

A large, black and white industrial machine with a yellow stripe, labeled 'Quattro'. It has a control panel on the right side with a screen and buttons. The machine is designed for automated material processing.

Component	Part Number	Lot Number
Calibrator Concentrate	105630	430514
Control 1 Concentrate	105632	430516
Control 2 Concentrate	105633	430517

Review/Approval

Dhwani Patel QA Specialist II DP / 11 Nov 2025

Name Title Signature/Date