



BECKMAN COULTER CARES AWARD



WE CARE THAT YOU CARE

PANEL OF JUDGES



Maurine M. Murtagh

CEO of the International Diagnostics Centre at the London School of Hygiene & Tropical Medicine

Maurine Murtagh is the CEO of the International Diagnostics Centre at the London School of Hygiene & Tropical Medicine. She is also a consultant specializing in diagnostics for global health in which capacity she has worked with the World Health Organization, the Bill & Melinda Gates Foundation and UNITAID, among others. Previously, Maurine was the Director of Diagnostic Services for the Clinton Health Access Initiative (CHAI), a position she held for almost five years. Prior to joining CHAI, Maurine was a corporate partner at the law firm of Howard Rice Nemerovski Canady Falk & Rabkin in San Francisco, California, where her focus was on mergers and acquisitions and securities law. She graduated from Georgetown University and holds an M.B.A. from the University of Texas at Austin, an M.Phil. in finance from Columbia University and a J.D. from Duke University.



Francois Venter

MD, FCP (SA), Mmed, DTM&H, Dip HIV Management (SA), Deputy Executive Director of the Wits Reproductive Health and HIV Institute at the University of the Witwatersrand.

Dr Venter is the Deputy Executive Director of the Wits Reproductive Health and HIV Institute at the University of the Witwatersrand. He leads multiple anti-retroviral treatment optimisation studies, and has an active interest in public sector access to HIV services. He is currently working on new first and second line antiretroviral options, patient linkage to care interventions, and self-testing projects. Previously, he led large PEPFAR-funded HIV programmes in South Africa, including one that focused on truckers and sex workers. He has been represented on South African and regional guidelines for over a decade, having done almost all his training within South Africa.

His major research currently focuses on low dose stavudine, low dose darunavir in second line treatment, and combinations of newer drugs to improve the resistance and potency of first line, as well as using patient information to drive improved linkage to care after diagnosis.



Ana Teasdale

Senior Adviser, Strategic Initiatives Pangaea Global AIDS

Ana Teasdale is the Senior Adviser for Strategic Initiatives. She provides senior counsel to Pangaea's CEO and Leadership Team on new and current programs and strategies, particularly in integrating HIV and broader women's health priorities, and liaison with the international community.

Ana has over 15 years of experience in international development. Prior to joining Pangaea Global AIDS, she worked at Susan G. Komen (Komen), Catholic Relief Services (CRS), and at the National Democratic Institute for International Affairs (NDI). At Komen Ana oversaw international community outreach grants and several global initiatives in Africa, Asia, Europe, and the Middle East to encourage breast cancer awareness, screening and treatment. She launched Komen's first public private partnership in The Bahamas (September 2012) and has played a key role in strengthening Komen's relationships with many NGO, government, and corporate partners worldwide. At CRS she supported CRS' field offices in the Middle East and Southeast Asia through strategic planning, training workshops, proposal development, and evaluations. At NDI she helped develop the Institute's political party, civil society, and parliamentary programs in Eastern Europe by collaborating with various political stakeholders.

A native of Serbia, Ana holds a B.A. in international relations and a M.A. in international communications with a focus on international development, both from the American University in Washington, D.C..



Lesley Scott

Professor Lesley Scott, PhD, Department of Molecular Medicine and Haematology at the University of the Witwatersrand, Johannesburg, South Africa.

Professor Scott is an applied scientist working in the field of laboratory medicine in the department of Molecular Medicine and Haematology at the University of the Witwatersrand, Johannesburg, South Africa. She is a WITS graduate, having performed all her qualifications at this University and for the past 18 years, focussed her research on designing, developing, evaluating and implementing laboratory quality diagnostics for HIV and TB, and 9 years prior in therapeutic vaccine research and manufacture. As head of research and development for the National Priority Program of the National Health Laboratory Service and Laboratory Director for the AIDS Clinical Trials Group, Lesley and her team have not only developed patented technologies that are used throughout the public sector laboratories in South Africa and other resource limited settings, but also developed a companion diagnostic that ensures quality testing in high throughput and point of care tuberculosis testing centres, that has also expanded into a global program now servicing 24 countries. Lesley continues to support the academic program at WITS University and supervisors scientists, registrars and biomedical engineering students in many exciting post graduate projects to improve patient care. Since 2004, Lesley has funded her group's activities through grants from NIH, Grand Challenges Canada, CDC, USAID/Peppar and others, and contributed to 78 publications and 130 conference abstracts resulting in several health care policy changes.



Francesca Celletti

M.D., Vice President of Innovation and New Technologies at the Elizabeth Glaser Pediatric AIDS Foundation.

Francesca Celletti is a global health expert, with a focus on innovations and health systems, global health initiatives and HIV/AIDS.

Francesca is the Vice President, Innovation and New Technology for EGPAF. Francesca manages the innovation to optimize program implementation efforts and expand access to HIV services, increase quality and improve the return on investments. Before this position, Francesca was Senior Director for Technical Leadership and Sustainability and she was overseeing the work of the Foundation on Technical Assistance and Health Systems. Before joining EGPAF, she was one of the Coordinator of the Health Systems cluster in WHO, where she led the WHO global effort to scale up the health workforce including mid-level cadres and community health workers in low and middle income countries. Before joining WHO, she worked in several countries in Africa and the Middle East to assist governments develop national health policies and implementing programmes aiming at strengthening health systems and optimize the service delivery system.

A native of Italy, Francesca received a MD degree from the University of Bologna, Italy. She trained in internal medicine and intensive care at University of Bologna and Harvard University, has a post-doctoral degree from Stanford University and PhD in International Relations from the Geneva School of Diplomacy.

