



Highlighted Celselect Publications

Cheng YH, Chen YC, Lin E, Brien R, Jung S, Chen YT, Lee W, Hao Z, Sahoo S, Min Kang H, Cong J, Burness M, Nagrath S, S Wicha M, Yoon E. Hydro-Seq enables contamination-free high-throughput single-cell RNA-sequencing for circulating tumor cells. *Nat Commun.* 2019 May 15;10(1):2163. doi: 10.1038/s41467-019-10122-2. PMID: 31092822; PMCID: PMC6520360.

Gogoi P, Sepehri S, Zhou Y, Gorin MA, Paolillo C, Capoluongo E, Gleason K, Payne A, Boniface B, Cristofanilli M, Morgan TM, Fortina P, Pienta KJ, Handique K, Wang Y. Development of an Automated and Sensitive Microfluidic Device for Capturing and Characterizing Circulating Tumor Cells (CTCs) from Clinical Blood Samples. *PLoS One.* 2016 Jan 25;11(1):e0147400. doi: 10.1371/journal.pone.0147400. PMID: 26808060; PMCID: PMC4726586.

Horimoto Y, Tokuda E, Murakami F, Uomori T, Himuro T, Nakai K, Orihata G, Iijima K, Togo S, Shimizu H, Saito M. Analysis of circulating tumour cell and the epithelial mesenchymal transition (EMT) status during eribulin-based treatment in 22 patients with metastatic breast cancer: a pilot study. *J Transl Med.* 2018 Oct 20;16(1):287. doi: 10.1186/s12967-018-1663-8. PMID: 30342534; PMCID: PMC6195982.

Konno N, Suzuki R, Takagi T, Sugimoto M, Asama H, Sato Y, Irie H, Hikichi T, Ohira H. Clinical utility of a newly developed microfluidic device for detecting circulating tumor cells in the blood of patients with pancreatico-biliary malignancies. *J Hepatobiliary Pancreat Sci.* 2021 Jan;28(1):115-124. doi: 10.1002/jhbp.850. Epub 2020 Nov 6. PMID: 33090657.

Kou R, Zhao J, Gogoi P, Carskadon S, Chow W, Hwang C, Palanisamy N, Leung C, Wang Y. Enrichment and mutation detection of circulating tumor cells from blood samples. *Oncol Rep.* 2018 Jun;39(6):2537-2544. doi: 10.3892/or.2018.6342. Epub 2018 Mar 30. PMID: 29620284; PMCID: PMC5983925.

Nakasato T, Kusaka C, Ota M, Hasebe Y, Ueda K, Unoki T, Oshinomi K, Morita J, Maeda Y, Shichijo T, Naoe M, Ogawa Y. A Novel, Circulating Tumor Cell Enrichment Method Reduces ARv7 False Positivity in Patients with Castration-Resistant Prostate Cancer. *Diagnostics (Basel).* 2020 Mar 11;10(3):151. doi: 10.3390/diagnostics10030151. PMID: 32168745; PMCID: PMC7151149.

Riahi R, Gogoi P, Sepehri S, Zhou Y, Handique K, Godsey J, Wang Y. A novel microchannel-based device to capture and analyze circulating tumor cells (CTCs) of breast cancer. *Int J Oncol.* 2014 Jun;44(6):1870-8. doi: 10.3892/ijo.2014.2353. Epub 2014 Mar 21. PMID: 24676558; PMCID: PMC4063542.

Ujiiie D, Matsumoto T, Endo E, Okayama H, Fujita S, Kanke Y, Watanabe Y, Hanayama H, Hayase S, Saze Z, Ohki S, Kono K. Circulating tumor cells after neoadjuvant chemotherapy are related with recurrence in esophageal squamous cell carcinoma. *Esophagus.* 2021 Jul;18(3):566-573. doi: 10.1007/s10388-021-00829-x. Epub 2021 Mar 4. PMID: 33661456.

Wang X, Jaimes M, Gu H, Shults K, Putta S, Sharma V, Chow W, Gogoi P, Handique K, Patterson BK. Cell by cell immuno- and cancer marker profiling of non-small cell lung cancer tissue: Checkpoint marker expression on CD103+, CD4+ T-cells predicts circulating tumor cells. *Transl Oncol.* 2021 Jan;14(1):100953. doi: 10.1016/j.tranon.2020.100953. Epub 2020 Nov 18. PMID: 33217647; PMCID: PMC7683336.



**Bio-Rad
Laboratories, Inc.**