

Focus	Article	Authors	Journal
2024			
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			
Extracellular Vesicles	Addressing heterogeneity in direct analysis of extracellular vesicles and analogues using membrane-sensing peptides as pan-affinity probes	A Gori et al.	bioRxiv. 2023 Dec
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
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
Neurology	Plasma extracellular vesicle tau isoform ratios and TDP-43 inform about molecular pathology in frontotemporal dementia and amyotrophic lateral sclerosis	A Schneider et al.	Preprint, Research Square. 2023 Jul
Technology	Ultrasensitive protein detection technologies for extracellular vesicle measurements	A Shami-Shah et al.	Mol Cell Proteomics. 2023 Jun

Focus	Article	Authors	Journal
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	Morales RT and Ko J	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
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

Focus	Article	Authors	Journal
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 July
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 June
Neurology	Synaptic and complement markers in extracellular vesicles in multiple sclerosis	P Bhargava et al.	Mult Scler. 2021 Apr
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

Focus	Article	Authors	Journal
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
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

Focus	Article	Authors	Journal
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