

## qHTS for Identifying the Cell Membrane Permeable Inositol Monophosphatase Inhibitors with HTRF IP-One Reagents

Wei Zheng

National Institutes of Health, National Human Genome Research Institute, NIH Chemical Genomics Center

Lithium has been used for the treatment of Bipolar Disorder in last 50 years with a narrow therapeutic index. Despite its severe adverse effects, it remains a main therapeutic choice for the patients with acute mania. Inositol monophosphatase (IMPase) has been speculated as a target of lithium therapeutic action. The traditional enzyme-based assays have failed to discover the cell-membrane permeable and bioavailable IMPase inhibitors. Cell-based enzyme assay offers a new approach for identifying inhibitors of intracellular enzymes. We have optimized and miniaturized the HTRF inositol-1-phosphate (IP1) assay in 1536-well plate format and applied it for a screen of ~100, 000 compounds. A total of 30 potential IMPase inhibitors were identified from the primary screening. The confirmation and follow-up assays on these active compounds are currently carrying out. Our results indicated that this HTRF IP-One assay is cost effective and robust for high throughput screening.