

Compounds: A 1208 LOPAC library is tested in HTS condition. All the compounds are titrated in 15-concentrations.

The IC50 values of 30 known cholinergic antagonists determined from both screens are correlated (figure).



Conclusion

IC50 values of cholinergic antagonists tested are similar on FLIPR and IP-One assay.

Wei Sheng - Current Chemical Genomics, 1, 70-78 2008