

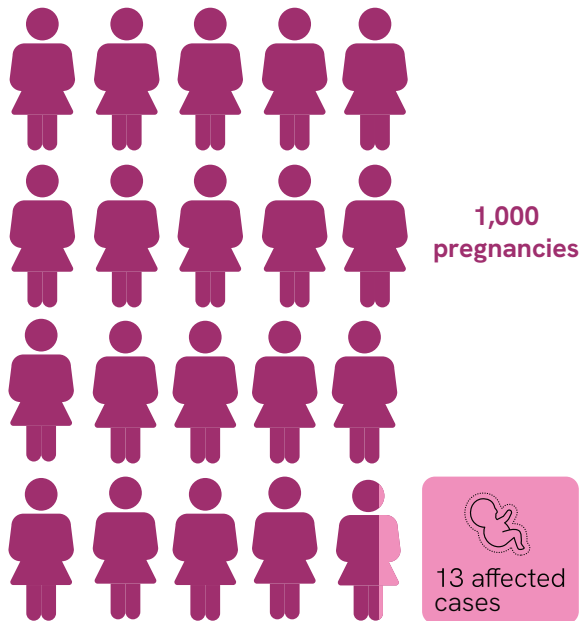


# NIPT: why no-call rates matter.

In clinical contexts, a noninvasive prenatal test (NIPT) with a low no-call rate delivers more results to women, reducing their chances of anxiety, financial stress<sup>1</sup>, or potential pregnancy complications.<sup>2</sup>

## Example

A hypothetical NIPT screening is conducted on **1,000 pregnancies** to detect a condition with 1.3% prevalence. Which test would you choose?



## Test A has a 4% no-call rate



40 no-call results

While 39/40 or 98% were unaffected, they still may experience



Anxiety



Financial burden<sup>1</sup>



Pregnancy loss<sup>2</sup>

96% of women receive a result



12 affected cases reported



948 unaffected cases reported

## Test B has a 0% no-call rate

100% of women receive a result



13 affected cases reported



987 unaffected cases reported

NIPTs with lower no-call rates deliver more results to more women, minimizing the chances a patient may experience anxiety, financial strain<sup>1</sup>, or be at risk for pregnancy loss.<sup>2</sup>

Revvity, Inc. does not endorse or make recommendations with respect to research, medication, or treatments. All information presented is for informational purposes only and is not intended as medical advice.

\*Akolekar, R., et al. (2015) Ultrasound Obstet Gynecol, 45: 16-26. doi.org/10.1002/uog.14636

1. In the form of follow-up consultations and further testing
2. Due to invasive procedures.\*