



Lab Water Customer Request Form

Sales Rep.: _____ Date: _____

Customer: _____

Address: _____ Room: _____

Phone-Number: _____ Fax-Number: _____ E-mail: _____

Application(s):

- HPLC
- GC-MS, AAS, ICP-MS
- Nutrient media for cell culture
- Photometry
- ELISA, RIA
- Ion chromatography
- Immunocytochemistry
- TOC analysis
- Histology
- Production of monoclonal antibodies
- Microbiological media & reagents
- Solutions for chemical analysis & synthesis
- Electrophoresis
- Endotoxin analysis
- PCR
- Buffers and pH solutions
- Feed water for lab. devices humidifier | autoclave dishwasher

Feed Water source:

- RO
- EDI
- TAP
- DI
- Distilled

Other application

System should be hooked up to:

- Dishwasher
- Autoclave
- Humidifier
- Ultrapure water system
- Others: _____
- Should not be hooked up (stand-alone system)

System should be:

- New system:
- Replacement system

Currently used system: _____

Water Quantity

per hour: min. max. _____

per day: min. max. _____

per week: min. max. _____

Additional Information

Available Budget: _____

Expected delivery date: _____

Comments: _____

