

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

Contact Name
 E-Mail
 Phone
 Date
 Project

Vessel Specification (please attach vessel drawing)

Vessel ☐ new/design phase ☐ existing

Vessel orientation ☐ horizontal ☐ vertical ☐ other

Vessel inner diameter ☐ mm ☐ inch

Wall thickness ☐ mm ☐ inch

Vessel material

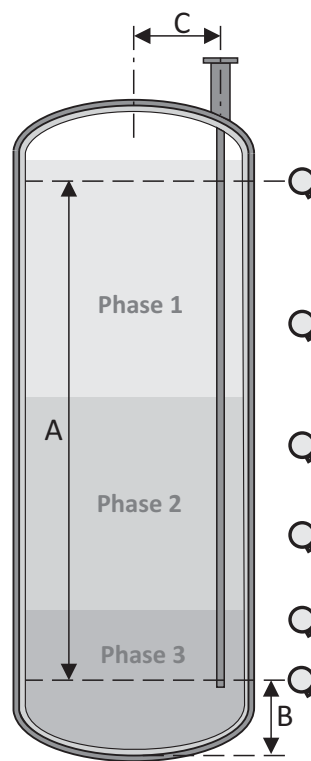
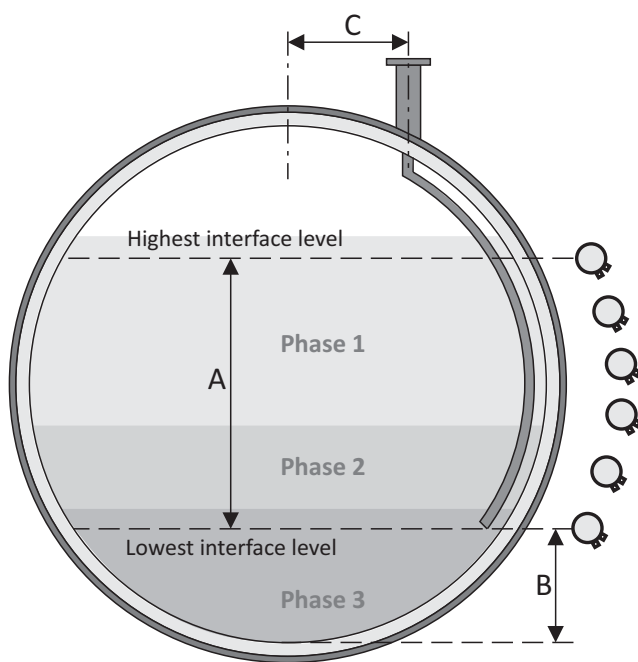
Insulation thickness ☐ mm ☐ inch

Insulation material

Additional cladding/heating jacket (thickness & material)

Measurement range (A) ☐ mm ☐ inch

Offset (B) ☐ mm ☐ inch



Installation Details (please attach drawing)

Dip pipe ☐ new/design phase ☐ existing

Thickness

Nozzle

Tag Size Pressure Diameter Thickness

Position - distance from center line (C):

☐ Obstruction i.e. electrical grids, heating (please attach drawing)

Measured Material Specifications

		Density range		Level range (from vessel bottom)	
		Unit g/cm ³ (if other, please specify)		☐ mm ☐ inch	
		min.	max.	min.	max.
top	Phase 1 Name				
	Phase 2 Name				
	Phase 3 Name				
	Phase 4 Name				
bottom	Phase 5 Name				

Emulsion layer ☐ no ☐ yes Expected layer thickness ☐ mm ☐ inch

Build-ups on vessel wall ☐ no ☐ yes Expected thickness ☐ mm ☐ inch

Process Conditions

	Unit (if other, please specify)	normal	min.	max.
Process temperature	°C			
Ambient temperature	°C			
Process pressure	bar			

Output Value

- ☐ Density: Number of measuring points (resolution)
- ☐ Density & Level of each phase

Detectors

Power supply ☐ 100-240 VAC ☐ 24 VDC

Exproof requested ☐ No ☐ Yes Type

Functional safety ☐ none ☐ SIL 2 ☐ SIL 3

Housing material ☐ SS304 (standard) ☐ SS316L (e.g. offshore) ☐ others Type

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.