

## Bottled at the Source



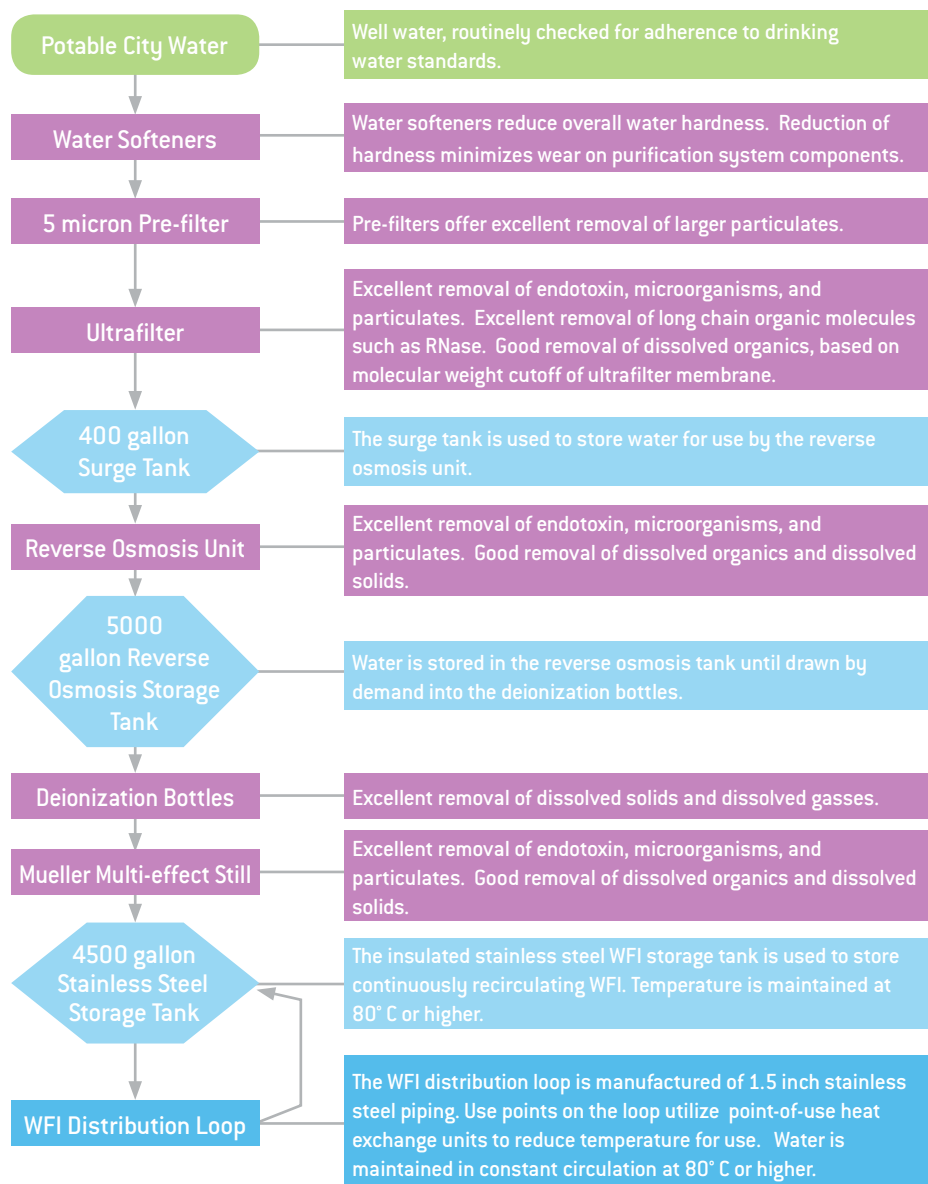
## WFI Quality Water for every need

At Lonza, we manufacture and test our Water For Injection (WFI) products to strict industry standards. Prepared through a process of ultrafiltration, reverse osmosis, deionization, distillation and sterile filtration, we meet or exceed USP and EP requirements.

Our off-the-shelf WFI products are available in 500 ml or 1 L bottles, 20 L carboys and 20 L, 50 L, 100 L and 200 L Platinum UltraPAK™ bags for your convenience. We can also provide made-to-order WFI in custom packaging options to fit your application.

When you are looking for high quality water for cell culture work, in start up mode with your bioprocessor, or for cleaning your core, we offer ready to use, cost effective WFI options.

## WFI Process Flow



## Contact Information

### North America

Customer Service: 800-638-8174  
 Technical Service: 800-521-0390  
 E-mail: [biotechserv@lonza.com](mailto:biotechserv@lonza.com)  
 Online Ordering: [www.lonza.com](http://www.lonza.com)

### Europe

Customer Service: 32 (0) 87 321 611  
 Technical Service: 32 (0) 87 321 651  
 (German Speaking)  
 32 (0) 87 321 626  
 (French Speaking)  
 E-mail: [techsup.europe@lonza.com](mailto:techsup.europe@lonza.com)  
 or [techsup.uk@lonza.com](mailto:techsup.uk@lonza.com)  
 Online Ordering: [www.lonza.com](http://www.lonza.com)

### International

Contact your local Lonza Distributor  
 Customer Service: 301-898-7025, ext. 2322  
 Fax: 301-845-8291  
 E-mail: [biotechserv@lonza.com](mailto:biotechserv@lonza.com)

### International Offices

Australia	61 3 9550 0883
Austria	0800 201 538
Belgium	32 (0) 87 321 611
Denmark	45 43 56 74 07
France	0800 91 19 81
Germany	0800 182 52 87
India	91 22 6697 2883
Italy	39 0363 35 14 70
The Netherlands	0800 022 4525
Spain	34 902 531 366
Switzerland	0800 83 86 20
United Kingdom	44 118 979 5234

## Ordering Information

Cat. No.	Description	Size
17-724F	Water for Cell Culture	500 ml bottle
17-724Q	Water for Cell Culture	1 L bottle
04-585P	Water for Cell Culture	20 L carboy
08-199F/02-038P20	Water for Injection*	20 L UltraPAK™
08-199P	Water for Injection*	50 L UltraPAK™
08-199D	Water for Injection*	100 L UltraPAK™
08-199E/02-038P200	Water for Injection*	200 L UltraPAK™

\*Bulk packaged for commercial use. Not for parenteral administration.

Lonza Walkersville, Inc.  
 Walkersville, MD 21793

For further manufacturing and laboratory use. Not for use in diagnostic procedures.

Unless otherwise noted, all trademarks herein are marks of the Lonza Group or its affiliates.

© Copyright 2007, Lonza Walkersville, Inc. All rights reserved.  
 FL-WFI-1 08/07

# Lonza