

| Focus | Article | Authors | Journal |
|------------------------|---|--------------------------|--------------------------------------|
| 2024 | | | |
| Neurology | Comparison of plasma soluble and extracellular vesicles-associated biomarkers in Alzheimer's Disease patients and cognitively normal | E Boyer et al. | Alzheimers Res Ther. 2024 Jun |
| Neurology | Plasma extracellular vesicle tau and TDP-43 as diagnostic biomarkers in FTD and ALS | M Chatterjee et al. | Nat Med. 2024 Jun |
| Extracellular Vesicles | Before Translating Extracellular Vesicles into Personalized Diagnostics and Therapeutics: What We Could Do | CA Cheng | Mol Pharm. 2024 Jun |
| Extracellular Vesicles | Addressing Heterogeneity in Direct Analysis of Extracellular Vesicles and Their Analogs by Membrane Sensing Peptides as Pan-Vesicular Affinity | A Gori et al. | Adv. Sci. 2024 May |
| Extracellular Vesicles | 2023 White Paper on Recent Issues in Bioanalysis (PART 2) | O Kholmanskikh et al. | Bioanalysis. 2024 May |
| Neurology | Single-extracellular vesicle (EV) analyses validate the use of L1 Cell Adhesion Molecule (L1CAM) as a reliable biomarker of neuron-derived EVs | C Noguerras-Ortiz et al. | J Extracell Vesicles. 2024 May |
| Oncology | Measuring EGFr on extracellular vesicles in human plasma samples | DA Olsen et al. | AACR poster, 2024 Apr |
| Neurology | Specific binding of Alzheimer's A β peptides to extracellular vesicles | C Coughlan et al. | Int. J. Mol. Sci. 2024 Mar |
| Neurology | Endo-lysosomal dysfunction in neurodegenerative diseases: opinion on current progress and future direction in the use of exosomes as biomarkers | M Herman et al. | Phil. Trans. R. Soc. 2024 Feb |
| Extracellular Vesicles | Exosome subpopulations: the isolation and the functions in diseases | X Ma et al. | Gene. 2024 Jan |
| Oncology | Ultrasensitive and high-resolution protein spatially decoding framework for tumor extracellular vesicles | CA Cheng et al. | Adv. Sci. 2024 Jan |
| 2023 | | | |
| Neurology | Proteomic comparison between non-purified cerebrospinal fluid and cerebrospinal fluid-derived extracellular vesicles from patients with Alzheimer's, Parkinson's and Lewy body dementia | Y Hirschberg et al. | J Extracell Vesicles. 2023 Dec |
| Neurology | ATP1A3 as a target for isolating neuron-specific extracellular vesicles from human brain and biofluids | Y You et al. | Sci. Adv. 2023 Sept |
| Extracellular Vesicles | Extracellular vesicles: techniques and biomedical applications related to single vesicle analysis | J Zhang et al. | American Chemical Society. 2023 Sept |

| Focus | Article | Authors | Journal |
|------------------------|---|------------------------|--|
| Neurology | Assessment of risk factors for postoperative delirium in older adults who underwent spinal surgery and identifying associated biomarkers using exosomal protein | W Baek et al. | J Korean Acad Nurs. 2023 Aug |
| Extracellular Vesicles | Proof of concept of using a membrane-sensing peptide for sEVs affinity-based isolation | B Benayas et al. | Front. Bioeng. Biotechnol. 2023 Aug |
| Technology | Ultrasensitive protein detection technologies for extracellular vesicle measurements | A Shami-Shah et al. | Mol Cell Proteomics. 2023 Jun |
| Technology | Improved isolation of extracellular vesicles by removal of both free proteins and lipoproteins | D Ter-Ovanesyan et al. | eLife. 2023 May |
| Neurology | An update on peripheral blood extracellular vesicles as biomarkers for Parkinson's disease diagnosis | K Wang et al. | Neuroscience. 2023 Feb |
| Cardiology | Rapid activation of neuroinflammation in stroke: plasma and extracellular vesicles obtained on a mobile stroke unit | RG Kowalski et al. | Stroke. 2023 Feb |
| 2022 | | | |
| Technology | Comparing digital detection platforms in high sensitivity immune-phenotyping of extracellular vesicles | R Frigerio et al. | Journal of Extracellular Biology. 2022 Aug |
| Technology | Future of digital assays to resolve clinical heterogeneity of single extracellular vesicles | RT Morales et al. | ACS nano. 2022 Jul |
| Neurology | Pathological (dis)similarities in neuronal exosome-derived synaptic and organellar marker levels between Alzheimer's disease and frontotemporal dementia | G Krishna et al. | Journal of Alzheimer's disease. 2022 Jul |
| Neurology | β -Amyloid in blood neuronal-derived extracellular vesicles is elevated in cognitively normal adults at risk of Alzheimer's disease and predicts cerebral amyloidosis | TR Li et al. | Alzheimers Res Ther. 2022 May |
| Neurology | KIBRA regulates amyloid β metabolism by controlling extracellular vesicles secretion | X Han et al. | EBioMedicine. 2022 Mar |

| Focus | Article | Authors | Journal |
|-------------|---|------------------------|------------------------------------|
| Neurology | Extracellular vesicle neurofilament light is elevated within the first 12-months following traumatic brain injury in a U.S military population | VA Guedes et al. | Sci Rep. 2022 Mar |
| Oncology | Fast quantification of extracellular vesicles levels in early breast cancer patients by single molecule detection array (SIMOA) | C Morasso et al. | Breast Cancer Res Treat. 2022 Feb |
| Neurology | Poor sleep quality is linked to elevated extracellular vesicle-associated inflammatory cytokines in warfighters with chronic mild traumatic brain injuries | JL Gottshall et al. | Front Pharmacol. 2022 Jan |
| 2021 | | | |
| Technology | Framework for rapid comparison of extracellular vesicle isolation methods | D Ter-Ovanesyan et al. | eLife. 2021 Nov |
| Technology | Advances in the analysis of single extracellular vesicles: a critical review | SH Hilton et al. | Sensors and Actuators Reports. |
| Oncology | Universal extracellular vesicles and PD-L1+ extracellular vesicles detected by single molecule array technology as circulating biomarkers for diffuse large B cell lymphoma | JW Li et al. | Oncoimmunology. 2021 Oct |
| Neurology | Neuronally-derived tau is increased in experienced breachers and is associated with neurobehavioral symptoms | KA Edwards et al. | Sci Rep. 2021 Sept |
| Neurology | Spinal fluid myeloid microvesicles predict disease course in multiple sclerosis | S Gelibter et al. | Ann Neurol. 2021 Aug |
| Neurology | Extracellular vesicles as distinct biomarker reservoirs for mild traumatic brain injury diagnosis | K Beard et al. | Brain Commun. 2021 Jul |
| Oncology | PD-L1 detection on circulating tumor-derived extracellular vesicles (T-EVs) from patients with lung cancer | F Wu et al. | Transl Lung Cancer Res. 2021 Jun |
| Technology | Microfluidics-based technologies for the analysis of extracellular vesicles at the single-cell level and single-vesicle level | F Zhu et al. | Chinese Chemical Letters. 2021 Jun |
| Neurology | L1CAM is not associated with extracellular vesicles in human cerebrospinal fluid or plasma | M Norman et al. | Nat Methods. 2021 Jun |
| Neurology | Synaptic and complement markers in extracellular vesicles in multiple sclerosis | P Bhargava et al. | Mult Scler. 2021 Apr |

| Focus | Article | Authors | Journal |
|-------------|--|----------------------|--------------------------------|
| Oncology | Enrichment of circulating tumor-derived extracellular vesicles from human plasma | KE Yoh et al. | J Immunol Methods. 2021 Mar |
| Neurology | Extracellular vesicle concentrations of glial fibrillary acidic protein and neurofilament light measured 1 year after traumatic brain injury | S Flynn et al. | Sci Rep. 2021 Feb |
| 2020 | | | |
| Oncology | Clinical significance of circulating exosomal PD-L1 and soluble PD-L1 in extranodal NK/T-cell lymphoma, nasal-type | JW Li et al. | Am J Cancer Res. 2020 Dec |
| Neurology | Functional and structural neuroimaging correlates of repetitive low-level blast exposure in career breachers | JR Stone et al. | J Neurotrauma. 2020 Dec |
| Neurology | Blood biomarkers of traumatic brain injury and cognitive impairment in older veterans | CB Peltz et al. | Neurology. 2020 Sep |
| Oncology | Plasma extracellular vesicles detected by single molecule array technology as a liquid biopsy for colorectal cancer | P Wei et al. | J Extracell Vesicles. 2020 Aug |
| Neurology | Proteomic profiling of extracellular vesicles derived from cerebrospinal fluid of Alzheimer's disease patients: a pilot study | S Muraoka et al. | Cells. 2020 Aug |
| Neurology | Long noncoding RNA POU3F3 and alpha-synuclein in plasma L1CAM exosomes combined with beta-glucocerebrosidase activity: potential predictors of Parkinson's disease | J Zou et al. | Neurotherapeutics. 2020 Jul |
| Neurology | Exosomal neurofilament light: a prognostic biomarker for remote symptoms after mild traumatic brain injury? | VA Guedes et al. | Neurology. 2020 Jun |
| Neurology | Circulating brain injury exosomal proteins following moderate-to-severe traumatic brain injury: temporal profile, outcome prediction and therapy implications | S Mondello et al. | Cells. 2020 Apr |
| 2019 | | | |
| Neurology | Association of extracellular vesicle biomarkers with Alzheimer's disease in the Baltimore longitudinal study of aging | D Kapogiannis et al. | JAMA Neurol. 2019 Nov |

| Focus | Article | Authors | Journal |
|-------------|---|--------------------|-----------------------------|
| Neurology | Assessing neuronal and astrocyte derived exosomes from individuals with mild traumatic brain injury for markers of neurodegeneration and cytotoxic activity | CN Winston et al. | Front Neurosci. 2019 Oct |
| Neurology | New windows into the brain: central nervous system-derived extracellular vesicles in blood | M Shi et al. | Prog Neurobiol. 2019 Apr |
| Neurology | Novel tau fragments in cerebrospinal fluid: relation to tangle pathology and cognitive decline in Alzheimer's disease | C Cicognola et al. | Acta Neuropathol. 2019 Feb |
| 2018 | | | |
| Neurology | Diagnosis of traumatic brain injury using miRNA signatures in nanomagnetically isolated brain-derived extracellular vesicles | J Ko et al. | Lab Chip. 2018 Dec |
| Neurology | Higher exosomal tau, amyloid-beta 42 and IL-10 are associated with mild TBIs and chronic symptoms in military personnel | J Gill et al. | Brain Inj. 2018 Jun |
| Neurology | Higher exosomal phosphorylated tau and total tau among veterans with combat-related repetitive chronic mild traumatic brain injury | K Kenney et al. | Brain Inj. 2018 Jun |
| Neurology | Detection of aggregation-competent tau in neuron-derived extracellular vesicles | FX Guix et al. | Int J Mol Sci. 2018 Feb |
| 2016 | | | |
| Neurology | CNS tau efflux via exosomes is likely increased in Parkinson's disease but not in Alzheimer's disease | M Shi et al. | Alzheimers Dement. 2016 Nov |