



Analyze chromosomes 21, 18, 13, X and Y with ease.

Vanadis® cfDNA platform

Revvity is proud to introduce the easy-to-use and cost-effective **Vanadis cfDNA platform**. Analyzing the common chromosome abnormalities (21, 18, 13, X & Y), Vanadis automates all critical steps, from primary tube to final data analysis, without the use of PCR or sequencing.

Easy-To-Use

Determinations are available in as soon as 72 hours*. Without PCR or sequencing, Vanadis does not require bioinformatics or genetic expertise to operate.

Cost-Effective

Requires only one lab technician to operate. Vanadis avoids costly IT infrastructure and or pre/post PCR room requirements, common across other cfDNA platforms.

Precise

Built to minimize the failure rate. Vanadis captures fragments of target chromosomes to sense the smallest chromosome elevations, bringing more determinations to your lab.

For research use only. Not for use in diagnostic procedures.
*Dependent on sample volume.



Vanadis View®

Not complex – just high tech.

Taking the complexity out of cfDNA analysis, our mission is to make cfDNA accessible to any laboratory. Wherever you're starting from, Vanadis can assist.

Learn more at www.revvity.com.