

DAY 2

Agenda- Wednesday 12th

Redefining Newborn Screening Practices

Start time: 1:00 pm EEST

Improving the Lab Efficiencies

- 1:00 pm | Effective and Efficient Newborn Screening for X-linked Adrenoleukodystrophy:
One Extraction with Two Injections
Mei Baker, USA
- 1:20 pm | Tennessee's Progress in Implementing Steroid Profiling for Congenital Adrenal Hyperplasia
Lauren Tyler, USA
- 1:40 pm | Transitioning from Lab Developed Tests to NeoBase 2
Sarah Smith, UK
- 2:00 pm | Software Integration and Workflow Improvements in the NBS Laboratory
Brendan Reilly, USA
- 2:20 pm | Q&A
- 2:50 pm | Coffee Break (30 min)

DAY 2

Agenda- Wednesday 12th

Redefining Newborn Screening Practices

New Disorders in Consideration

- 3:20 pm | Newborn Screening for MLD: Methodology and Results of Pilot Studies Across 4 Countries
Mike Gelb, USA
- 3:40 pm | MLD: a Treatable Disease, the Perspective of a Neuropediatrician
Francesca Fumagalli, Italy
- 4:00 pm | Michigan's Experience Adding Guanidinoacetate and Creatine for GAMT Deficiency to the NeoBase™ 2 Assay
Krystyna Weiss-Pawlak, USA
- 4:20 pm | Implementation of MPS II Testing in a High Throughput Newborn Screening Laboratory
Patricia Ryland
- 4:40 pm | Q&A

End time: 5:10 pm EEST