

Title	Session	Author(s)	Date	Session Time	Location
GENOME-WIDE ASSOCIATION STUDIES OF BLOOD-BASED BIOMARKERS OF ALZHEIMER'S DISEASE IMPLICATE APOE AND TMEM106B	2730 - ANCESTRY AND GENETICS OF NEURODEGENERATION (I D 15)	Michael J. Rigby (United States of America)	Wed, 06.03.2024	08:40 - 10:40	Auditorium V
DEVELOPMENT OF A PROTOTYPE POINT OF CARE TEST WITH FOPPR TECHNOLOGY FOR ULTRASENSITIVE DETECTION OF NEUROFILAMENT LIGHT CHAIN IN PLASMA	2890 - FLUID BIOMARKERS, AND IMAGING AND DIGITAL TOOLS (ID 31)	Madison I. Honey (Netherlands)	Wed, 06.03.2024	17:00 - 19:00	Auditorium VI+VII
COMPARISON OF TWO PLASMA P-TAU217 ASSAYS TO DETECT AND MONITOR ALZHEIMER'S PATHOLOGY	2890 - FLUID BIOMARKERS, AND IMAGING AND DIGITAL TOOLS (ID 31)	Joseph Therriault (Canada)	Wed, 06.03.2024	17:00 - 19:00	Auditorium VI+VII
BRAINS IN FOCUS: UNCOVERING TAU AGGREGATE CHANGES ACROSS ALZHEIMER'S DISEASE WITH SINGLE-MOLECULE IMAGING	3630 - TAU MECHANISMS 02 (ID 43)	Elizabeth A. English (United Kingdom)	Thu, 07.03.2024	08:40 - 10:40	Auditorium III + IV
AGREEMENTS BETWEEN COGNITIVE STATUS AND PLASMA BIOMARKERS POSITIVITY IN PRECLINICAL AND PRODROMAL ALZHEIMER'S DISEASE	3710 - FLUID BIOMARKERS (ID 51)	Bjørn-Eivind Kirsebom (Norway)	Thu, 07.03.2024	13:50 - 15:50	Auditorium II
RENAL FUNCTION AND PLASMA ATN BIOMARKERS OF ALZHEIMER'S DISEASE AMONG DIVERSE MIDDLE-AGED AND OLDER HISPANICS/LATINOS: SOL-INCA (HCHS/SOL) RESULTS.	3710 - FLUID BIOMARKERS (ID 51)	Hector Gonzalez (United States of America)	Thu, 07.03.2024	13:50 - 15:50	Auditorium II
TARGETING MULTIPHOSPHORYLATED TAU: PLASMA TAU SIMULTANEOUSLY PHOSPHORYLATED AT T217 AND T231 SURPASSES PERFORMANCE OF P-TAU217	3710 - FLUID BIOMARKERS (ID 51)	Anna Lidia Woidala (Italy)	Thu, 07.03.2024	13:50 - 15:50	Auditorium II
THE ASSOCIATION BETWEEN PLASMA BIOMARKERS, ROUTINE CSF AND BLOOD BIOMARKERS, AND COMORBIDITIES IN A MIXED MEMORY CLINIC COHORT	3710 - FLUID BIOMARKERS (ID 51)	Helena Sophia Gleerup (Denmark)	Thu, 07.03.2024	13:50 - 15:50	Auditorium II
THE EFFECT OF APOE4 CARRIER STATUS ON ALZHEIMER'S DISEASE PLASMA BIOMARKER CONCENTRATIONS IN HEALTHY SENIORS	3710 - FLUID BIOMARKERS (ID 51)	Jennifer G. Cooper (Canada)	Thu, 07.03.2024	13:50 - 15:50	Auditorium II
PLASMA PTAU-217 SUPERIORITY COMPARED TO OTHER PLASMA BIOMARKERS IN PREDICTION OF AMYLOID PET POSITIVITY IN TERTIARY MEMORY CLINIC PATIENTS	6190 - FLUID BIOMARKERS AND IMAGING 01 (ID 73)	Marco Buccì (Sweden)	Fri, 08.03.2024	08:40 - 10:40	Auditorium VIII
RACIAL DISCRIMINATION DURING MIDDLE AGE PREDICTS HIGHER SERUM PHOSPHORYLATED TAU AND NEUROFILAMENT LIGHT CHAIN LEVELS A DECADE LATER	6190 - FLUID BIOMARKERS AND IMAGING 01 (ID 73)	Michelle Mielke (United States of America)	Fri, 08.03.2024	08:40 - 10:40	Auditorium VIII
PLASMA P-TAU212 IS ASSOCIATED WITH COGNITIVE DECLINE AND DISEASE PROGRESSION IN COGNITIVELY UNIMPAIRED PEOPLE.	6250 - FLUID BIOMARKERS AND IMAGING 02 (ID 79)	Przemysław R. Kac (Sweden)	Fri, 08.03.2024	13:50 - 15:50	Auditorium I
RELATIONSHIP OF PLASMA NFL, GFAP, AB1-42/AB1-40 AND PTAU181 WITH SYNAPTIC DENSITY AND ALZHEIMER'S DISEASE HALLMARKS IN NON-DEMENTED OLDER ADULTS	6250 - FLUID BIOMARKERS AND IMAGING 02 (ID 79)	Steffi De Meyer (Belgium)	Fri, 08.03.2024	13:50 - 15:50	Auditorium I
EVALUATING SEX-RELATED DISPARITIES IN THE PROGNOSTIC PROPERTIES OF PLASMA BIOMARKERS FOR ALZHEIMER'S DISEASE	6270 - SEX SIMILARITIES AND DIFFERENCES IN BIOMARKERS & PATHOPHYSIOLOGY (ID 81)	Marta Milà Alomà (United States of America)	Fri, 08.03.2024	13:50 - 15:50	Auditorium II

PLASMA P-TAU181, NFL AND GFAP ARE ASSOCIATED WITH COGNITIVE, FUNCTIONAL, AND MRI MARKERS IN A LONGITUDINAL COHORT OF NEURODEGENERATIVE DISEASES	6350 - MRI (ID 90)	Erlan Sanchez (Canada)	Fri, 08.03.2024	16:20 - 18:30	Auditorium V
IMPROVING DIFFERENTIAL DIAGNOSIS OF AD AND OTHER TAUOPATHIES IN BLOOD WITH A NEW BIOMARKER OF INFLAMMASOME FORMATION: ASC SPECKS	6360 - FLUID BIOMARKERS AND IMAGING 03 (ID 91)	Evgeniia Lobanova (United Kingdom)	Fri, 08.03.2024	16:20 - 18:30	Auditorium VIII
PLASMA BIOMARKERS IN FORMER PROFESSIONAL BOXERS: A 7-YEAR LONGITUDINAL STUDY	6360 - FLUID BIOMARKERS AND IMAGING 03 (ID 91)	Suzie Kamps (Netherlands)	Fri, 08.03.2024	16:20 - 18:30	Auditorium VIII
SIMOA CSF P-TAU205 ASSAY TRACKS IN VIVO TAU PATHOLOGY DEPOSITION, NEURODEGENERATION AND COGNITIVE DECLINE IN ALZHEIMER'S DISEASE PATIENTS.	6360 - FLUID BIOMARKERS AND IMAGING 03 (ID 91)	Juan Lantero-Rodriguez (Sweden)	Fri, 08.03.2024	16:20 - 18:30	Auditorium VIII
DEVELOPMENT OF POTENT, ORALLY BIOAVAILABLE LRRK2 PROTAC® DEGRADER MOLECULES AS POTENTIAL DISEASE MODIFYING THERAPEUTICS FOR PARKINSON'S DISEASE	6361 - LRRK2, ALPHA – SYNUCLEIN, DIAGNOSIS AND THERAPEUTIC TARGETS (ID 74)	Jere Meredith (United States of America)	Fri, 08.03.2024	16:20 - 18:30	Auditorium III + IV
USING BLOOD BIOMARKERS TO IMPROVE DIAGNOSIS OF BRAIN INJURY AND INFORM ON BRAIN AGING AND DEMENTIA RISK IN SURVIVORS OF INTIMATE PARTNER VIOLENCE	6720 - BIOMARKERS IN CLINICAL USE ACROSS DEMENTIAS (ID 103)	Cheryl Wellington (Canada)	Sat, 09.03.2024	08:40 - 10:40	Auditorium VI
104-WEEK EFFECTS OF ALZ-801/VALILTRAMIPROSATE ON PLASMA, MRI BIOMARKERS AND COGNITION SUPPORT DISEASE MODIFICATION IN APOE4 CARRIERS WITH EARLY ALZHEIMER'S DISEASE	6730 - TRANSLATIONAL TREATMENT STRATEGIES AND NEW TARGETS (ID 105)	John Hey (United States of America)	Sat, 09.03.2024	08:40 - 10:40	Auditorium V
DEFICIENCY OF THE SORL1 AND VPS26B-RETROMER COMPONENT PHENOCOPY EACH OTHER AT THE ANATOMICAL, BEHAVIORAL, AND BIOMARKER LEVELS.	6780 - SORL1 (ID 429)	Sabrina Simoes (United States of America)	Sat, 09.03.2024	11:10 - 13:10	Auditorium III + IV
CEREBROSPINAL FLUID PHOSPHO-TAU181, 217, AND 231 ARE INCREASED IN AUTOPSY-VERIFIED CREUTZFELDT-JAKOB DISEASE AND ASSOCIATE WITH NEURODEGENERATION AND AMYLOID NEUROPATHOLOGY	6800 - BIOMARKERS AND MULTIMODAL IMAGING IN AD, PD, STROKE AND CREUTZFELDT-JAKOB DISEASE (ID 111)	Andreja Emersic (Slovenia)	Sat, 09.03.2024	11:10 - 13:10	Auditorium VI
ASSOCIATION OF EGFR WITH PLASMA ALZHEIMER'S DISEASE BIOMARKERS IN AMYLOID-NEGATIVE INDIVIDUALS – A PILOT STUDY	6900 - FLUID BIOMARKERS AND IMAGING 04 (ID 122)	Burak Arslan (Sweden)	Sat, 09.03.2024	14:10 - 16:10	Auditorium III + IV
DIAGNOSTIC ACCURACY OF TWO PLASMA PTAU217 IMMUNOASSAYS AND A MASS SPECTOMETRY PTAU217 OCCUPANCY RATIO ASSAY FOR PREDICTION OF AMYLOID PATHOLOGY.	6900 - FLUID BIOMARKERS AND IMAGING 04 (ID 122)	Alicia Algeciras-Schimmich (United States of America)	Sat, 09.03.2024	14:10 - 16:10	Auditorium III + IV
POOR SLEEP EFFICIENCY IS LINKED TO ELEVATED BLOOD-BASED BIOMARKERS OF ALZHEIMER'S DISEASE AND NEURODEGENERATION IN OLDER ADULTS LIVING WITH SUBJECTIVE COGNITIVE DECLINE	6950 - EEG, SLEEP, AND VIRTUAL ASSESSMENT TOOLS (ID 126)	Johannes C. Michaelian (Australia)	Sat, 09.03.2024	16:40 - 18:40	Auditorium V
EFFECTS OF TAU DELETION AND TAU INDUCTION ON NEURODEGENERATION IN A MOUSE MODEL OF CEREBRAL BETA-AMYLOIDOSIS	6970 - NEURODEGENERATION: DIVERSE MECHANISMS (ID 128)	Carina Bergmann (Germany)	Sat, 09.03.2024	16:40 - 18:40	Auditorium II
ALZPATH PTAU217 PERFORMANCE IDENTIFYING ALZHEIMER'S PATHOLOGY IN THE BRAIN AND BODY DONATION PROGRAM (BBDP) AT BANNER SUN HEALTH RESEARCH INSTITUTE	6980 - NEURODEGENERATION: DIVERSE MECHANISMS FLUID BIOMARKERS AND IMAGING (ID 431)	Lauren Chaby (United States of America)	Sat, 09.03.2024	16:40 - 18:40	Auditorium III + IV
EFFECT OF KIDNEY FUNCTION ON PLASMA PTAU217 CONCENTRATIONS USING DIFFERENT PTAU217 ASSAYS.	6980 - NEURODEGENERATION: DIVERSE MECHANISMS	Joshua A. Bornhorst (United States of America)	Sat, 09.03.2024	16:40 - 18:40	Auditorium III + IV

	<u>FLUID BIOMARKERS AND IMAGING (ID 431)</u>				
PLASMA P-TAU217 DEMONSTRATES INCREASED ASSOCIATION WITH REGIONAL AMYLOID AND TAU PET LOAD COMPARED TO AB42/40 RATIO OR PTAU181 IN AD	VIRTUAL ORAL (VO039 / #2970) : THEME A: B-AMYLOID DISEASES / A04.C. IMAGING, BIOMARKERS, DIAGNOSTICS: PET - AMYLOID	<u>Han Yin1</u> , <u>Pallavi Sachdev2</u> , <u>Arnaud Charil3</u> , <u>David Verbel1</u> , <u>Michael Irizarry4</u> , <u>Larisa Reyderman3</u> , <u>Shobha Dhadda5</u> , <u>Joel Braunstein6</u> , <u>Philip Vergheze6</u> , <u>Tim West6</u>			
CEREBROSPINAL FLUID PHOSPHO-TAU181, 217, AND 231 ARE INCREASED IN AUTOPSYVERIFIED CREUTZFELDT-JAKOB DISEASE AND ASSOCIATE WITH NEURODEGENERATION AND AMYLOID NEUROPATHOLOGY	VIRTUAL ORAL (VO130 / #2946) : THEME F: PRION DISEASES / F03. IMAGING, BIOMARKERS, DIAGNOSTICS	<u>Andreja Emersic1</u> , <u>Nicholas Ashton2,3,4,5</u> , <u>Agathe Vrillon6,7</u> , <u>Juan Lantero-Rodriguez4</u> , <u>Mlakar Jernei8</u> , <u>Milica Gregoric Kramberger1,9</u> , <u>Fernando Gonzalez-Ortiz4</u> , <u>Przemyslaw Kac4</u> , <u>Maciej Dulewicz4</u> , <u>Jörg Hanrieder4,10</u> , <u>Eugeen Vanmechelen11</u> , <u>Cucnik Saša1,12,13</u> , <u>Uroš Rot1,9</u> , <u>Henrik Zetterberg4,14,15,16,17,18</u> , <u>Thomas Karikari4,19</u> , <u>Kaj Blennow4,14</u>			