

Company		Contact Name	
Street/P.O.Box		E-Mail	
City		Phone	
Postal code		Date	
Country / State		Project	

Process Specifications

Measuring tag	
Application/Process	
Product	

	Unit (if other, please specify)	normal	min.	max.
Moisture (weight fraction)	%			
Bulk density	kg/m ³			
Particle size	mm			
Product operating temperature	°C			

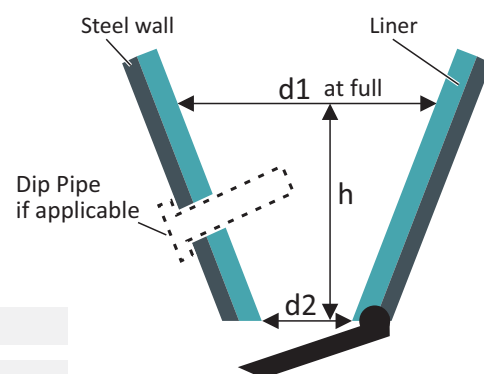
Crystal water ☐ No ☐ Yes If yes, how much?

Material composition is varying? ☐ No ☐ Yes

Accuracy % (absolute value in weight %)

Vessel Specification

	Dimensions ☐ mm ☐ inch	Material	Density g/cm ³
Inner distance d1 (at full)			
Inner distance d2 (at outlet)			
Product level h			
Steel wall thickness			
Liner wall thickness			
Others			



Vessel drawing and a vessel cross section drawing are mandatory!

Is it possible to install a dip pipe? ☐ No ☐ Yes

☐ Batch measurement

☐ Continuous product filling level

Instrumentation

	min.	max.	Unit (if other, please specify)
Ambient temperature at measuring point			°C
Power supply	☐ 90-250V AC/DC	☐ 24V AC/DC	