

11th of May 2016

Laboratory Training on Transporter Assays

FINAL PROGRAM

Meeting place and time: In front of the Continental Hotel Budapest, 42-44 Dohány street, 1074 Budapest, at 9:15 AM.

10:00 Registration and Coffee break

10:30-11:30 **Presentation:**

Timing and application of transporter assays during drug development through case studies

Herédi Szabó Krisztina, Head of Service Operations, SOLVO Biotechnology

11:30-14:30 **Laboratory demonstration and lunch** in groups

*Vesicular Transporter Assay: *Emese Kis, Head of Production, SOLVO Biotechnology*

*Uptake Transporter Assay: *Beáta Tóth, Scientist, SOLVO Biotechnology*

*Transwell Assay: *Károly Mogyorósi, Study Director, SOLVO Biotechnology*

*Sandwich cultured hepatocytes: *Erzsébet Beéry, Head of Laboratory, SOLVO Biotechnology*

*HepatoPac tox: *Zsuzsanna Gáborik, Scientific Development Manager, SOLVO Biotechnology;*

Jack McGeehan, Vice President of Operations, Ascendance

Coffe Break

14:30-15:30 Data evaluation and interpretation

Anna Klukovits, Study Manager, SOLVO Biotechnology

15:30- Transportation back to Continental Hotel Budapest