

# Krios G4

Designed for true atomic-resolution cryo-EM and speed

**Autoloader**  
Samples can be retrieved and transferred contamination free to another autoloader equipped microscope for re-investigation – optimal tool connectivity

**E-CFEG**  
Cold field emission gun optimized for low energy spread ( $\leq 0.3$  eV) enables higher contrast at highest resolutions ( $\leq 2.0$  Å) in a shorter time

**Selectris X Imaging Filter**  
Stable, zero-loss energy filtering for atomic-resolution cryo-EM

**Falcon 4i Direct Electron Detector**  
Speed, productivity, and efficient data compression without compromising image quality