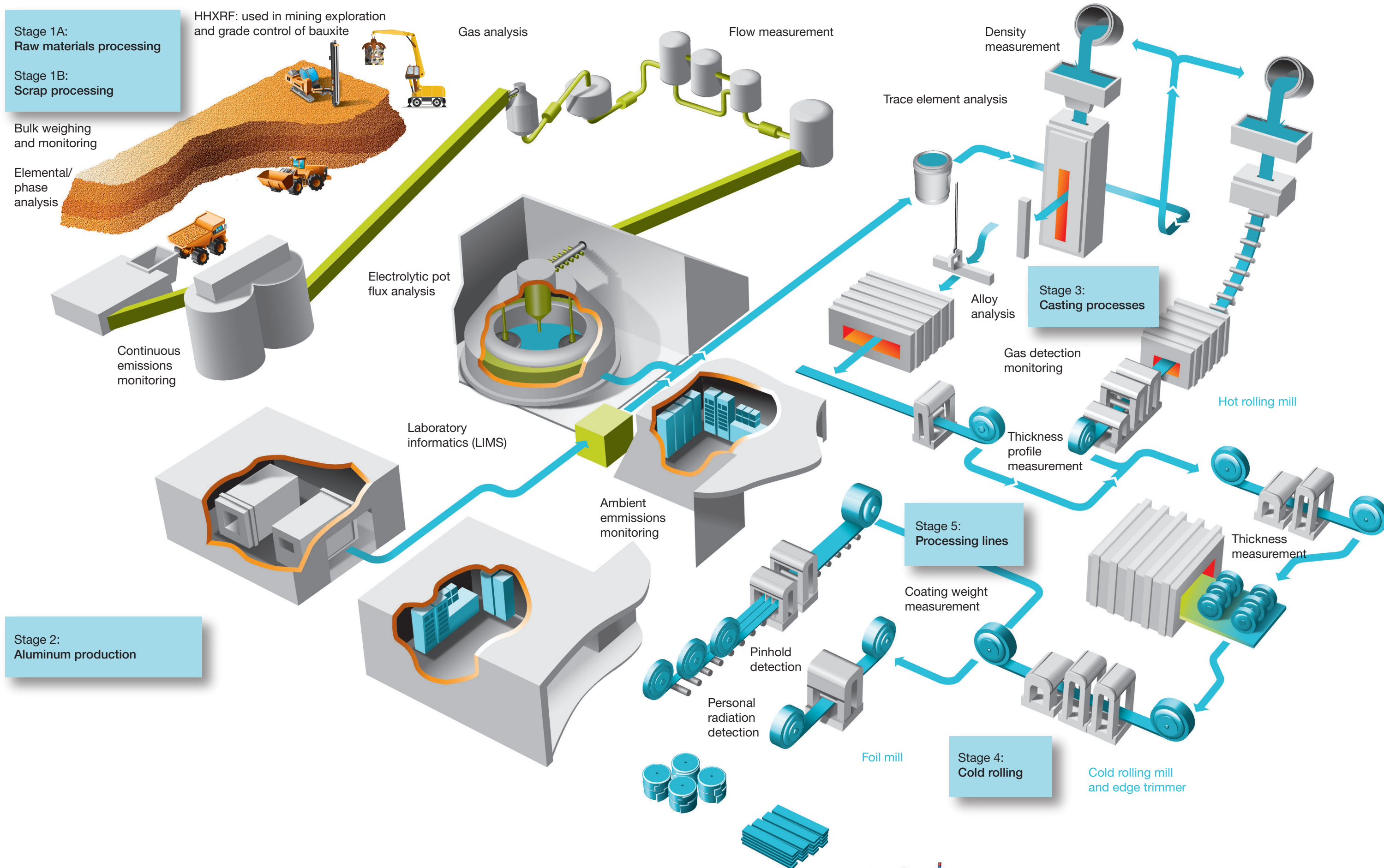


Aluminum manufacturing process map



Stage 2: Aluminum production

Stage 1B, 2, & 5
 Thermo Scientific™
 XL5 Plus Handheld XRF Analyzer

Stage 1A
 Thermo Scientific™
 ARL™ X'TRA Companion

Stage 2
 Thermo Scientific™
 ARL™ 9900 IntelliPower™

Stage 3
 Thermo Scientific™
 ARL™ iSpark™ Plus

Stage 4
 Thermo Scientific™
 ARL™ SMS-3300

Stage 4
 Thermo Scientific™
 RM 210 AS Aluminum Thickness Gauge

Stage 5
 Thermo Scientific™
 RM 300 EL Coating Weight Gauge

Learn more at thermofisher.com/aluminum