

Performance and Design for any Environment



HFC Scan - Hand & Foot Contamination Monitor

- Very high sensitivity
- Highly robust for use in industrial applications
- Integrated transport wheels for ease of mobility
- Easy decontamination
- One detector type for hands, feet and clothing
- Detachable hand detector
- Intuitive touch display

SZ345	Background count rate [s ⁻¹]	Efficiency [%] Foot/Hand*	MDA [Bq/cm ²] Foot/Hand*
¹⁴ C	< 15 (β)	8/14	0.76/0.42
³⁶ Cl		31/51	0.10/0.06
²⁴¹ Am	< 0.1 (α)	25/48	0.03/0.02

*ISO 7503