

60-second research spotlight: prenatal screening for autosomal trisomies



Study:

Article title: Cell-Free DNA Screening For Common Autosomal Trisomies Using Rolling- Circle Replication In Twin Pregnancies

Publication: PRENATAL DIAGNOSIS

Author: Vivanti, AJ *et al*

Publication date: May 27, 2024

Summary

Prospective 2-center study of 862 twin pregnancies. Compared no-call rate and performance against 16,382 singleton pregnancies. Twin samples analyzed on Revvity's Vanadis™ NIPT System showed great performance. While Vanadis™ NIPT System twin samples reported a higher no-call rate than singletons, its twin no-call rate was better than those of competing cfDNA technologies.

[Read here](#)

Products may not be licensed in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability

The views and opinions expressed herein are those of those of the study and do not necessarily reflect the views of Revvity, Inc.

Key takeaways:

1. The study found that using cell-free DNA (cfDNA) as a first-line screening test for common aneuploidies in both singletons and twins is better than biochemical screening.
2. The no-call rate, at first attempt, for common aneuploidy in twins is lower with Vanadis technology (2.0%) compared to other cfDNA technologies (2.9 - 13.2%).
3. In twin pregnancies, the Vanadis™ NIPT System performed strongly with a sensitivity of 100% (25 out of 25 pregnancies) and no false negatives recorded. Note: the number of affected samples in this study was limited, so the performance observed may not be typical.
4. IVF increased the probability of a no-call result in twin pregnancies, but maternal weight or gestational age did not have an impact.
5. The study suggests that the Vanadis™ technology NIPT System is less technically and analytically complex, offers a more attractive cost/benefit ratio when compared to sequencing technologies, and may be a viable alternative to other cfDNA technologies due to recent worldwide reagent shortages.