

Company	<input type="text"/>	Contact Name	<input type="text"/>
Street/P.O.Box	<input type="text"/>	E-Mail	<input type="text"/>
City	<input type="text"/>	Phone	<input type="text"/>
Postal code	<input type="text"/>	Date	<input type="text"/>
Country / State	<input type="text"/>	Project	<input type="text"/>

**Process Specifications**

Measuring tag

Application/Process

Product

**Measurement Task** (Please mark if density measurement or concentration measurement is required)

In case of a density measurement (e.g. g/cm<sup>3</sup>, SG):

**Density min/max**

	Unit (if other, please specify)	normal	min.	max.
Density	g/cm <sup>3</sup>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Product operating temperature	°C	<input type="text"/>	<input type="text"/>	<input type="text"/>
Do you know the chemical formula? <input type="text"/>				
Accuracy	Unit (if other, please specify)			
<input type="text"/>	%	<input type="text"/>		

In case of concentration measurement (e.g. %):

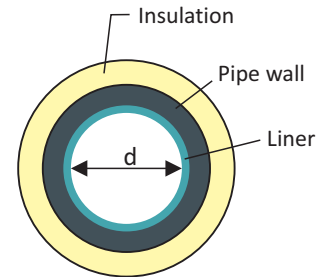
**Concentration (composits of 2 products)**

Name of product 1 (i.e. carrier liquid)

	Unit (if other, please specify)	normal	min.	max.
Density of product 1	g/cm <sup>3</sup>	<input type="text"/>	<input type="text"/>	<input type="text"/>
Name of product 2 (i.e. solids) <input type="text"/>				
	Unit (if other, please specify)	normal	min.	max.
Density of product 2	g/cm <sup>3</sup>	<input type="text"/>	<input type="text"/>	<input type="text"/>
	Unit (if other, please specify)	normal	min.	max.
Concentration (product 2 in product 1)	%	<input type="text"/>	<input type="text"/>	<input type="text"/>
Accuracy	Unit (if other, please specify)			
<input type="text"/>	%	<input type="text"/>		

**Pipe Specification**

	Thickness <input type="checkbox"/> mm <input type="checkbox"/> inch	Material	Density g/cm <sup>3</sup>
Inner diameter (d)	<input type="text"/>		
Pipe wall	<input type="text"/>	<input type="text"/>	<input type="text"/>
Liner wall	<input type="text"/>	<input type="text"/>	<input type="text"/>
Insulation wall	<input type="text"/>	<input type="text"/>	<input type="text"/>
Others	<input type="text"/>	<input type="text"/>	<input type="text"/>



Operating pressure in pipe     constant  bar  
     variable from  to  bar

Measurement on     horizontal pipe  
                           vertical pipe  
                           not on a pipe (e.g. vessel, tank or chute)  
                          Please attach drawing of the geometry, wall thickness and dimensions.

**Instrumentation**

	min.	max.	Unit (if other, please specify)
Ambient temperature at measuring point	<input type="text"/>	<input type="text"/>	°C <input type="text"/>
Power supply	<input type="radio"/> 90-250VAC <input type="radio"/> 24VDC		
Exproof requested	<input type="radio"/> No <input type="radio"/> Yes    Type <input type="text"/>		
Process signal:	<input type="radio"/> 4 ... 20 mA <input type="radio"/> HART <input type="radio"/> FF <input type="radio"/> 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     Unit (if other, please specify) Mbq

Type of isotope   

Supplier of source   

**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.