

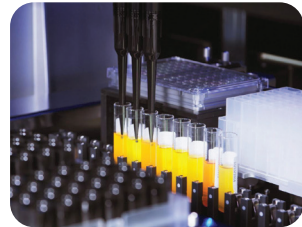
## Easy to learn, easy to use.

With a sleek, simple, flexible user interface and a monitoring dashboard that enables access from any current Web browser, these technologies provide the reliability and support laboratories can rely on.

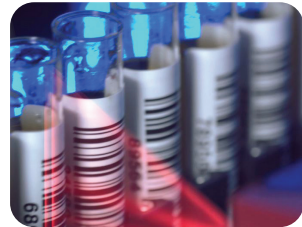
To start your run, simply load your instrument, following the intuitive guides in the software. The run starts automatically after the instrument detects that all samples, reagents, and consumables are in place. The integrated system software guides and monitors your workflow - including sample identification and processing, demographic data entry, sample classifications and quality assessment.

The performance dashboard enables monitoring of markers and assay performance, so you can check:

- Images of tubes, reagents, and consumables used
- Positioning of plates
- Number of samples per batch
- Assay performance statistics
- Quality control and review



Automated plasma pipetting helps avoid contamination.



Reagents and samples are barcoded to enable complete traceability.



Intuitive software guides you through the process.

# Analyze cfDNA samples in-house.

The Revvity logo, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

1246252

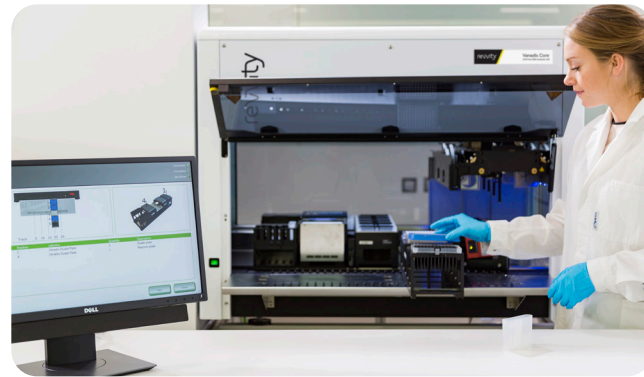
[www.revvity.com](http://www.revvity.com)

The Revvity logo, featuring the word "revvity" in a lowercase, sans-serif font.

Vanadis® cfDNA platform  
For research use only. Not for use in diagnostic procedures.

# Vanadis cfDNA platform

Vanadis® is the turnkey solution for analyzing the common chromosome abnormalities (21, 18, 13, X & Y). Eliminating PCR amplification and gene sequencing, Vanadis makes cfDNA analysis accessible to all labs. Better yet - Vanadis is equipped with walkaway automation, needing only one lab technician to analyze up to 14,000 samples per year.



## What makes the Vanadis platform unique?

**Breakthrough technology,** eliminating PCR and sequencing to provide cfDNA analysis to any laboratory

**Walkaway automation,** enabling a streamlined workflow for greater efficiency

**More precise assay** designed to minimize the failure rate

**A scalable platform,** allowing a single technician to run up to 14,000 samples per year

**Easy-to-use system** with full traceability of both samples and reagents

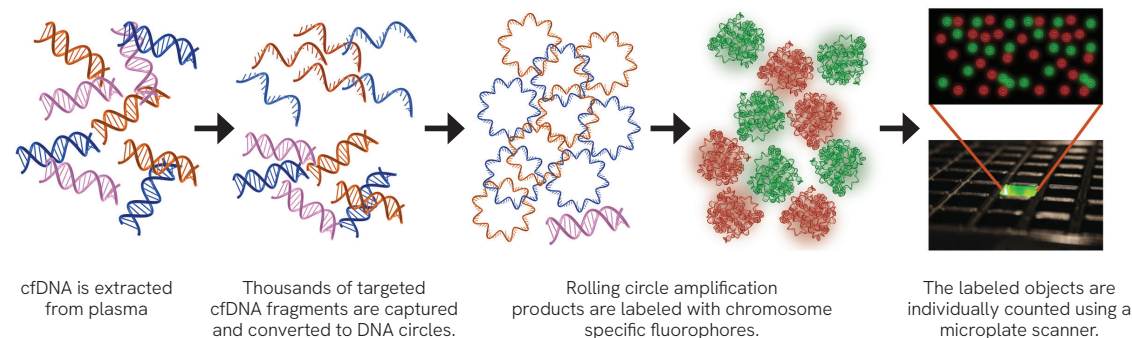
**Integrated system software** for remote workflow monitoring

## Groundbreaking cell-free DNA technology

Unlike complex, expensive sequencing-based platforms, Vanadis enables cost-efficient cell-free DNA analysis using standard microplates and automated sample processing - from primary tube to final data analysis. The precise Vanadis platform converts target chromosomes into rolling circle amplification products that simplify data analysis.

### No sequencing, no PCR

The Vanadis platform enables targeted cfDNA analysis without PCR, instead directly capturing target fragments of chromosomes 21, 18, and 13 and labeling them for counting. This eliminates the need for expensive and data-intensive steps such as DNA sequencing, microarrays, and microfluidics.

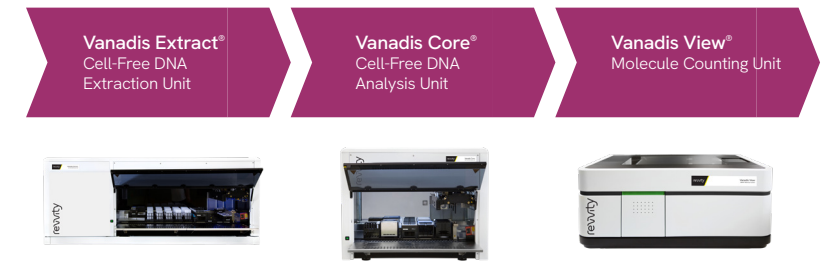


## Walkaway automation for greater efficiency

With the Vanadis technology, all critical steps are automated, starting with pipetting steps to reduce manual errors. Plasma volumes are monitored by camera to avoid contamination by the buffy coat, and samples & reagents are barcoded for complete tracking throughout the workflow. Software guides you through the process, with user-friendly touchscreens, and remote workflow management for ease of use and access.

## Automated workflow, from primary tube to final result

Results available in 72 hours\*



\* For 48 samples

## Minimum hands-on time

Apart from sample centrifugation steps, hands-on time can be as little as 40 minutes, depending on the number of samples being analyzed.

	Vanadis Extract®	Vanadis Core®	Vanadis View®	Total
48 samples	20 minutes	~20 minutes	3 minutes	~40 minutes
84 samples	40 minutes	~20 minutes	3 minutes	~60 minutes

## Scalable capacity to support evolving needs

The Vanadis platform was developed to offer maximum flexibility to laboratories whose workload encompasses from 2,000 to 14,000 determinations per year. What's more, with no minimum number of samples required to start extraction, labs can start using Vanadis right away.

## Precise assay for minimum failure rate

Assay precision describes how much the measured ratio for normal samples varies, and is highly influenced by the design of the assay (PCR, counting statistics, level of automation...).

The precision of the Vanadis assay greatly reduces variation and tightens the Chromosomal Ratio distribution. As shown in the figure to the right, Vanadis enables the detection of small Chr21 elevations and minimizes the failure rate.

