

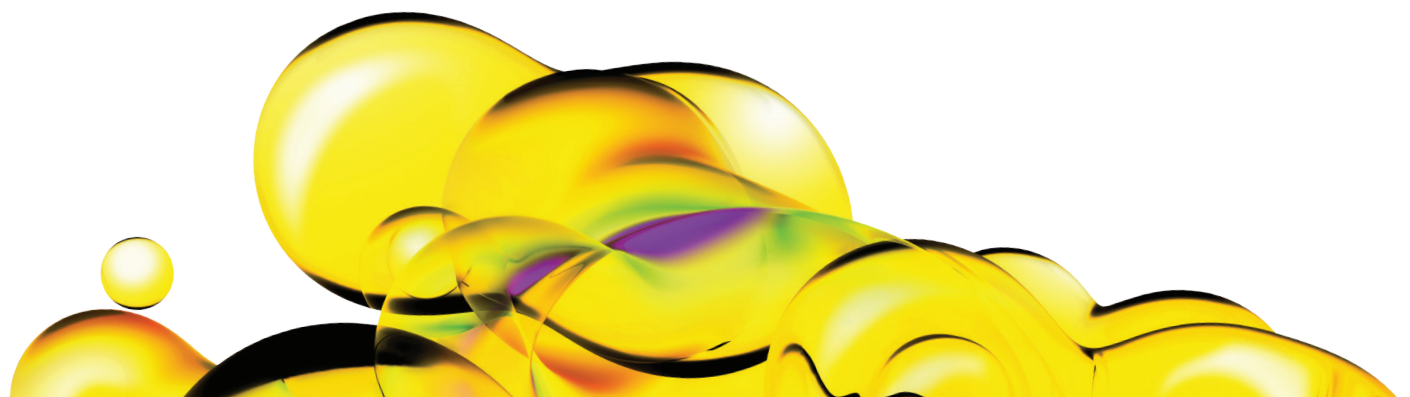
📅 Friday, 24th April 2026

📍 The Heritage Resort, Manesar

Reporting Time: 09:00 am to 10:30 am
(Registrations & Kits Distribution)

AGENDA

Time	Sessions	Speakers
10:30 am - 10:50 am	The Next Growth Chapter: Revvity's Vision for India & South Asia	Shripad Joshi <i>President India, South Asia, Thailand and Vietnam</i>
10:50 am - 11:10 am	Life Sciences & Applied Genomics: Unlocking the Next Wave of Discovery	Mayank Srivastava <i>Director - Life Sciences and Applied Genomics</i>
11:10 am - 11:30 am	Transforming Diagnostics: Innovations That Impact Lives	Chetan Patil <i>General Manager Sales - Reproductive Health</i>
11:30 am - 11:40 am	Growing Together: Financial Strength for Mutual Success	Alok Kumar <i>Finance Controller & Director</i>
11:40 am - 11:50 am	Built on Trust: Strengthening Compliance for Sustainable Growth	Suvajit Chakraborty <i>Associate General Counsel & Legal Director</i>
11:50 am - 12:00 pm	Delivering Excellence: Creating a Stronger Operations Ecosystem	Srinivas Rao K <i>General Manager - Customer Operations</i>
12:00 pm - 12:15 pm	Simplifying Success: Channel Policies	Hardik Sheth <i>Channel Operations Manager</i>
12:15 pm - 12:30 pm	Partners for Purpose: Strategic Marketing Vision	Himanshu Vashishth <i>Marketing Head</i>
12:30 pm - 01:00 pm	Q&A	
01:00 pm - 02:30 pm	Lunch & Networking	
02:30 pm - 04:00 pm	Breakout Sessions: Innovation Circle	
04:00 pm - 06:00 pm	Serenity Walk & High Tea	
07:00 pm - 09:00 pm	Celebrating Excellence Night	
09:00 pm onwards	Dinner and Cocktails	



📅 Friday, 24th April 2026

📍 The Heritage Resort, Manesar

Reporting Time: 09:00 am to 10:30 am



Vision 2026: The road ahead

Step into the future with an exclusive deep dive into Revvity's Vision 2026—where we are headed, how the market is evolving, and the role you will play in shaping that journey.



Strategic projects & market drivers

Discover the key growth engines—emerging technologies, funding trends, and market drivers—that will define success in the coming years.



How we win-together

A focused session on winning strategies, co-creation models, and how we can strengthen execution on the ground to accelerate mutual growth.



Fun, food & networking

An unforgettable evening filled with entertainment, music, and meaningful conversations, strengthening bonds beyond business.



Awards night

Recognizing excellence, performance, and partnership—celebrating those who have truly made an impact.

Important notes

- Please confirm your participation in advance to ensure seamless arrangements.
- Travel and accommodation details will be shared separately.
- Please adhere to the event schedule to make the most of each session.
- Active participation is encouraged—this is a collaborative platform.
- Business formals for the day & traditional evening attire for the awards night.
- Room bookings will be arranged based on your confirmation, and the shared details will be used for your delegate badge. We request you to please share your information at the earliest.

For any enquiries, please reach out to:

Himanshu Vashishth

☎ 96433 29402

✉ himanshu.vashishth@revvity.com

Anushka Sharma

☎ 77108 57343

✉ anushka.sharma@revvity.com