



# PREDEASY™ ATPase Assay Kit

**SOLVO Biotechnology introduces the revolutionary PREDEASY™ ATPase Assay Kit.**

A unique product for measuring the ATPase-dependent Drug/ABC membrane transporter interaction in a convenient, "ready to assay" package!

For the ultimate in convenience, **PREDEASY™** kits contain all you need to conduct ATPase activation and inhibition assays, including all necessary reagents and solutions. Each kit also comes with a CD that includes the assay protocol, data sheet, MSDS, electronic data processing and evaluation templates and example data. The recommended layout allows activation and inhibition data in duplicate for 8 concentrations of the test article with all of the relevant controls.

The Kit delivers major technical improvements over the conventional ATPase products, such as:

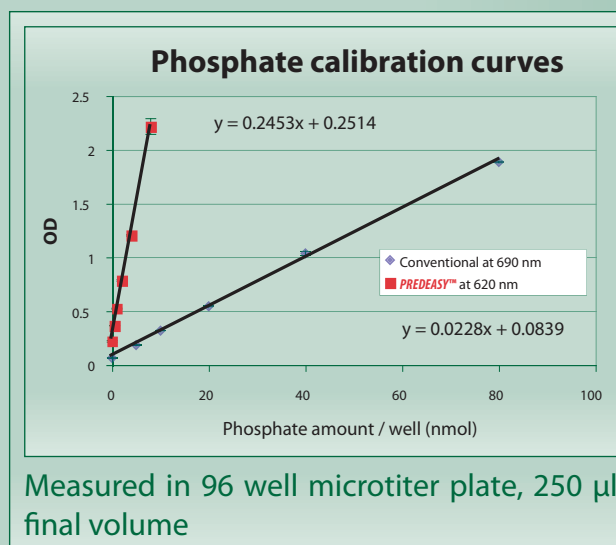
- Higher sensitivity
- More robust
- Easy to use
- More reliable
- All-in-one
- Easier automation, Higher Throughput
- Lower cost for each Data Point generated



## Benefits of PREDEASY™ ATPase Assay Kit over the conventional method

### Higher sensitivity

Ten times more sensitive phosphate detection. Shorter incubation times.



### More robust

Color stability of phosphate detection lasts **more than 3 hours**.

