

# The Role of Multiplex Immunoassays in Early Cancer Detection

## Meeting the Demand for Sensitivity, Reliability, and Efficiency

Nearly 30% of bladder cancer patients are diagnosed at stage 2 or higher, where five-year survival falls to around 50%.<sup>1</sup> The reason? Current diagnostics cannot reliably detect the disease before clinical presentation. This gap represents more than a technical challenge—it reflects a fundamental limitation of single-biomarker approaches in a disease defined by complexity.

Cancer progression and therapeutic response rarely depend on a single molecular signal. Instead, they reflect the interplay of cytokines, growth factors, and signaling networks within the tumor microenvironment. To capture this complexity, researchers are increasingly turning to multiplex immunoassays—platforms capable of measuring dozens of analytes from small-volume samples.

By conserving patient specimens, increasing throughput, and revealing systems-level insights, multiplexing provides a more comprehensive view of tumor biology than single-analyte approaches can offer. But realizing this potential demands assays that deliver more than high throughput. They must achieve

ultra-high sensitivity, reproducibility, and low cross-reactivity, while performing reliably in complex matrices such as serum or plasma.

Meeting these standards requires antibody expertise, rigorous validation, and panel designs flexible enough to address diverse research needs. **R&D Systems Luminex® Assays Instruments** combine broad multiplexing capacity with validated antibodies, optimized diluents, and one of the industry's largest mix-and-match analyte menus, providing the sensitivity, reliability, and flexibility essential for advancing cancer detection and therapeutic development.



## Connecting the Dots in Cancer Biology

Cancer biology cannot be reduced to a single molecular signal. Cytokines such as IL-6, TNF- $\alpha$ , and IL-10 play overlapping yet distinct roles in the tumor microenvironment. Measuring one analyte provides only a fragment of the picture, whereas multiplexing offers a panoramic view of immune–tumor interactions. This broader perspective enables the identification of biomarker patterns that correlate with disease stage, predict therapeutic response, or signal recurrence.

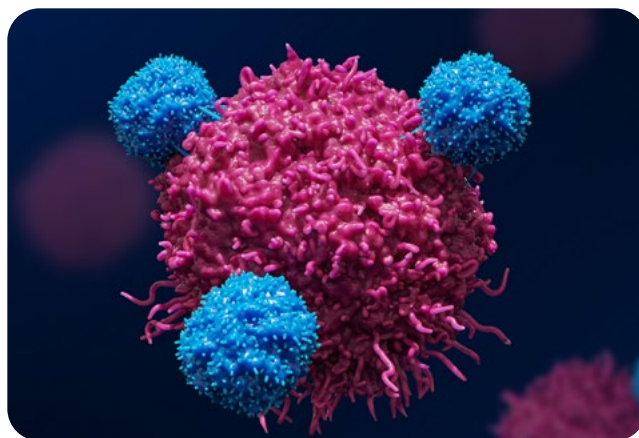
Traditional single-analyte assays fall short in this setting. They may overlook subtle but clinically important changes in low-abundance cytokines or fail to capture the interplay between opposing immune factors. For instance, a rise in a pro-inflammatory cytokine could be counterbalanced by an anti-inflammatory mediator—a dynamic invisible when measured in isolation.

Multiplex immunoassays overcome these limitations by simultaneously quantifying dozens of cytokines and soluble factors from a single plasma or serum sample. This dual capture of tumor-derived signals and systemic immune activation has already revealed cytokine signatures linked to prognosis and treatment response. In studies of gastrointestinal cancers, patients with high baseline levels of two cytokine patterns—one associated with checkpoint activity and another with T-cell trafficking—showed better prognosis and higher overall response rates to checkpoint blockade. Critically, multiplex profiling enabled these signatures to be identified before clinical outcomes were evident.<sup>2</sup>

Longitudinal multiplex analysis extends this capability further, tracking immune shifts during therapy and distinguishing emerging responders from those developing resistance or toxicity. Such insights not only refine patient selection but also give clinicians the opportunity to adapt treatment strategies earlier and more effectively.

## Accuracy Starts with the Right Antibodies

Detecting low-abundance biomarkers—such as early-stage tumor cytokines—requires assays with both exceptional sensitivity and specificity. In multiplex settings, this challenge is even greater: the assay must remain sensitive enough to detect subtle changes while avoiding cross-reactivity among closely related proteins.



The key lies in antibody quality. High-performance systems rely on rigorously selected antibody pairs with proven specificity. Extensive validation against potential cross-reactivity ensures that even large, complex panels produce accurate, reproducible results. This attention to antibody design allows multiplex assays to detect analytes at picogram-per-milliliter concentrations—levels that can serve as early warning signs before clinical symptoms emerge.

Recent advances in multiplex technology have pushed sensitivity into new territory. Modern platforms such as the **Luminex xMAP® System** can now achieve detection limits as low as 0.06 pg/mL. Widely applied for protein assays—including cytokines, chemokines, and growth factors—as well as for gene expression analysis, Luminex technology exemplifies how technical innovation directly expands research capability.



These advances translate into meaningful clinical insight. In a lung cancer biomarker study, researchers used an **R&D Systems Human XL Cytokine Luminex® Performance Assay** to measure both VEGF and PD-L1 in patient serum.<sup>3</sup> The panel delivered ELISA-comparable sensitivity while enabling simultaneous detection, revealing co-expression patterns linked to tumor angiogenesis and immune evasion. This dual readout provided a more complete picture of pathway interactions than single-biomarker approaches could achieve.

By analyzing VEGF and PD-L1 together, the study identified prognostic indicators of poor survival in lung adenocarcinoma while highlighting the therapeutic potential of targeting angiogenesis and immune evasion in tandem. Multiplex profiling exposed complex biological signatures within the tumor microenvironment that would remain invisible in isolation—underscoring how integrated biomarker analysis sharpens both prognostic insight and treatment strategy.

## Turning Complexity into Clarity

Cancer samples such as serum, plasma, urine, and tumor lysates present a uniquely difficult testing environment. They are rich in high-abundance proteins, interfering lipids, and heterophilic antibodies—factors collectively known as matrix effects. Left unaddressed, these interferences can mask true analyte signals, leading to false negatives, exaggerated positives, or loss of sensitivity.

One of the most common challenges lies in the imbalance between high- and low-abundance proteins. For example, IL-8 may be present at very high levels while IL-17 exists only in trace amounts. Without careful dilution and assay design, the stronger IL-8 signal can overwhelm IL-17, making accurate detection nearly impossible.

R&D Systems Multiplex Assays are engineered to overcome these barriers. Optimized diluents and buffers are specifically formulated to minimize nonspecific binding, reduce cross-reactivity, and control variability from factors such as pH, viscosity, or salt concentration. This ensures accurate spike recovery and linear dilution across a wide range of analytes—even in the most complex biological samples. Signal integrity is preserved for both high- and low-concentration proteins, delivering enhanced precision and lot-to-lot reliability across panels measuring dozens of biomarkers.

## Adapting Panels for Discovery and Validation

Cancer is not a static disease. As tumors evolve, so too does the need to monitor shifting patterns of protein expression. Because no two cancer studies are alike, assay flexibility is essential. R&D Systems addresses this need with two complementary formats:

- Discovery Assays offer customizable panels with access to more than 450 analytes, ideal for broad exploratory research where researchers cast a wide net to identify candidate biomarkers.
- High Performance Assays provide pre-validated panels optimized for sensitivity and reproducibility on par with ELISA, designed for translational and clinical applications where refinement and validation are critical.

This two-tiered model provides a natural progression: researchers begin with Discovery Panels to explore the biological landscape, then transition to High Performance Panels for focused validation as studies advance toward clinical endpoints.

Building on this foundation, R&D Systems recently introduced the Human Tumor Biomarker Performance Panel—a rigorously validated 28-plex assay optimized for plasma samples collected in preservative blood collection tubes. Leveraging Luminex xMAP technology, the panel encompasses tumor markers relevant to liver, ovarian, breast, pancreatic, neuroendocrine, and thyroid cancers. Each analyte has undergone stringent validation to ensure optimal performance in plasma-based applications.

Pairing validated multiplex panels with specialized sample collection methods further strengthens early detection workflows. Preservative tubes such as Streck® Cell-Free DNA BCT® and Protein Plus BCT™ maintain sample integrity by preventing genomic DNA release and preserving circulating tumor components critical for downstream analysis. Together, these advances bridge genomic and proteomic tools—bringing multi-cancer early detection and diagnostic development within closer reach.

# From Research to Clinical Impact: The Oncuria Story

The path from biomarker discovery to clinical diagnostics is notoriously difficult. It requires not only scientific insight but also platforms that combine sensitivity, reproducibility, and validated reagents capable of meeting regulatory standards. Bladder cancer diagnostics illustrate both the challenge and the opportunity.

Although relatively uncommon, bladder cancer has one of the highest recurrence rates among cancers. Current diagnostics are unable to reliably detect the disease before clinical presentation; nearly 30% of patients are first diagnosed at stage 2 or higher, where the five-year survival rate falls to around 50%. The clinical need for earlier, more accurate detection is clear.

To address this gap, Nonagen Bioscience developed Oncuria®, a first-of-its-kind multiplex protein test built with high-quality reagents from R&D Systems. Oncuria measures 10 proteins associated with bladder cancer in urine samples and has been validated in more than 4,300 patients. Clinical studies demonstrated 93% sensitivity and 95% specificity—giving urologists greater diagnostic confidence than existing urine-based methods.<sup>4</sup> (See Figure 1)

FIGURE 1.

	AUROC	Sensitivity	Specificity
Overall	0.95	0.93	0.93
Low-grade tumors	0.94	0.89	0.93
High-grade tumors	0.95	0.94	0.93
Low-grade tumors (NMIBC)	0.93	0.93	0.93
High-grade tumors (MIBC)	0.97	0.94	0.93

Figure 1. Diagnostic performance of Oncuria in patients with no personal bladder cancer history: presenting to a urology outpatient clinic for bladder cancer evaluation. 46 patients with de novo bladder cancer and 316 controls. Patients with no history of bladder cancer; n=362.

# Diagnostic Performance of Oncuria

This case underscores a broader reality: advancing from research use only (RUO) assays to laboratory-developed tests (LDTs) requires platforms that combine sensitivity, reproducibility, and validated reagents. Cancer research often spans years, which makes reproducibility across assay lots absolutely critical. Even small inconsistencies can compromise longitudinal data, undermining confidence in insights related to disease progression, relapse, or therapeutic response.

To ensure consistency, every R&D Systems assay panel undergoes rigorous validation—particularly important in longitudinal oncology trials, such as tracking cytokine release in patients undergoing immunotherapy. Because these studies often rely on limited clinical specimens, sample conservation is essential. Multiplex assays meet both requirements, enabling broad biomarker profiling from as little as 25–50 µL of serum or plasma, making repeated measurements possible even in constrained patient cohorts.

Trusted suppliers with deep immunoassay expertise and regulatory-grade manufacturing serve as essential partners in this process. Custom panel design further accelerates the RUO-to-LDT pathway, enabling researchers to save time, mitigate risk, and build assays capable of meeting clinical standards. The Oncuria example demonstrates how multiplex assays can transition from research tools to clinically actionable diagnostics—transforming biomarker discovery into real-world diagnostic impact.

## The Path Forward: Integration and Precision

One of the most promising aspects of protein biomarkers is their detectability in minimally invasive samples such as blood, urine, or saliva. This opens the door to routine monitoring and population-wide screening with far greater patient comfort and compliance. By reducing barriers to participation, non-invasive sampling makes it possible to follow high-risk individuals more closely and detect disease earlier.

Multiplex immunoassays extend this advantage by enabling sensitive, simultaneous measurement of multiple biomarkers from very small sample volumes. Dozens of analytes can be quantified in just hours using standard laboratory instrumentation—work that would otherwise take days if run as single-analyte assays. The result is faster data generation, streamlined workflows, and more timely decision-making in both clinical trials and translational research.

Early cancer detection demands more than a single molecular signal. Protein biomarkers and tumor-derived nucleic acids each provide valuable insights, but together—integrated within multi-omics approaches—they deliver a more complete view of cancer biology. Multiplex immunoassays sit at the center of this evolution, enabling systems-level analysis that supports earlier detection, prediction of therapeutic response, and personalized treatment strategies.

The hurdles of multiplexing—cross-reactivity, matrix effects, dilution challenges, and reproducibility—have always been part of the equation. Advances in antibody quality, assay design, and validation now make it possible to generate reliable, reproducible data even from complex clinical samples. With high sensitivity, broad analyte coverage, flexible panel formats, and decades of immunoassay expertise, R&D Systems Assays provide researchers with tools that are both rigorous and practical.

As cancer research enters a new era defined by precision and integration, multiplex immunoassays provide more than just technical solutions—they lay the foundation for earlier detection, more effective therapies, and ultimately, better patient outcomes.

## Key Considerations When Selecting Multiplex Assays

- **Antibody validation:** Build on well-characterized antibodies tested for cross-reactivity and proven for multiplex performance.
- **Panel-wide validation:** Choose panels validated as integrated systems, not just collections of single-analyte assays.
- **Matrix optimization:** Account for high- and low-abundance targets with proper grouping and optimized diluents.
- **Supplier expertise:** Partner with suppliers who combine antibody expertise and assay development experience to reduce risk and ensure reliable results.

## R&D Systems Luminex Assay Capabilities

- **Flexible formats:** Discovery and High Performance Panels designed for different research stages.
- **Validated quality:** Antibodies tested for specificity, panels validated for recombinant and natural proteins.
- **Efficient workflows:** Generate results in 3–3.5 hours using  $\leq 50 \mu\text{L}$  of sample.
- **Proven reproducibility:** Stringent lot-to-lot QC ensures consistent performance over time.

## REFERENCES

1. Charles J Rosser, MD, MBA, Hideki Furuya, Cedars-Sinai Medical Center, A Novel Multiplex Immunoassay for the Early Detection of Bladder Cancer, IIT2021-19-Furuya-EarlyDx, 2022 04-29
2. Yarahmadi A, Afkhami H. The role of microbiomes in gastrointestinal cancers: new insights. Front Oncol. 2024 Feb 1;13:1344328. doi: 10.3389/fonc.2023.1344328.
3. Journal, Cancer Research, Publisher: American Association for Cancer Research. PMID: 37074042; PMC Id: 10236157; DOI: 10.1158/0008-5472.CAN-21-4425
4. Furuya H, Tabula L, Lee R, Kralovec P, Ramsden M, Wong R, Rosser CJ. Analytical validation of ONCURIA™ a multiplex bead-based immunoassay for the non-invasive bladder cancer detection. Pract Lab Med. 2020 Nov 13;22:e00189. doi: 10.1016/j.plabm.2020.e00189.

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners.  
18300256523\_1125

**biotechne®** / Global Developer, Manufacturer, and Supplier of High-Quality Reagents, Analytical Instruments, and Precision Diagnostics.

INCLUDES

R&D Systems™

Novus Biologicals™

Tocris Bioscience™

ProteinSimple™

Asuragen™

ACD™

Lunaphore™