

Concentration & Moisture Measurement with Microwave Systems

Please fill out this questionnaire as completely as possible for project planning.

Project: _____ Date: _____
Company: _____ Name: _____
Street / P.O. Box: _____
Postal Code / City: _____ Country / State: _____
Phone: _____ Fax: _____
Email: _____

Product description _____
Chemical composition _____
Measured variable %DS Weight-% _____
Measuring range min.: _____ nom.: _____ max.: _____
Requested accuracy _____ Standard deviation _____
Consistence liquid pasty lumpy
Particle size refined coarse approx. non-varying
Grain size distribution % _____ mm % _____ mm % _____ mm
Abrasion low medium high
(Bulk)density [g/cm³] min.: _____ nom.: _____ max.: _____
Product temperature [°C] min.: _____ nom.: _____ max.: _____
Ambient temperature [°C] min.: _____ nom.: _____ max.: _____
Electrically conductive Yes _____ mS/cm No
Salt content constant _____ Weight.-%
Power supply 100-240 V AC 24 V DC
Ex-classification Yes Zone/Class: _____ No

Arrangement Pipeline

Nominal Width, Flange mm _____ Inch _____ Flange _____
Pressure range [bar] min.: _____ nom.: _____ max.: _____
Flow horizontal vertical

Arrangement Belt

A sketch with belt and material profile necessary!

Belt data Width: _____ Belt material: _____
Belt reinforced (steel) Yes No If yes, please provide sketch.
Belt speed [m/s] min.: _____ max.: _____ constant: _____
Dumping height [mm] min.: _____ nom.: _____ max.: _____
 smooth Scraper Leveling

Arrangement Other

A sketch of the arrangement is very helpful to clarify the project.

Appendix

Comments: _____

