

2020 HEPATOPAC® User Group Meeting
Thursday, October 22, 2020

Time	Presentation
10:00 AM EST	Opening Remarks <i>Speaker: Chris Black, Ph.D., Senior Vice President, ADME-Tox, BioIVT</i>
10:05 AM EST	HEPATOPAC: Past, Present, and Future Applications in Drug Development <i>Speaker: Dr. Salman Khetani, Associate Professor, University of Illinois at Chicago</i>
11:00 AM EST	Development and Application of a Transcriptomic Signature of Bioactivation in the HEPATOPAC Model to Reduce Drug-induced Liver Injury Risk Early in Small Molecule Drug Development <i>Speaker: Wen Kang, Ph.D., Senior Principal Scientist, Merck</i>
12:00 PM EST	HEPATOPAC use at Boehringer Ingelheim Pharmaceuticals: A Decade of Learning <i>Speaker: Tom Chan, Ph.D., Senior Principal Scientist, Boehringer Ingelheim Pharmaceuticals</i>
1:00 PM EST	Live cell imaging of HEPATOPAC to characterize LNP delivery and mRNA expression to Hepatocytes <i>Speaker: Joe Sarkis, Ph.D., Principal Scientist, Moderna Therapeutics</i>
1:40 PM EST	Incorporation of Kupffer Cells in Long Term Hepatocyte Cultures in the HEPATOMUNE Model <i>Speaker: Scott Heyward, M.S., Director, R&D, Scientific Affairs, BioIVT</i>
2:20 PM EST	Closing Remarks <i>Speaker: Scott Heyward, M.S.</i>