

## **HTS solutions for protein kinases**

Jason Brown, Ph.D.

*Upstate Biotech*

While significant improvements in automation and miniaturization technologies have – to a degree – increased screening throughput, many companies face a bottleneck at the assay development and the subsequent candidate-profiling phases. Technologies and services that can facilitate these processes for protein kinases, the second largest target group for HTS labs after GPCRs, will be discussed.

The IMAP fluorescence polarization protein kinase kits provide a homogeneous HTS compatible solution that circumvents the assay development step. Furthermore all assays in the IMAP range follow very similar protocols thus further reducing tool-up and familiarisation time from kit delivery to screen. The utility of IMAP in HTS configurations will be discussed with reference to specific examples. In addition a new fluorescence ELISA assay for the SapKs will be presented.

Finally the application of KinaseProfiler™ for accelerating candidate-profiling by screening hits against a panel of 60 kinases will be reviewed.