

DAY 1

Agenda- Tuesday 11th

Activating Newborn Screening Around the World

Start time: 9:00 am EEST

9:00 am Welcome & Opening Remarks

Petra Furu, Finland

9:10 am Greetings from International Society for Neonatal Screening

Jim Bonham, UK

9:30 am WHO Progress on Maternal and Newborn Health Globally

Anshu Banerjee, WHO, Switzerland

9:50 am MLD Family Journey 'Nala and Teddi'

Ally Wallace, UK

Emerging Programs in LMICs

10:50 am Convince Governments to Act

Carmencita Padilla, Philippines

11:10 am Partnerships and Advocacy to Address the Sickle Cell Funding Paradox

Hon. Jane Ruth Aceng, Minister of Health, Uganda

11:30 am Leveraging HIV DBS Infrastructure for NBS in Uganda

Charles Kiyaga, Uganda

11:50 am Unitaid: Saving Lives Through Health Innovation

Kelsey Barrett, Switzerland

1:30 pm Round Table: How to Accelerate Access of Newborn Health innovations in LMICs?

Panelists: Hon. Jane Ruth Aceng (MOH, Uganda), Anshu Banerjee (WHO),

Carmencita Padilla (Philippines), Kelsey Barrett (UNITAID)

Moderator: Marika Kase

End time: 2:15 pm EEST