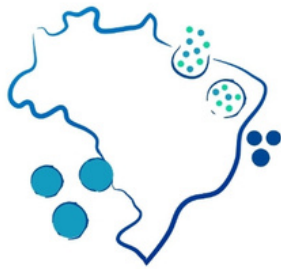


1ST Brazilian NanoScale Workshop: Exploring Extracellular Vesicles

APRIL, 09TH

08H - 09H	ACCREDITATION WITH OPENING COFFEE	
09H - 09H30	OPENING: WHAT TO EXPECT FROM THE 1ST BRAZILIAN WORKSHOP OF EXTRACELLULAR VESICLES?	DR ANDREA ALICE DA SILVA (UFF)
09H30 - 10H30	KEYNOTE: BIOFLUIDS AS A SOURCE OF EVS AND NON-EV COMPONENTS; CONSIDERATIONS FOR SAMPLE COLLECTION AND PROCESSING	DR JOHN TIGGES (BETH ISRAEL MEDICAL CENTER)
10H30 - 12H	SESSION: EXTRACELLULAR VESICLES AND PATHOGENS I	
10H30 - 11H	THE INFLUENCE OF EXTRACELLULAR VESICLES ON PARASITE-HOST INTERACTION AND THE IMMUNE SYSTEM	DR PATRÍCIA XANDER BATISTA (UNIFESP)
11H - 11H30	INVOLVEMENT OF EXTRACELLULAR VESICLES IN THE TROPISM AND PERSISTENCE OF TCRUZI DURING HOST INFECTION IN THE CHRONIC PHASE OF CHAGAS DISEASE	DR MARCEL RAMIREZ (FIOCRUZ-PR)
11H30 - 12H	EXTRACELLULAR MICROVESICLES CONTRIBUTE TO THE PATHOPHYSIOGENESIS OF HUMAN CUTANEOUS LEISHMANIASIS: A NANO-FLOW CYTOMETRIC APPROACH IN SAMPLES FROM PATIENTS BEFORE AND UNDER TREATMENT	DR ÁLVARO BERTHO (FIOCRUZ - RJ)
12H - 13H30	LUNCH BREAK	
13H30 - 15H	SESSION: EXTRACELLULAR VESICLES AND CANCER	
13H30 - 14H	THE PRO-TUMOR AND PRO-THROMBOTIC ROLE OF EXTRACELLULAR VESICLES	DR ROBSON DE QUEIROZ MONTEIRO (UFRJ)
14H - 14H30	IMPORTANCE OF TUMOR EXTRACELLULAR VESICLES IN NEUTROPHIL POLARIZATION: NEW THERAPEUTIC APPROACHES	DR JOÃO ALFREDO DE MORAES (UFRJ)
14H30 - 15H	EXTRACELLULAR VESICLES AS PROGNOSTIC BIOMARKERS AND CLINICAL MONITORING IN CANCER	DR ANDREA TEIXEIRA (UFMG)
15H - 16H	KEYNOTE: MISEV GUIDELINES UPDATES: INTERNATIONAL SOCIETY OF EXTRACELLULAR VESICLES (ISEV)	DR ANA CLÁUDIA TORRECILHAS (UNIFESP/ISEV)
16H - 17H	COFFEE BREAK & POSTER SESSION	
17H - 18H	KEYNOTE: THE THERAPEUTIC POTENTIAL OF EXTRACELLULAR VESICLES IN CANCER	DR PILAR MARTÍN (UNIVERSIDAD DE ZARAGOZA)

Organization:



1ST Brazilian NanoScale Workshop: Exploring Extracellular Vesicles

APRIL, 10TH

09H - 10H	KEYNOTE: URINARY EXTRACELLULAR VESICLES AS BIOMARKERS IN VASCULAR AND RENAL DISEASES	DR DYLAN BURGER (OTTAWA HOSPITAL RESEARCH INSTITUTE)
10H - 11H	SESSION: EXTRACELLULAR VESICLES AND APPLICATIONS IN THE CLINIC	
10H - 10H30	MESENCHYMAL EXTRACELLULAR CELL AND VESICLE THERAPY FOR CHRONIC LUNG DISEASES: EXPERIMENTAL AND CLINICAL EVIDENCE	DR PATRICIA RIEKEN MACEDO ROCCO (UFRJ)
10H30 - 11H	PROFILE OF CIRCULATING MICROVESICLES ASSOCIATED WITH CLINICAL-LABORATORY PARAMETERS AS A PREDICTIVE TOOL FOR THE SEVERITY OF COVID-19	DR ANDRÉA TEIXEIRA DE CARVALHO (FIOCRUZ-MG)
11H - 12H	SESSION: EXTRACELLULAR VESICLES, INFLAMMATION AND REPAIR	
11H - 11H30	EXTRACELLULAR MATRIX VESICLES AND THEIR IMPORTANT ROLE IN THE BIOMINERALIZATION PROCESS	DR PIETRO CIANCAGLINI (USP)
11H30 - 12H	EVS IN INFLAMMATION AND THROMBOSIS	DR PATRÍCIA BOZZA (FIOCRUZ)
12H - 13H30	LUNCH BREAK	
13H30 - 16H	SESSION: EXTRACELLULAR VESICLES AND PATHOGENS II	
13H30 - 14H	EXTRACELLULAR VESICLES AS A FUNGAL VIRULENCE FACTOR	DR FAUSTO DOS REIS ALMEIDA (FMRP/USP)
14H - 14H30	IMMUNE-STIMULATORY EFFECT OF MEMBRANE VESICLES RELEASED BY MYCOBACTERIUM SPP-INFECTED MACROPHAGES	DR TONYA AZEVEDO DUARTE (UFBA)
14H30 - 15H	EXTRACELLULAR VESICLES FROM NEUROPEPTIDE-ACTIVATED HUMAN MACROPHAGES MEDIATE ANTI-SARS-COV-2 PROTECTION IN MONOCYTES.	DR DUMITH CHEQUER BOU HABIB (FIOCRUZ-RJ)
15H - 15H30	INVOLVEMENT OF EXTRACELLULAR VESICLES FROM SPOROTHRIX BRASILIENSIS YEAST IN THE IMMUNE RESPONSE AGAINST SPOROTRICHOSIS	DR KAREN SPADARI FERREIRA (UNIFESP)
15H30 - 16H	MALARIA'S COMMUNICATORS: EXTRACELLULAR VESICLES IN SURVIVAL STRATEGIES	DR NETA REGEV RUDZKI (WEIZMANN INSTITUTE OF SCIENCE, ISRAEL)
16H - 17H	COFFEE BREAK & POSTER SESSION	
17H - 18H	PRESENTING THE NEW CYTOFLEX NANO	BECKMAN COULTER LIFE SCIENCES
18H - 18H30	WRAP-UP & AWARDS	COMISSÃO ORGANIZADORA

Organization:

