

# Simplifying failure analysis with plasma-FIB

The right tool for the right job

Cut precision without compromising throughput

Allowing efficient bulk milling, controlled end-pointing, and high-quality polishing, resulting in maximal throughput and success rate.

FEATURE DENSITY

Mechanical polish

*Ex situ* laser

*In situ* laser

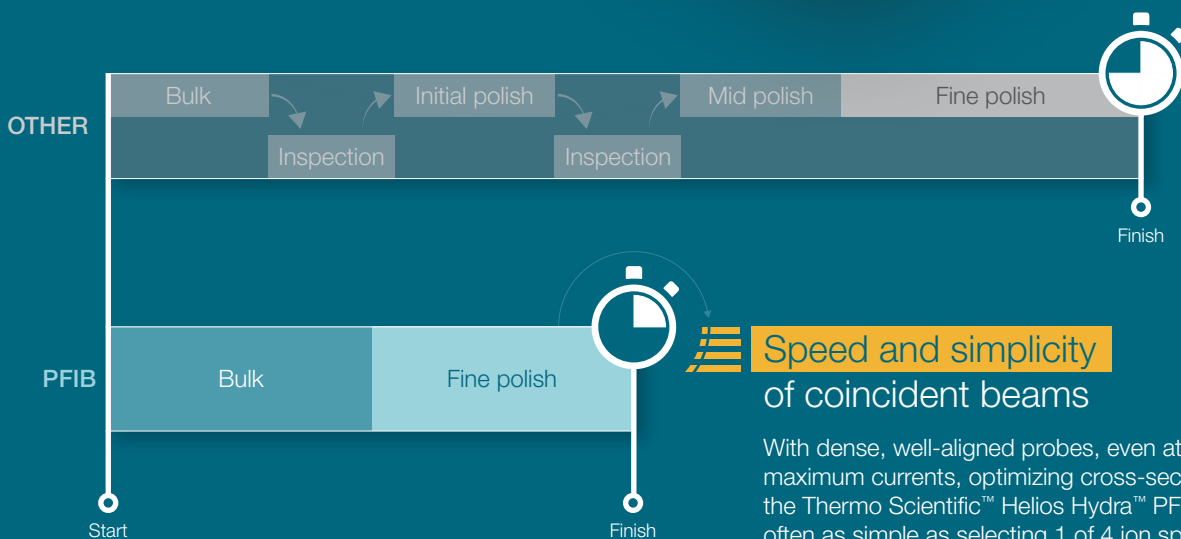
Coincident laser

PFIB

Pinpoint easy-to-miss targets buried deep within complex 3D structures

For small defects or features, the precision of PFIB milling helps ensure that you don't completely miss or even destroy your target.

Faster results in fewer steps



Speed and simplicity of coincident beams

With dense, well-aligned probes, even at maximum currents, optimizing cross-sections with the Thermo Scientific™ Helios Hydra™ PFIB-SEM is often as simple as selecting 1 of 4 ion species.

Learn more at [thermofisher.com/helioshydra](https://thermofisher.com/helioshydra)