

DAY 3

Agenda- Thursday 13th

The Future Shape of Neonatal Screening

Start time: 8:30 am EEST

World Programs Expanding the Menu to SCID, SMA, and LSDs

- 8:30 am | Building Tomorrow: The Professional Journey of an SMA Scientist in the Newborn Screening Era
Vanessa Romanelli, Brazil
- 8:50 am | Implementation of the Eonis Q System for Newborn Screening for Severe Combined Immunodeficiency and Spinal Muscular Atrophy in Catalonia
Ana Argudo-Ramirez, Spain
- 9:10 am | Thailand Government National Mandate: Expanding to Cover IEMs and SMA
Khunton Wichajarn, Thailand
- 9:30 am | National Implementation of LSD NBS in Korea
Sung-Eun Cho, Korea
- 10:00 am | Q&A

Next Potential Expansion Candidates on Molecular NBS

- 11:00 am | The Case for Universal CMV Screening: How Minnesota Became the First State in the USA to Screen All Newborns for Congenital CMV Infection
Mark Schleiss, USA
- 11:20 am | One Year of Universal Screening for Cytomegalovirus in Minnesota
Carrie Wolf, USA
- 11:40 am | Toward Prospective Newborn Screening for Angelman, Prader-Willi, and Dup15q Syndromes; Validation of a DNA Methylation-Based Test
Kate Kucera, USA
- 12:00 pm | Q&A

DAY 3

Agenda- Thursday 13th

The Future Shape of Neonatal Screening

NBS Genomics

- 1:30 pm | Principal Clinician Newborn Genomes Programme
David Bick, UK
- 1:50 pm | Experience from a United States Statewide Newborn Sequencing Study: Early Check
Holly Peay, USA
- 2:10 pm | Improving Screening Performance with 2nd Tier NGS - Expanding newborn screening towards genomics
Asbjørg Stray-Pedersen, Norway
- 2:30 pm | Considerations of Ethics and Equity When Using Genomics in Newborns
Amy Gaviglio, USA
- 3:20 pm | Round Table: NBS Genomics
Panelists: *Speakers*
Moderator: *Madhuri Hegde*
- 4:05 pm | Closing Remarks
Prahlad Singh, CEO, USA

End time: 4:05 pm EEST