

NEXUS 2022

INFORMATICS EU 2022 USER CONFERENCE



October 11th

10:00 – 11:30	Workshops (W1 & W2 topics will be presented) W1: Automating Workflows using the Signals Platform W2: Readily Expand your Signals Notebook Implementation into Validated Workspaces
Noon	Registration & Lunch
13:00	Meeting start
	Keynotes
13:15 – 13:45	Keynote: The PerkinElmer Informatics Signals Platform for Research Development and Clinical: The Power of the Suite David Gosalvez <i>Executive Director, Strategy & Informatics Portfolio, PerkinElmer Informatics</i>
13:45 – 14:30	Keynote: Digital Transformation across the R&D Value Chain David Drake <i>Drug Discovery IT Capability Lead, AstraZeneca</i>
14:30 – 15:00	Lightening Intro to Exhibition
15:00 – 15:30	Break & Exhibition
15:30 – 16:15	Keynote: Analytics in Time and Space Michael O'Connell, <i>Chief Analytics Officer, TIBCO</i>
16:15 – 16:55	Scaling Global Adoption of Signals Notebook Jen Kraljevich <i>Product Manager - Research Laboratory Methods Management Informatics - Science & Research F. Hoffmann-La Roche Ltd</i>
16:55 – 17:30	Compound-Substance-Product. A New Structure to Our Data at AZ Wolfgang Klute <i>Associate Director, Pharmaceutical Sciences, AstraZeneca</i>
17:30 – 17:45	Day Wrap up
18:00 – 22:00	Evening Social Networking Event at the Universeum

October 12th			
9:00 – 10:30	Workshops (W1 –W4 topics will be presented) W1: Automating Workflows using the Signals Platform W2: Readily Expand your Signals Notebook Implementation into Validated Workspaces W3: Streamlining Formulation Research in the Cloud W4: Three Enhancements for Medical Review within Spotfire		
10:30 – 11:00	Break & Exhibition		
	Research Track	Clinical Track	ELN Track
11:00 – 11:30	There and Back again - a Signals Research Suite Walk-through from Material Creation through <i>In Vitro</i> assay and Back to SAR Analysis <i>Dan Weaver, Sr Product Leader Screening & Lead Discovery PerkinElmer Informatics</i>	Enhancing Clinical Quality-by-Design (QbD) with Analytics <i>Muneeb Kahn Clinical Trial Spotfire Analytics Lead, PerkinElmer Informatics</i>	The Future of Signals Notebook: Partnering Together <i>Josh Bond Sr. Director, Product Portfolio PerkinElmer Informatics</i>
11:30 – 12:00	Streamlined Data Integration of CETSA Data at Pelago Bioscience <i>Rikard Alm, PhD Senior Scientist, Pelago Biosciences</i>	Identifying Anomalies and Outliers in Clinical Trial Data <i>Pekka Tiikkainen Principal Clinical Data Scientist, Bayer Pharmaceuticals</i>	R&D Digitalization Through Signals Notebook Implementation at Novéal <i>Pierrick Ondet, PhD Senior Scientist & ELN Project Leader Novéal</i>
12:00 – 12:30	Image Data Management in the Context of New Applications Like Cell Painting, AI/ML Derived Knowledge <i>Dr. Karsten Kottig Vice President, Imaging & Analysis Evotec SE</i>	Adding Value to Clinical Data: Perspectives from the Implementation of FAIR Principles <i>Giovanni Nisato Project Manager / FAIR implementation Pistoia Alliance</i>	First Impact of SNB Rollout and Beyond <i>Andreas Muheim Head of Research Information Systems Givaudan</i>
12:30 – 13:45	Lunch		
13:45 – 14:15	Selection of De Novo Molecules from REINVENT with Spotfire - from Large Scale Property Analysis to Synthesis Decisions <i>Eva Nittinger Associate Principle Scientist in Computational Chemistry AstraZeneca</i>	Visual & Integrated Support of Translational Analysis (VISTA) – to Shed Some Light on the Support Concept Behind <i>Ute Burkard Global Team Lead Descriptive Analysis Boehringer Ingelheim</i>	Capturing Data, a lot of Data, for a Quality by Design Approach to Product Development <i>Michele Ghirardi Formulation Development Manager Custom Pharma Services</i>
14:15 – 14:45	Use of Spotfire in the Development of Lipid Nanoparticles for Vaccine and Targeted Genome Editing <i>Michael Munson, PhD Senior Research Scientist AstraZeneca</i>	Industry Presentation	Experiences with the Introduction of Signals Notebook, Present Use and Future Ventures at Ajinomoto Bio-pharma Services <i>Stan Coene Lab Technician Process R&D Ajinomoto Bio Pharma Services</i>
14:45 – 15:00	Break & Exhibition		
15:00 – 15:30	TIBCO Spotfire What's New and What's Next <i>Niklas Amberntsson, Director, Spotfire Product Management TIBCO</i>		
15:30 – 16:00	Spotfire as an Enabler for AI-Augmented Drug Discovery and Development <i>Johan Ulander, Principal Scientist, Data Science and Modelling Justin Morley, Cheminformatics Data Scientist AstraZeneca</i>		
16:00 – 16:30	Signal Inventa's Cutting Edge: Integrated Matched Molecular Pair and Reaction Pair Analysis Reveal Complex Structure-activity Relationships <i>Dan Weaver, Sr Product Leader, Screening & Lead Discovery PerkinElmer Informatics</i>		
16:30 – 16:45	Conference Wrap & Close		

REGISTER / LOG IN NOW

<https://info.perkinelmer.com/INF-GNRL-NEXUS-Europe-2022-Conference-Register>

PerkinElmer, Inc.
940 Winter Street
Waltham, MA
02451 USAP:
(800) 762 –4000
or
(+1) 203 –925 –4602
<https://perkinelmerinformatics.com/>

