

International Women's Day



revvity

Editorial

When science meets strength

Celebrating the women shaping healthier futures

Every breakthrough has a story.

Sometimes it begins in a laboratory — under the quiet hum of instruments, where a scientist studies patterns invisible to the human eye. Sometimes it begins in a neonatal ward, where a clinician makes a timely decision that changes the course of a newborn's life. Sometimes it begins in a research center, where a question refuses to be ignored.

This International Women's Day, at Revvity, we celebrate the women who turn these beginnings into impact.

Across laboratories, hospitals, universities, and public health programs, women scientists and clinicians are reshaping the future of healthcare. They are decoding complex biology, refining diagnostics, advancing cancer research, strengthening newborn screening, and building systems that reach the most vulnerable communities. But beyond their titles and designations, what truly defines them is purpose.

Purpose to detect disease earlier.

Purpose to treat more precisely.

Purpose to ensure that no patient is unseen and no possibility unexplored.



Behind every dataset is dedication. Behind every clinical protocol is compassion. Behind every published paper is perseverance. Many of these journeys have required navigating structural barriers, limited resources, and the subtle expectations placed on women in demanding professions. Yet what stands out is not the difficulty — it is the determination.

We see women leading diagnostic laboratories with precision and integrity. We see researchers mentoring young minds, encouraging curiosity where doubt once lived. We see clinicians balancing evidence with empathy, science with humanity.

Their leadership is not loud — it is impactful. It transforms institutions, inspires teams, and builds cultures rooted in excellence.

In India and around the world, women are accelerating the adoption of advanced technologies, integrating AI into research, strengthening quality standards, and expanding access to critical diagnostics. They are not waiting for the future to arrive — they are designing it.

At Revvity, we are privileged to collaborate with these changemakers. Our mission has always been to enable earlier insights and smarter decisions in healthcare. But technology alone does not create transformation. People do. And time and again, it is the resilience, clarity, and vision of women in science that convert potential into progress.

When a newborn receives an early diagnosis that safeguards their life, when a cancer patient benefits from more precise monitoring, when a research breakthrough opens a new therapeutic pathway — we are reminded that impact is deeply human.

This International Women's Day, we honor the women who bring both rigor

and heart to science. The ones who ask harder questions. The ones who stay curious. The ones who lead with integrity. The ones who quietly, consistently change the trajectory of lives.

Your work is not just advancing science.
It is strengthening societies.
It is building healthier generations.

And it is inspiring the next young girl who will step into a laboratory one day believing that she belongs there.

At Revvity, we celebrate you — not only today, but every day we work alongside you to create a future where discovery knows no limits.

Happy International Women's Day.



Dr. Anubama Rajan

Organoid Technologies & Translational Research
IIT Madras



Background

Assistant Professor in the Department of Medical Sciences & Technology at IIT Madras. Ph.D. and postdoctoral training from Baylor College of Medicine (USA), where she pioneered organoid (“mini-organ”) technology for disease modeling. Her lab (Orgamed) develops patient-derived stem-cell organoids to model human diseases.

Key Contributions

Dr. Rajan’s team uses organoids as surrogates for human subjects, bridging bench research and clinical testing. She was awarded the prestigious Wellcome Trust-India Alliance Early Career Fellowship (2023) and a DBT Ramalingaswami Re-entry Fellowship (2022) to establish her translational program. Her work has already identified novel biomarkers and therapeutic targets for infectious diseases.

Dr. Rajan’s journey shows how cutting-edge stem-cell science can be harnessed for real-world impact. As a young leader in bioengineering and microbiology, she exemplifies the power of curiosity and persistence. Her success inspires scientists to push the boundaries of technology — proving that even “mini-organs” on a dish can make a big difference in early diagnosis and personalized treatment.

Inspiration

Dr. Rajan’s work demonstrates how innovative thinking and scientific perseverance can transform fundamental research into meaningful healthcare solutions. By advancing organoid technology and translational science, she is helping redefine how diseases are studied and treated. Her journey reflects the power of curiosity, resilience, and interdisciplinary collaboration in modern biomedical research. As a young leader in bioengineering and microbiology, she inspires emerging scientists—especially women in science—to pursue bold ideas and push the boundaries of technology in order to create impactful solutions for global health challenges.

Dr. Sherine Sonia Cubelio

Marine Biodiversity & Genetic Barcoding

CMLRE, Kochi



Background

Scientist F at India's Centre for Marine Living Resources & Ecology (CMLRE), under the Ministry of Earth Sciences. Holds MFSc and Ph.D with specialization in marine biosciences.

Key Contributions

Dr. Cubelio leads research on deep-sea biology and the Census of Marine Life. She applies DNA barcoding to catalog ocean organisms, identifying species from the deep ocean. Her projects include large-scale surveys and genetic profiling of marine organisms to improve environmental diagnostics and resource management.

Dr. Cubelio's career reminds us that diagnostics isn't limited to hospitals - it extends to our planet. By mapping ocean biodiversity, she strengthens our understanding of marine ecosystems and climate change. Her passion for exploring uncharted waters - scientific and literal - encourages other women to dive into STEM fields where there are few role models. She proves that dedicated research in extreme environments can yield discoveries with global impact.

Inspiration

Dr. Sherine Sonia Cubelio's work highlights how scientific exploration can expand our understanding of the natural world and its impact on human life. By studying deep-sea biodiversity and applying advanced genetic tools such as DNA barcoding, she contributes to protecting marine ecosystems and improving environmental knowledge. Her journey reflects courage, curiosity, and dedication to discovery, inspiring young scientists—especially women in STEM—to explore new scientific frontiers and contribute to research that benefits both science and the planet.

Dr. Anupama Kumari

M.V.Sc., B.V.Sc. & AH

Veterinary Translational Scientist
Laboratory Animal Science Leader



Background

Dr. Anupama Kumari is a Veterinary Translational Scientist and Laboratory Animal Science specialist with over a decade of experience in preclinical research and translational hepatology. She currently serves as Veterinary Officer at the Centre for Comparative Medicine, Institute of Liver and Biliary Sciences (ILBS), New Delhi, where she works on integrating advanced animal models with clinical hepatology research to better understand liver diseases and therapeutic strategies.

Key Contributions

Dr. Kumari's research focuses on the mechanisms of infectious, metabolic, and immune-mediated liver diseases using advanced preclinical models. She has served as Principal Investigator and Co-Principal Investigator on several nationally funded projects supported by agencies such as DBT, BIRAC, DST, and SERB. Her work spans areas including cancer immunotherapy, hepatitis virus pathogenesis, metabolic liver diseases, and translational drug discovery. In recognition of her scientific contributions, she received the ICLAS-AFLAS Youth Progress Award (2025) from the International Council for Laboratory Animal Science and the Asia Federation of Laboratory Animal Science Associations.

Inspiration

Dr. Kumari's journey reflects the powerful role of translational research in bridging laboratory science with clinical solutions. Her dedication to scientific discovery, mentorship, and capacity building in laboratory animal sciences highlights the importance of collaborative research in addressing complex health challenges. Her achievements and international recognition continue to inspire young scientists—especially women in biomedical research—to pursue innovation and contribute to advancing global healthcare.

Dr. Akila Prashant

Professor & Head

Department of Biochemistry, JSS Medical College
JSS Academy of Higher Education and Research
Molecular Diagnostics, Genomics & Translational
Laboratory Medicine



Background

Dr. Akila Prashant is Professor and Head of Biochemistry at JSS Medical College and Hospital, JSS Academy of Higher Education & Research, Mysuru. With training in medicine, biochemistry, cancer biology, and genomic medicine, her work bridges clinical laboratory diagnostics and molecular research, focusing on rare genetic disorders, cancer biology, and precision medicine.

Key Contributions

Dr. Prashant has played a pivotal role in strengthening advanced diagnostics and translational research within academic healthcare settings. She has been instrumental in establishing molecular biology, cytogenetics, and genomic diagnostic facilities, enabling clinicians to integrate genetic insights into routine patient care. Her work spans molecular biomarkers, inflammatory pathways, and genomic approaches to complex diseases including cancer, sepsis, and rare genetic disorders.

With over 200 peer-reviewed publications, including contributions to global collaborations such as the Global Burden of Disease studies, her research has appeared in leading journals including The Lancet. Beyond research, she is deeply committed to mentorship and academic leadership, guiding doctoral scholars and early-career researchers while fostering collaborative, interdisciplinary scientific environments.

Inspiration

Dr. Prashant's journey reflects the power of curiosity-driven science combined with clinical purpose. By advancing molecular diagnostics and nurturing future scientists, she continues to help shape a healthcare landscape where earlier diagnosis, precision medicine, and scientific collaboration lead to healthier generations.

Dr. K. Aparna Sharma

Maternal-Fetal Medicine

AIIMS, New Delhi



Background

Professor, Department of Obstetrics & Gynaecology, All India Institute of Medical Sciences (AIIMS), New Delhi. Dr. K. Aparna Sharma is a leading expert in high-risk pregnancy and fetal imaging.

Key Contributions

She heads the Maternal-Fetal Medicine division at AIIMS, overseeing diagnostics for complex pregnancies. Her clinical and research work involves advanced obstetric ultrasound, prenatal interventions (e.g. amniocentesis, fetal therapy), and maternal health innovation. She trains fellows in fetal medicine and has developed protocols for managing twin pregnancies and birth defects. As faculty, she also contributes to national guidelines in obstetric care.

Dr. Aparna Sharma exemplifies how women can lead in life-saving diagnostics. By managing high-risk pregnancies and pioneering fetal therapies, she helps bring healthy babies into the world. Her leadership at India's premier medical institute encourages women to enter and excel in obstetrics and gynecology research. Her commitment to patient-centered care inspires the next generation of female clinicians to advance maternal and child health.

Inspiration

Dr. K. Aparna Sharma's work reflects the profound impact that expertise and compassion can have in shaping maternal and fetal healthcare. By advancing diagnostic techniques and pioneering interventions for high-risk pregnancies, she contributes to safeguarding the health of both mothers and unborn children. Her leadership in maternal-fetal medicine demonstrates how dedication to science and patient care can transform outcomes for families. Through her achievements and mentorship, she inspires the next generation of clinicians and researchers—especially women in medicine—to pursue excellence in advancing maternal and child health.

Dr. Manisha Kumar

Fetal Medicine & Genetics

Lady Hardinge Medical College, Delhi



Background

Professor of Obstetrics & Gynecology and head of Fetal Medicine at Lady Hardinge Medical College (LHMC), Delhi. Chairperson of the AOGD fetal medicine subcommittee.

Key Contributions

Dr. Kumar leads prenatal diagnostic services at LHMC. Her interests include antenatal ultrasound, fetal autopsy and genetic counseling. She has won several awards (e.g. FOGSI Genetic & Fetal Medicine Award, Rajat Ray Award) for her contributions. Dr. Kumar has been Principal Investigator on major projects (ICMR preeclampsia screening, WHO stillbirth initiative) and established the NIDAN genetic lab at LHMC. She has over 80 publications in indexed journals.

Dr. Kumar's work highlights the importance of research-driven policy in prenatal care. By integrating genetics and ethics into obstetrics, she improves how India screens and treats fetal disorders. Her active role in national projects and publications encourages women researchers to take on leadership in public health. She shows that clinical excellence paired with scientific vision can transform diagnostic protocols for mothers and babies.

Inspiration

Dr. Manisha Kumar's work reflects the power of combining clinical expertise with research-driven innovation to improve maternal and fetal healthcare. By advancing prenatal diagnostics and integrating genetics into obstetric practice, she contributes to more informed and ethical approaches to managing fetal conditions. Her leadership in national research initiatives and academic contributions highlights the importance of evidence-based medicine in shaping healthcare policies. Through her dedication to science, education, and patient care, she inspires the next generation of clinicians and researchers—especially women—to lead advancements in fetal medicine and public health.

Dr. Mabi Jaleel

Drug Discovery & Translational Biology

Sai Life Sciences, Hyderabad



Background

Vice President and Head of Biology (India) at Sai Life Sciences, Hyderabad – a leading drug discovery services company. Ph.D. in Biochemistry from the Indian Institute of Science; postdoctoral fellowship in drug signalling pathways at the University of Dundee.

Key Contributions

Dr. Mabi Jaleel brings over two decades of experience in Drug Discovery R&D, with a career that spans academia, biotech, and global CRO leadership. She began her scientific journey at the Indian Institute of Science (IISc), after which she held senior research positions at Dr. Reddy's Laboratories and Aurigene Discovery Technologies.

During her tenure at Aurigene, Dr. Jaleel served as Project Lead, driving multiple end-to-end drug discovery programs. Under her leadership, seven preclinical candidates successfully advanced into clinical development, marking a significant contribution to the company's therapeutic pipeline.

In 2020, Dr. Jaleel joined Sai Life Sciences, and by 2024 progressed to the role of Vice President, Biology. She currently oversees the *in vitro*, *in vivo*, and translational biology groups, enabling global pharma and biotech partners to advance their programs from bench to clinic. Her work spans diverse therapeutic areas, including Oncology, Immuno-oncology, Inflammation & Immunology, CNS, and rare diseases.

Dr. Jaleel's career exemplifies the profound impact women bring to the biotech and CRO sectors. Rising through competitive industrial research environments, she has demonstrated that Indian women can lead major drug discovery initiatives at the highest levels. Her leadership, mentorship, and cross-functional expertise continue to inspire the next generation of women scientists pursuing careers in industrial R&D.

By bridging foundational biological science with cutting-edge pharmaceutical innovation, Dr. Jaleel embodies the forward-looking scientific leadership essential for tackling the diseases of the future.

Dr. Kalpana Singh

Department of Biochemistry,
King George's Medical University (KGMU)



Background

Dr. Kalpana Singh is a respected academic and researcher in the Department of Biochemistry at King George's Medical University (KGMU). Her work spans clinical biochemistry and endocrinology, fields that play a critical role in understanding metabolic disorders and hormonal regulation in human health.

Contribution & Impact

Dr. Singh's research in clinical biochemistry and endocrinology contributes to improving the biochemical understanding of diseases such as metabolic disorders, hormonal imbalances, and endocrine-related conditions. Through her academic research and laboratory work, she supports the development of diagnostic insights that help clinicians better understand disease mechanisms and improve patient management. Her work also contributes to the training of medical students and researchers, strengthening the scientific foundations of clinical diagnostics.

Inspiration

Dr. Singh's dedication to scientific research and academic mentorship highlights the vital role of biochemistry in modern medicine. By advancing biochemical research and nurturing future scientists, she contributes to building a stronger scientific ecosystem that drives healthcare innovation. Her journey stands as an inspiration for aspiring researchers who seek to translate biochemical discoveries into meaningful improvements in human health.

Dr. Neerja Gupta

Clinical Geneticist,
Department of Pediatrics AIIMS, Delhi



Background

Dr. Neerja Gupta (MD Pediatrics) is a faculty in charge in the Division of Genetics, Department of Pediatrics at the All India Institute of Medical Sciences (AIIMS), New Delhi. A distinguished clinician-scientist in medical genetics, she has been recognized with prestigious honors including the Dr. I.C. Verma Outstanding Researcher Award (2024) and the Dr. I.C. Verma Excellence in Research Award for Young Pediatricians (2020).

Contribution & Impact

Dr. Gupta has played a significant role in advancing clinical genetics and rare disease research in India. She serves as the nodal officer for the Centre of Excellence on Rare Disorders at AIIMS under the Ministry of Health and Family Welfare and coordinates the WHO Birth Defect Surveillance Program across the SEAR region. Her contributions extend to global and national scientific initiatives, including work with the Undiagnosed Disease Network (NIH) and leadership roles in professional societies focused on medical genetics and metabolic disorders. With more than 200 scientific publications, her research has significantly advanced the understanding of inherited metabolic disorders, genomic conditions, and rare pediatric diseases.

Inspiration

Dr. Gupta's work reflects the transformative power of genetics in improving diagnosis and care for rare and complex diseases. Through her clinical leadership, research excellence, and dedication to medical education, she continues to shape the future of genomic medicine in India.

Dr. Nimmi Kansal

MBBS, MD, Clinical Chemistry from University
College of Medical Sciences, Delhi

Director for Clinical Chemistry & Biochemical
Genetics, Working with Dr LalPath Labs since 2008



Background

Dr. Nimmi Kansal is an accomplished professional in laboratory medicine and diagnostics, associated with Lal PathLabs, one of India's most respected diagnostic institutions. With extensive experience in clinical diagnostics and laboratory science, she has contributed significantly to strengthening the quality and reliability of diagnostic services that support clinicians in delivering accurate medical care.

Contribution & Impact

Through her work in diagnostic medicine, Dr. Kansal has played an important role in ensuring high standards of laboratory testing, quality assurance, and scientific rigor. Her efforts support the development of reliable diagnostic systems that enable early disease detection and precise clinical decision-making. By strengthening laboratory processes and promoting quality-driven diagnostic practices, she contributes to improving patient outcomes and advancing the broader healthcare ecosystem.

Inspiration

Dr. Nimmi Kansal's work reflects the critical role of laboratory medicine in enabling accurate diagnosis and effective patient care. Through her commitment to quality, precision, and scientific excellence in clinical chemistry and biochemical genetics, she contributes to strengthening diagnostic standards that clinicians rely on every day. Her leadership in advancing reliable laboratory practices highlights how strong diagnostic systems form the backbone of modern healthcare. Her journey inspires young professionals—especially women in laboratory science—to pursue excellence in diagnostics and contribute to building a more robust and reliable healthcare ecosystem.

Dr. Aarti Nagpal

Associate Director and Head –
Biochemistry & Hematology
Pathkind Diagnostics Pvt. Ltd.



Background

Dr. Aarti Nagpal is a respected leader within the diagnostics ecosystem and is associated with Pathkind Labs, where she contributes to expanding and strengthening modern diagnostic services. Her professional journey reflects a deep commitment to advancing laboratory medicine and improving healthcare accessibility through reliable diagnostic solutions.

Contribution & Impact

Through her work, Dr. Nagpal has helped strengthen diagnostic infrastructure and improve laboratory systems that support clinicians in making timely and informed medical decisions. By promoting high standards of diagnostic accuracy and expanding access to quality testing services, she contributes to bridging gaps in healthcare delivery and enabling better disease management across communities.

Inspiration

Dr. Aarti Nagpal's leadership reflects the growing influence of women in shaping the future of diagnostics and healthcare innovation. With professional interests spanning allergy diagnostics, prenatal screening, newborn screening, and flow cytometry, she contributes to advancing specialized diagnostic areas that are essential for early disease detection and preventive healthcare. Her journey demonstrates how scientific expertise combined with visionary leadership can strengthen healthcare systems and improve patient outcomes. Through her work and dedication, she continues to inspire young professionals—especially women in science and medicine—to pursue impactful careers that combine knowledge, innovation, and purpose.

Dr. Debarati Mazumder Tagore

Scientific Director,
Bristol Myers Squibb, India



Background

Dr. Debarati Mazumder Tagore is a Scientific Director at Bristol Myers Squibb, leading the In vitro and Molecular Pharmacology team. Her work sits at the intersection of mechanistic rigor and translational relevance, advancing discovery programs across oncology, immunology, neuroscience and cardiovascular diseases. Through innovative, physiologically relevant assay platforms, Dr. Mazumder Tagore and her team drive high quality datasets to enable decisions, accelerate lead optimization, and strengthen therapeutic hypotheses. A passionate advocate for Women in STEM, she is committed to inclusive leadership, mentorship, and cultivating a strong scientific community grounded in purpose and impact.

Contribution & Impact

Under Debarati's leadership, the In vitro and Mechanistic Pharmacology team has successfully balanced automation and throughput with translational relevance, ensuring that scientific rigor remains at the core of discovery. As a highly accomplished woman leader, she has not only driven scientific excellence but also set a powerful example of resilience, clarity, and conviction. Dr. Debarati has led multiple discovery programs end to end and enabled several clinical candidate nominations that have advanced the portfolio. Her collaborative leadership style fosters strong cross functional partnerships, translating deep mechanistic insights into actionable strategies—and ultimately, into potential medicines that aim to improve patients' lives.

Inspiration

Dr. Mazumder Tagore's professional journey reflects her passion and purpose in advancing medicines that make a meaningful difference for patients. She firmly believes that a strong understanding of the mechanistic underpinnings of disease biology is key to enabling faster, smarter, and higher quality transitions from discovery to the clinic. Guided by this philosophy, she invests deeply in developing future scientific leaders—providing guidance, advocacy, and encouragement to those who aspire to pair scientific excellence with empathy and authentic leadership.

Dr. Prathima Radhakrishnan

Director and Consultant,
Bangalore Fetal Medicine Centre



Background

Dr. Prathima Radhakrishnan is a fetal medicine specialist and Director and Consultant at the Bangalore Fetal Medicine Centre (BFMC) with over 16 years of experience in fetal medicine. She holds prestigious qualifications including FRCOG, Diploma in Fetal Medicine (FMF, UK), and Diploma in Advanced Obstetric Ultrasound (RCOG, UK) and received advanced training under Prof. Kypros Nicolaides and Prof. Lindsay Allen in fetal medicine and fetal echocardiography.

Contribution & Impact

Dr. Radhakrishnan is recognized for pioneering contributions to fetal diagnostics and therapy in India. She was the first Indian to receive the FMF (UK) Diploma in Fetal Medicine and to perform therapeutic fetoscopic laser surgery for complicated monochorionic twins. Her clinical focus includes early fetal screening, prenatal diagnosis, and fetal therapy, and she actively contributes to medical education through lectures, workshops, and research presentations. She has received honors including the Arnimaal Award and Moninder Kochar Award for her academic contributions.

Inspiration

Dr. Radhakrishnan's work demonstrates how advanced prenatal diagnostics and specialized care can transform outcomes for both mothers and unborn babies. Through her pioneering clinical work, research, and dedication to training the next generation of fetal medicine specialists, she continues to expand the possibilities of prenatal healthcare. Her journey inspires doctors—especially women in medicine—to pursue specialized fields like fetal medicine and to push the boundaries of innovation in maternal and fetal healthcare.

The inspiring women featured in this report represent only a small glimpse of the many remarkable scientists, doctors, and healthcare professionals who continue to shape the future of diagnostics and medicine. Across laboratories, hospitals, research institutions, and diagnostic centers, countless women are advancing science, improving patient care, and strengthening healthcare systems through their dedication and expertise. While this report highlights a few exemplary leaders, it also stands as a tribute to the many others whose contributions quietly power progress every day. On this International Women's Day, we celebrate and salute all women in science and healthcare whose passion, perseverance, and commitment continue to inspire the next generation and move the world of life science and diagnostics forward.





www.revivity.com

revvity

Copyright ©2026, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

For research use only. Not for use in diagnostic procedures.