

Company

Street/P.O.Box

City

Postal code

Country / State

Contact Name

E-Mail

Phone

Date

Project

Process Specifications

Measuring tag Product

Application

Material composition

	Unit (if other, please specify)	normal	min.	max.
Bulk density	g/cm ³ <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Particle size	mm <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Product temperature	°C <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Ambient temperature	°C <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Conveying capacity	t/h <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Conveying speed	m/s <input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Speed signal	<input type="checkbox"/> constant speed <input type="checkbox"/> 4 ... 20 mA signal available <input type="checkbox"/> Tachometer required			
Accuracy	<input type="text"/> % of nominal load	t/h (+/-) <input type="text"/>	<input type="text"/> % at nominal load	

Belt Conveyor



Belt material

mm (if other, please specify)

Normal loading height (H)

Min. loading height (H)

Max. loading height (H)

Belt width (B)

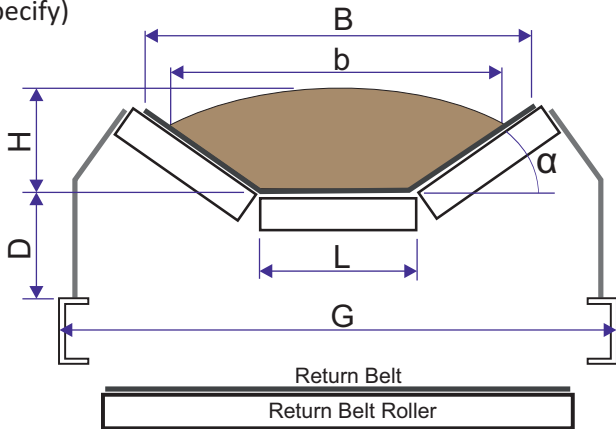
Typical loading width (b)

Distance belt to carrier (D)

Carrier distance outside (G)

Middle role width (L)

Pitch angle (α)



Please add drawing/photo of the belt conveyor with cross section and side view and indicate the possible installation place of the measuring frame

Trough Chain Conveyor

Material of scraper and trough

mm (if other, please specify)

Material height (h)

Trough width (B)

Trough height (H)

Scraper width (A)

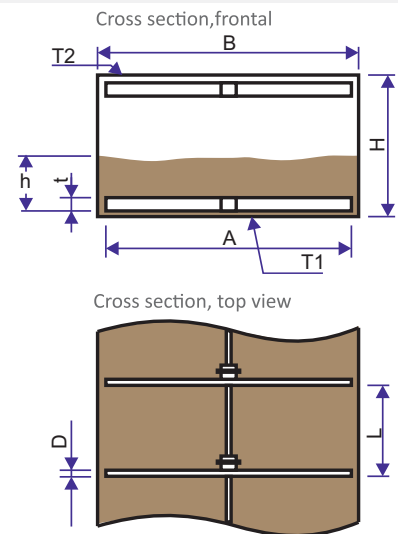
Scraper height (t)

Scraper thickness (D)

Sheet thickness bottom (T1)

Sheet thickness top (T2)

Distance between scrapers (L)



Screw Conveyor

Material of screw and pipe

mm (if other, please specify)

Internal pipe diameter (D1)

Shaft diameter (d)

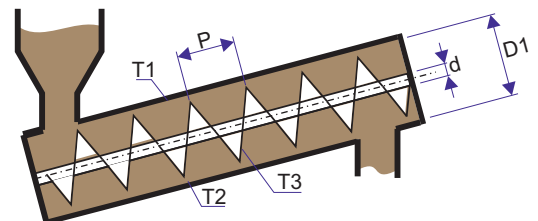
Housing thickness, top (T1)

Housing thickness, bottom (T2)

Blade thickness (T3)

Pitch (P)

RPM speed



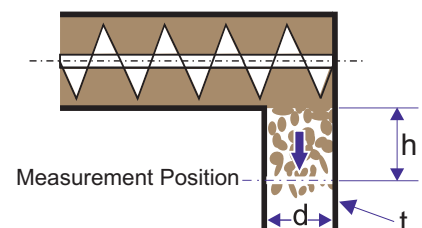
Free Fall Application

mm (if other, please specify)

Minimum distance to dropping point (h)

Pipe diameter or convey width (d)

Wall thickness (t) and material



Instrumentation

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

Exproof requested ☐ No ☐ Yes Type

Distance detector to evaluation unit m

What is more important? ☐ Price ☐ Accuracy

The products that Berthold Technologies offers are custom engineered systems. There are multiple family models and component options that are able to be selected based on the customer's process parameters. Also nuclear source sizes are calculated and selected for each individual system. These inputs are necessary to engineer a system that will meet the required needs and will function properly. Inaccuracies or omissions of the inputs could have a negative effect on the operation of the measurement. Berthold cannot be held accountable for the performance of their equipment if initial specifications were falsified or not presented fully.