

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

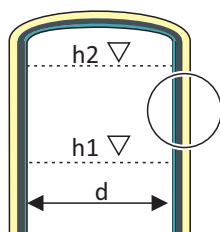
Contact Name   
 E-Mail   
 Phone   
 Date   
 Project

## Process Specifications

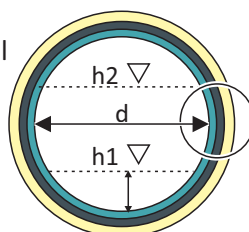
Measuring tag   
 Application/Process

## Vessel Specification

☐ Vertical



☐ Horizontal



☐ other (conical, conical-cylindrical, half conical, ...) **please attach drawing**

Measuring range

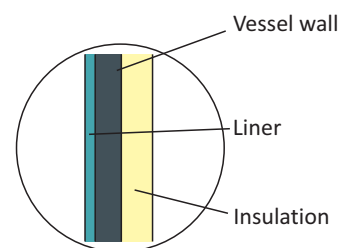
☐ mm ☐ inch

Product level at 0% (h1)

Product level at 100% (h2)

Inner diameter (d)

|                 | Thickness                                                 | Material             | Density                 |
|-----------------|-----------------------------------------------------------|----------------------|-------------------------|
|                 | <input type="checkbox"/> mm <input type="checkbox"/> inch |                      | <b>g/cm<sup>3</sup></b> |
| Vessel 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 (refractory layers, cladding, ...)

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

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

Are there additional radiometric measurements close by?

☐ No ☐ Yes if yes, please add vessel location plan

## Product

| <input type="radio"/> liquid <input type="radio"/> solid | Unit<br>(if other, please specify) | normal | min. | max. | Name |
|----------------------------------------------------------|------------------------------------|--------|------|------|------|
| Product density                                          | g/cm <sup>3</sup>                  |        |      |      |      |
| Gas, foam, second<br>liquid phase (if any)               | g/cm <sup>3</sup>                  |        |      |      |      |
|                                                          | g/cm <sup>3</sup>                  |        |      |      |      |
| Vessel pressure                                          | bar                                |        |      |      |      |
| Product temperature                                      | °C                                 |        |      |      |      |

## Instrumentation

|                                        | min.                                                                                                                                                           | max. | Unit<br>(if other, please specify) |
|----------------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------|------|------------------------------------|
| Ambient temperature at measuring point |                                                                                                                                                                |      | °C                                 |
| Power supply                           | <input type="radio"/> 90-250VAC <input type="radio"/> 24VDC                                                                                                    |      |                                    |
| 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"/> FF <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                  |                                                                     |
| Source Type                           | <input type="radio"/> Point Source <input type="radio"/> Rod Source |
| 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.