

## QUESTIONNAIRE

## Microwave Concentration & Moisture Measurement

Company Street/P.O.Box City Postal code Country / State				Contac E-Mail Phone Date Project	t Name			
Process Specifications								
	□ % dry substa	ance $\square$ %	water cont max.		Accuracy	,		Standard deviation
Consistence	□ liquid			☐ pasty		□ lun	npy	
Particle size	☐ refined	l		□ coarse		□ арј	orox. non-v	arying
Grain size distribution	%		mm	%	m	nm	%	mm
Abrasion	□ low	☐ medium	_					
(Bulk) density [g/cm³]	min.	nom.	max.	Electric	ally conduc	tive 🗌 Yes	n	nS/cm
Product temperature [°	[C]			Salt con	tont	☐ Yes		
Ambient temperature [	[°C]			Sait Coil	iterit	☐ No	V	Veight%
Installation								
☐ Conveyor Belt	Belt material		Ве	elt reinforced	11816611 —	Yes No		
Nom. dumping height (H) [mm]								
Min. dumping height (H) [mm]								
Max. dumping height (H		エ						
Typical dumping width (b) [mm]								
Belt width (B) [mm]	A							
Distance to return belt (A) [mm]								
Belt speed [m/s]	miı	n.	max.	со	nstant			
Belt occupancy	smooth	scrape	r available	☐ hei	ght sensor a	vailable		
☐ Pipeline								
Nominal width	□DN	□Inch			ection type	e: flange EN 1	1092-1/11	
Pressure range [bar] n	nin. no	om.	max.			IN 11853-1		
Flow	horizontal	☐ vertical		othe	r			
☐ Other Arrangement (sketch necessary!)								
Other								
Power supply	□100-240 V	AC 🗆 2	4 V DC					
Ex-proof needed								
Comments								