



April 18

9:00	Registration
9:30	Welcome Kapil Aggarwal, Director of Software Sales, Revvity Signals
9:45	Welcome to Revvity Signals NEXUS India 2024 Ralph Dioguardi, VP Software Sales, Revvity Signals
10:00	Keynote - Signals of Success - Empowering Scientific Innovation with Revvity Signals Platform David Gosalvez, Executive Director, Strategy & Informatics Portfolio, Revvity Signals
10:30	Collaborative Platforms: Enabling Seamless Communication & Knowledge Sharing in Research Atanu Roy, President & Group CIO, PI Industries
11:00	Streamlining Formulation R&D Using Revvity Signals Software Dr. Manish Khandagale, Senior Field Application Scientist, Revvity Signals
11:30	Tea / Coffee Break
11:45	Accelerating Laboratory Digital Transformation: Integration Signals Notebook with Scitara's DLX Technology Geoff Gerhardt, CTO, Scitara
12:15	Panel Discussion: Addressing Challenges in the Lab of the Future <ul style="list-style-type: none"> ➤ David Gosalvez, Executive Director, Strategy & Informatics Portfolio, Revvity Signals(Chair) ➤ Dr Chander Shekhar Devra, CIO, SRF ➤ Dr. Kamlesh Padiya, Director, Synvision Chem ➤ Yogendra Yadav, Director-IT, Sai LifeSciences ➤ Geoff Gerhardt, CTO, Scitara
13:00	Lunch
14:00	Navigating Our Digital Journey to Signals Notebook at Givaudan Andreas Muheim, Head of Research Information Systems, Givaudan
14:30	Signals ChemDraw: Taking Chemistry Communication to the Next Level Pierre Morieux, Ph.D., Chemistry Product Marketing Manager and ChemDraw Wizard, Revvity Signals
15:00	Industry Case Study: Aarti Industries Limited Vinod Shivanand Honmure, R&D Excellence Head, Aarti Industries Limited
15:30	Tea / Coffee Break

16:00	Value Generation through Digital: Mastering the Art of Harmonizing Business Objectives with Tech Initiatives Tushar Zade, Chief Information Officer, Aurigene Pharmaceutical Ltd.
16:30	Unleashing the Potential of Signal Notebook for Biological Workflows Dr. Tanvira Paul, Field Application Scientist, Revvity Signals
17:00	Bridging Efficiency and Compliance: The Expectations of the Generic Pharma Industry from Electronic Lab Notebooks Dr.Gopal Vaidyanathan, VP-Analytical R&D, Natco Pharma Ltd
17:30	Panel Discussion: The Role of AI in Digital Transformation <ul style="list-style-type: none"> ➤ Pierre Morieux, Chemistry Product Marketing Manager and ChemDraw Wizard, Revvity Signals (Chair) ➤ Rupali Aher, Head, R&D Digital Excellence, Cipla ➤ Ankit Aggarwal, Director, R&D IT, PI Industries ➤ Dr. Muthukumarasamy Karthikeyan, Chief Scientist, CSIR, National Chemical Laboratory ➤ Tushar Zade, CIO, Aurigene Pharmaceuticals
18:15	Conclude Kapil Aggarwal, Director Software Sales, Revvity Signals
18:30	Cocktails Followed by Dinner

REGISTER NOW

<https://info.revvity.com/NEXUS-India-2024-User-Conference-Registration>

Revvity Signals Software, Inc.
940 Winter Street
Waltham, MA
02451 USA
P: (800) 762-4000
or
(+1) 203-925-4602
<https://revvitysignals.com/>

revvity
signals

For a complete listing of our global offices, visit www.revvity.com/contact-us

Copyright ©2024, Revvity Signals Software, Inc. All rights reserved.