

Company		Contact Name	
Street/P.O.Box		E-Mail	
City		Phone	
Postal code		Date	
Country / State		Project	

Process Specifications

Measuring point		Measuring range		Unit ☐ K ₂ O ☐ KCL other
Material condition	☐ bulk ☐ liquid			

	Unit (if other, please specify)	normal	min.	max.
Bulk density	g/cm ³			
Particle size	mm			
Product temperature	°C			
Ambient temperature	°C			

Vessel Specifications

☐ Measurement inside the vessel (Dip tube)	Vessel form		Wall thickness	
☐ Measurement outside the vessel	External dimensions (please attach drawings)		Measuring position (please attach drawings)	

Belt Conveyor Specifications

	Unit (if other, please specify)	normal	min.	max.
Conveyor speed	m/s			
Loading height (H)	mm			

Smoothing with scraper to get a constant bulk height possible? ☐ no ☐ yes on mm (if other unit, please specify)

Speed signal ☐ constant speed ☐ 4 ... 20 mA signal available ☐ Tachometer required

Accuracy %

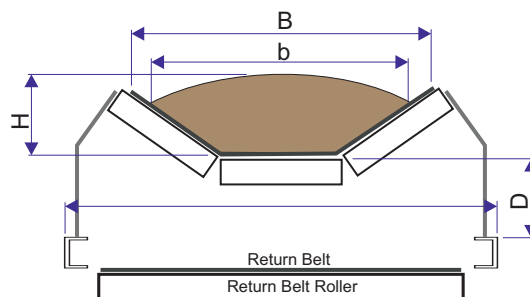
Belt material

Belt thickness

Belt width (B)

Typical loading width (b)

Distance belt to carrier (D)



Instrumentation

Power supply ☐ 90-250V AC/DC ☐ 24V AC/DC

Current output: ☐ Active/Source ☐ Passive/Sink