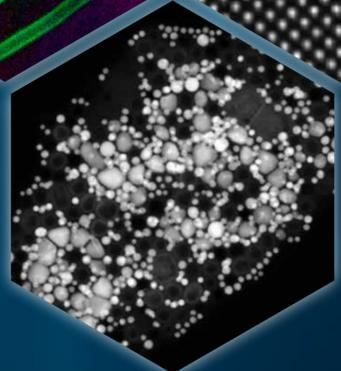
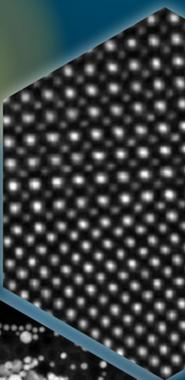
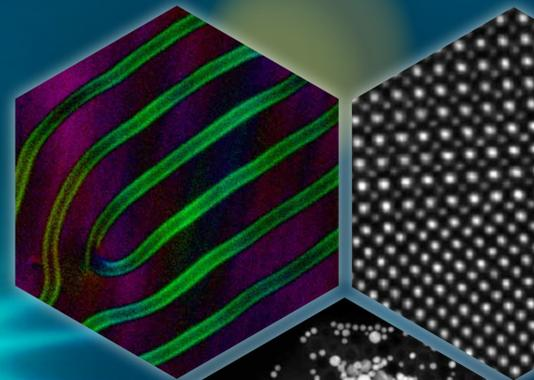
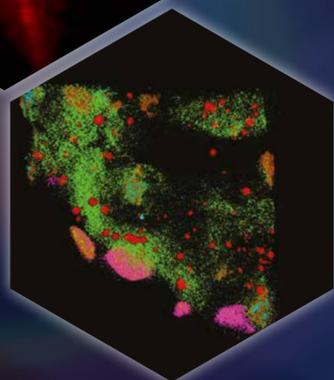
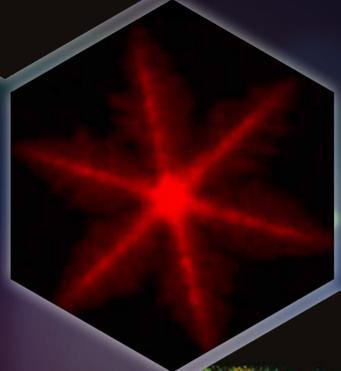
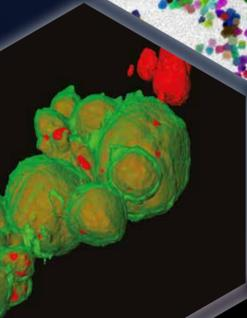
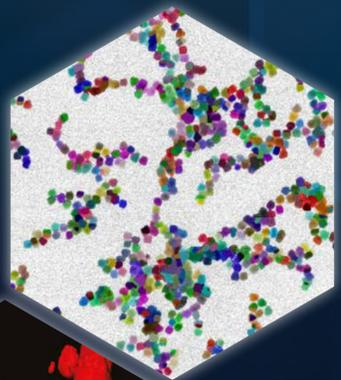
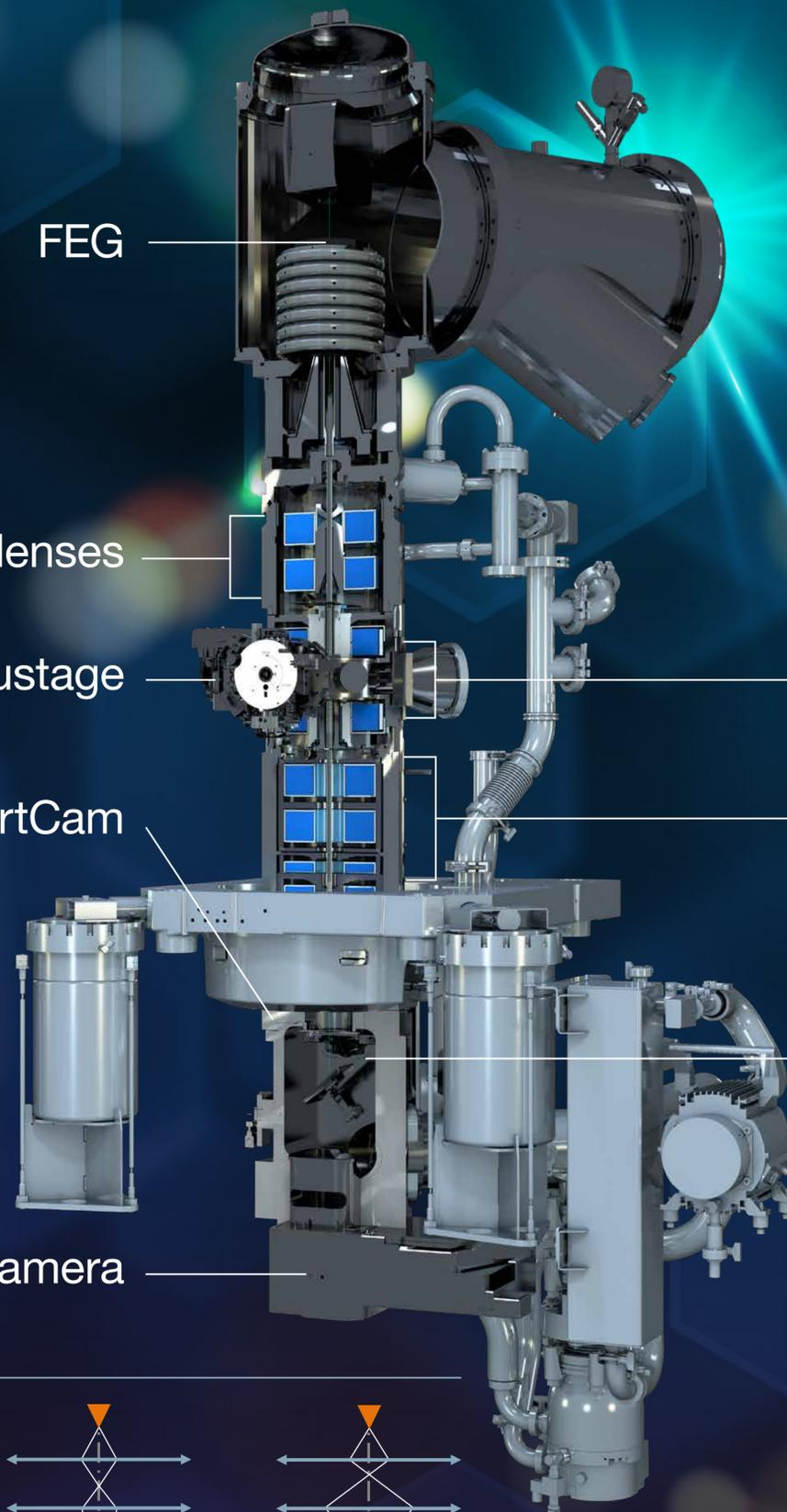


Thermo Scientific™ Talos™ Scanning Transmission Electron Microscope

Elevate Your Analysis Without Compromise

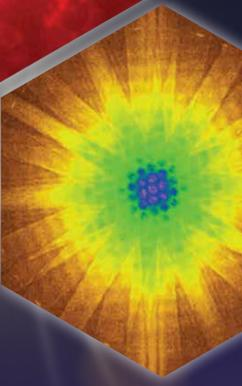
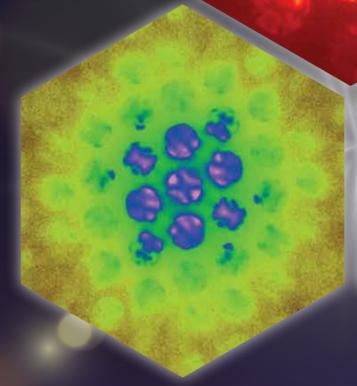
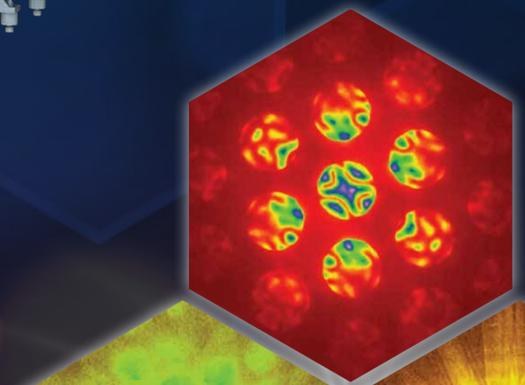
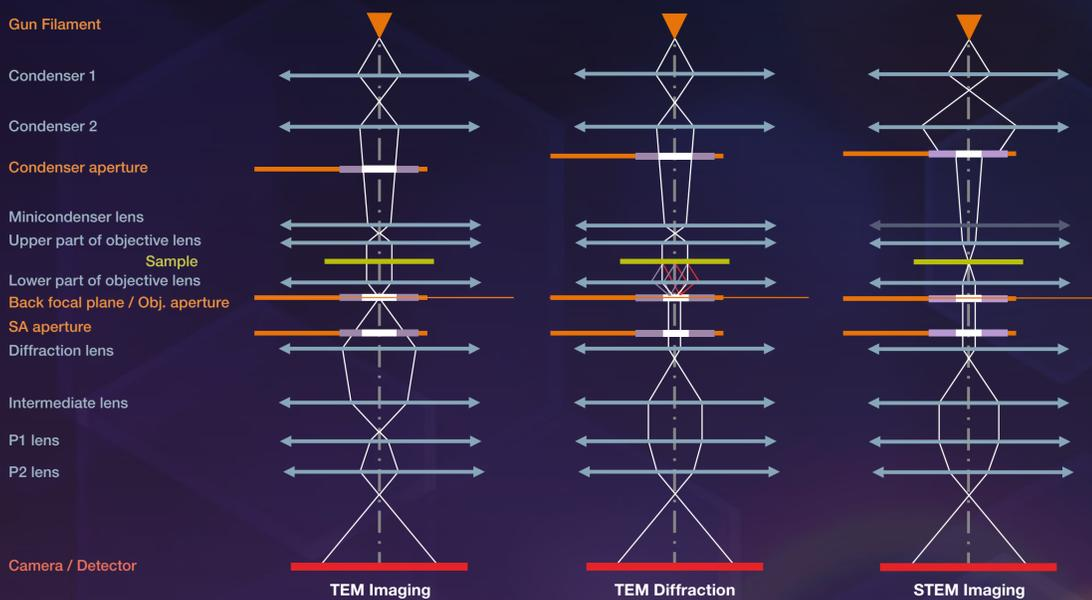


Constant power objective lenses

Projector system

Flu-screen

TEM modes



Learn more at thermofisher.com/tem

thermo scientific