

Company
 Street/P.O.Box
 City
 Postal code
 Country / State

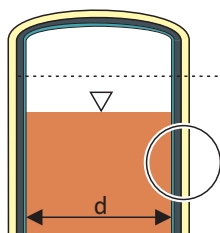
Contact Name
 E-Mail
 Phone
 Date
 Project

Process Specifications

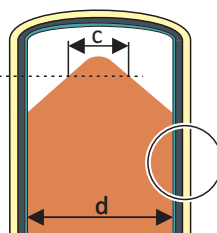
Measuring tag
 Application/Process

Vessel Specification

☐ Fluid level in vertical cylindrical vessel



☐ Bulk cone in vertical cylindrical vessel



☐ other
(please attach drawing)

Point of
measurement

Measuring range

☐ mm ☐ inch

Inner diameter (d) at point of measurement

Bulk cone diameter (c)

Thickness

☐ mm ☐ inch

Material

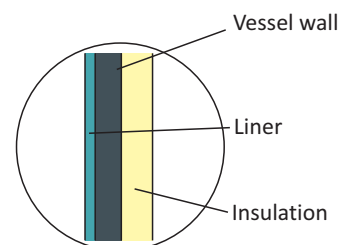
Density

g/cm³

Vessel wall

Liner wall

Insulation wall



Others (refractory layers, cladding, ...)

Structural Obstructions (agitator, collar, ...) ☐ No ☐ Yes if yes, please add drawing

Build Ups ☐ No ☐ Yes if yes, approximate thickness and density

Switching alarm function ☐ high ☐ low

Are there additional radiometric measurements close by?

☐ No ☐ Yes if yes, please add location plan

Product

<input type="radio"/> fluid <input type="radio"/> solid	Unit (if other, please specify)	normal	min.	max.	Name
Product/Bulk density	g/cm³				
Gas, foam, second fluid (if any)	g/cm³				
	g/cm³				
Vessel pressure	bar				
Product temperature	°C				

Instrumentation

Response time (s)		min.	max.	Unit (if other, please specify)
Ambient temperature at measuring point				°C
Power supply	<input type="radio"/> 90-250 V AC <input type="radio"/> 24V AC/DC			
Exproof requested	<input type="radio"/> No <input type="radio"/> Yes	Type		
Process signal:	<input type="radio"/> 4 ... 20 mA <input type="radio"/> HART <input type="radio"/> Relay <input type="radio"/> OFF <input type="radio"/> PA (FF = Foundation Fieldbus, PA = Profibus PA)			
Functional safety:	<input type="radio"/> none <input type="radio"/> SIL 2 <input type="radio"/> SIL 3			
Current output:	<input type="radio"/> Active/Source <input type="radio"/> Passive/Sink			

Retrofit (with existing source)

Original source date	
Original source activity	<input type="radio"/> Mbq <input type="radio"/> mCi
Type of isotope	
Radiation angle of shielding (degree)	
Supplier of source	

Please add drawing or at least a sketch of the existing installation with side and top view.

Comments / Special Requirements

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.