

SLAS Tutorials & Posters

Tutorials

“Screening Epigenetic Targets with the Transcreener® HTS Assay Platform”
February 6, 2 – 2:50 pm

“Transcreener® HTS Enzyme Assay Development Service”
February 7, 8 – 8:50 am

Posters

- Interrogating GTPases, GAPs and GEFs with the Transcreener GDP assay , MP113
- Validation of Transcreener® EPIGEN Assays for screening chromatin modifying enzymes: Methylation and Acetylation, MP115
- An Innovative Multiwell Plate that Enables Immunocytochemistry with Non-Adherent Cells, TP148
- Validation of Transcreener® Fluorescent Polarization Assays using BMG LABTECH`s PHERAstar PLUS , TP132
- Automation of a Generic Fluorescence Methyltransferase Activity Assay , MP77

Discover more about enzyme detection and HTS at Booth 128!!

