

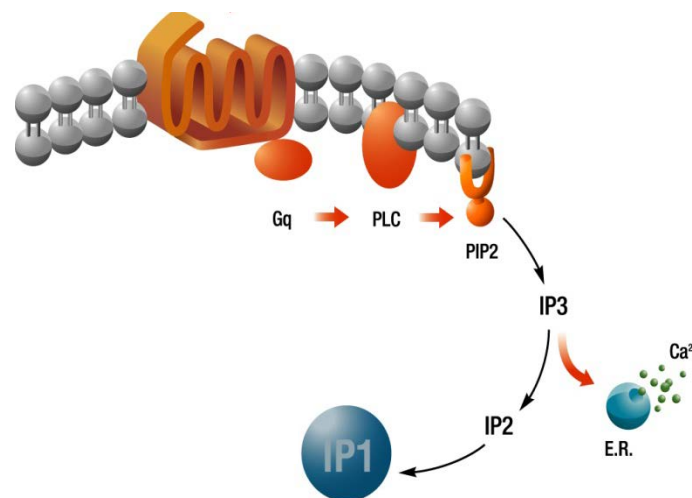
# A comparison of IP-One and FLIPR (calcium) assays. IC<sub>50</sub> & ranking of a selection anti-histaminic blockbuster.

## Assay description

**Cellular model:** Recombinant histamine H1 receptor expressed in CHOK1 cells.

Cells were seeded in 384 well plates at 10,000 cell / well and 30,000 cells/well respectively for the calcium and the IP-one assays and were incubated for 24 hours.

**Compounds:** Several marketed anti-H1 histamine blockbusters were tested in the two assays against 100 nM histamine.



	IC <sub>50</sub> (nM)	
	IP-One	FLIPR
Ketotifen	11	6.5
Mirtazapine	10	10
Doxepin	10	23
Clemastine	31	36
Cetirizine	127	1188
Astemizole	620	459
Loratadine	787	3130

## Results

Similar IC<sub>50</sub>'s and ranking are obtained for most of the blockbusters except in one case, **cetirizine**, where the data are in favor of the IP-One assay.