



Every sample matters.

Extract analytes from tough samples with ease.

Sample homogenization is a critical and universal step that starts diverse pathogenic organism or virus identification workflows. A robust, automation-enabled solution lends itself to analytical reproducibility and reduction-of-ambiguity in data results.



Tough & complex sample matrix processing

- Sample homogenization
- Cell and viral lysis

DNA/RNA extraction

- Automated nucleic acid isolation

Automated workflows & molecular assays

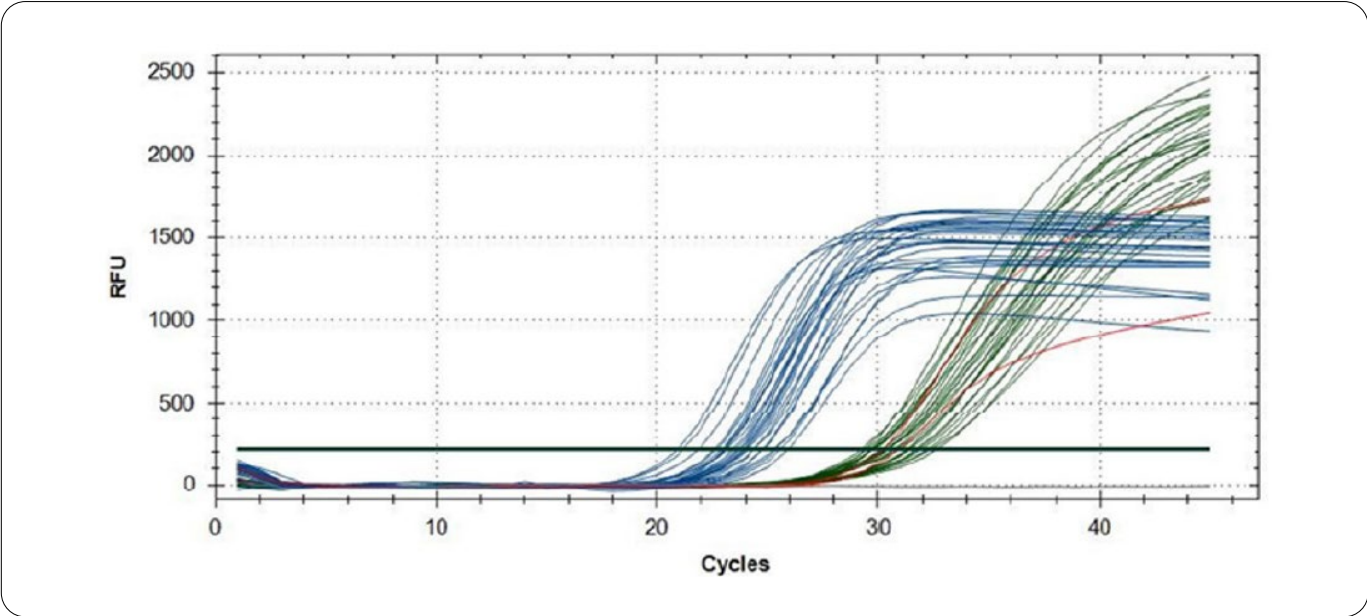
- Microarrays
- Multiplex PCR assays
- Next-gen sequencing (NGS)



Routine sample matrices

- Nail shavings
- Skin
- Hair stubs
- Cultured material
- Liquid or tissue biopsy samples
- Histological sections
- Wastewater filters

Leverage our BR cryo-cooling unit (Cat # 19-8005) with the Omni Bead Ruptor Elite™ bead mill homogenizer (Cat # 19-042E) for temperature-sensitive biological analytes.



RT-PCR method using saliva samples into a bead mill homogenization workflow compared to standard naso-pharyngeal swab PCR methods. Results demonstrated 100% agreement between methodologies. Blue: SARS-CoV-2 N1/N2 gene, Green: RNase P (internal control), Gray: negative control, Red: positive control. (n = 60)

Avoid potential headaches and invest in lab fundamentals

- Future-proof scalability with measurable operational savings
- Path to standardization to maximize efficiency and throughput

Additional resources

APPLICATION NOTE

Molecular identification of dermatophytes in dermatomycosis

APPLICATION NOTE

Direct-to-PCR method for influenza A virus detection

APPLICATION NOTE

Epigenetic analysis of tissues via ChIP-seq

Rely on our Omni Bead Ruptor series of bead mill homogenizers to streamline workflows and help optimize protocols. Explore more at www.revivity.com/category/bead-mill-homogenizers

