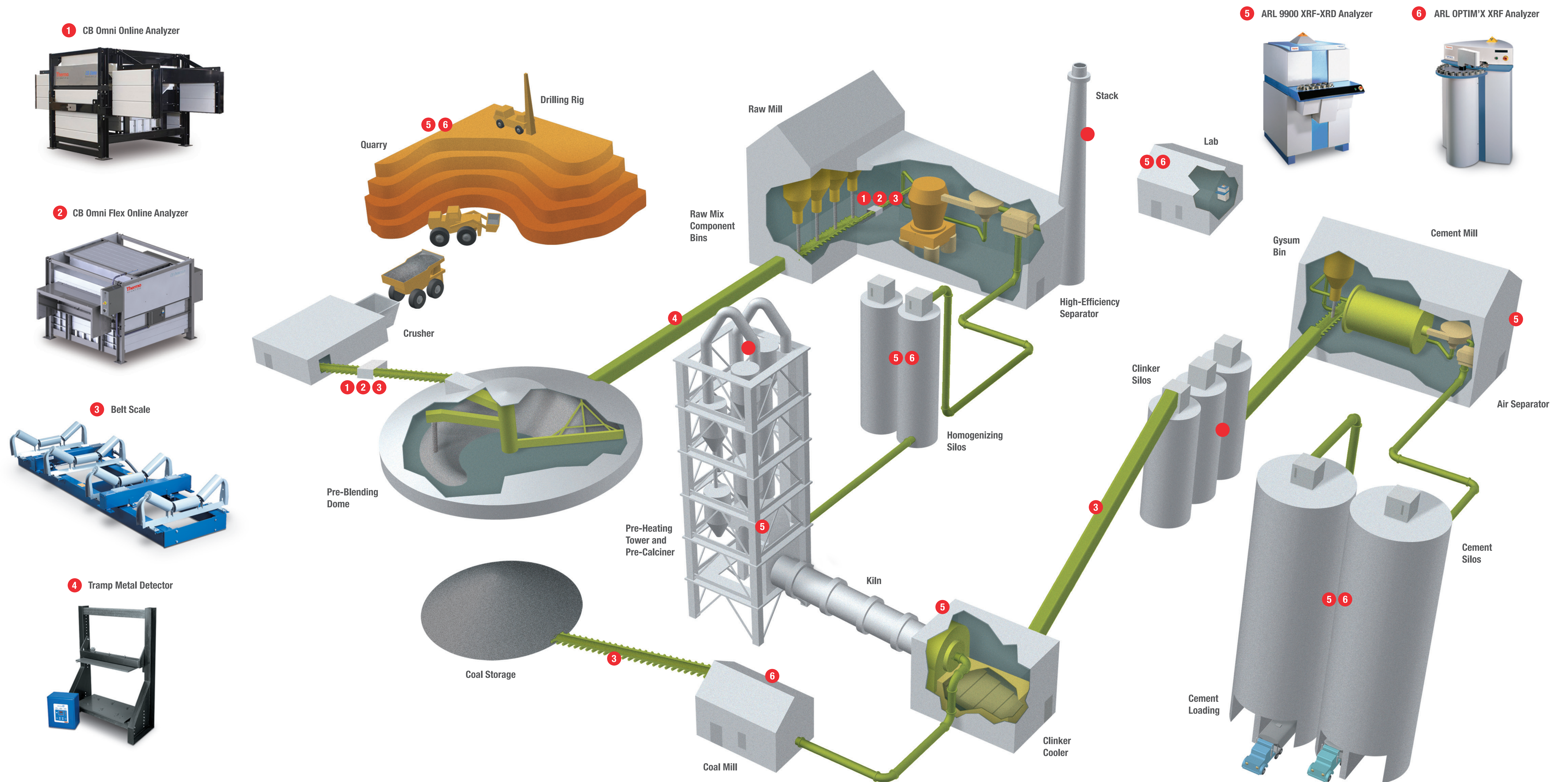


Thermo Scientific Products in the Cement Manufacturing Processing



Innovative Products for Your Cement Process

We continue to lead the way with innovative and unique product solutions that help you meet your profitability targets. We design, manufacture and service highly reliable specialized equipment for your industry and are dedicated to delivering equipment to meet your needs. The large portfolio of products includes elemental online analyzers, X-ray analyzers, belt scales, weigh belt feeders, level sensors and indicators, flow detectors, impact weighers, stack emission gas detectors, material storage tracking software and more.

Experience first-hand why the Thermo Scientific Brand represents the world's leading analytical instruments.

● Thermo Fisher Products are Useful at this Stage