



# Improving sample tracking in PGT-A.

Effective sample tracking and contamination detection are essential to prevent embryo misclassification. Sample tracking mistakes can lead to the discarding of viable or transferring of aneuploid embryos. Recent studies have highlighted that contamination from external sources effects on average 0.4% of biopsy samples\*.

## Take traceability to the next level

When combined with the innovative [Cap2™ tubes](#) from Azenta® Life Sciences, the [PG-Seq™ Rapid v2 kit](#), provides an efficient and error-minimizing solution for PGT-A. With robust sample tracking and a streamlined workflow, this workflow takes traceability of embryos to the next level.

\*Clark, G., et al. (2023). *Human Reproduction*, 38(Supplement\_1), dead093.337

Scan the QR codes to learn more:



The Azenta Cap2™



PG-Seq™ Rapid v2 kit

### Key features:

#### Azenta® Cap2™ tubes

- Sample collection and PCR in the same tube
- 2D matrix code and human readable ID
- Compatible with automated capping/decapping



#### Revvity PG-Seq™ Rapid kit v2

- Fast and streamlined workflow with minimal tube openings
- Enables the monitoring of external DNA contamination and sample swapping by analyzing mitochondrial DNA information from data obtained during the standard workflow
- Flexible throughput and resolution

