

Thermo Fisher Scientific (Bremen) GmbH, Hanna-Kunath-Str. 11, 28199 Bremen, Germany

Learn more at thermofisher.com/orbitrapastralzoom