

## QC Data Management Solutions

Key Points		Est. \$\$ Pkg	
<b>UnityWeb™ 2.0 UW</b> <i>(Web-based)</i>	<ul style="list-style-type: none"> <li>• Quick and easy implementation (no IT involvement)</li> <li>• Includes Westgard Rules, Multi Test Data Entry, Bench Review, Actions/Comments, Qual Data</li> </ul>	<ul style="list-style-type: none"> <li>• Simple point data entry (includes Instant QC) or monthly summary data entry</li> <li>• Best for those who only want to submit data for peer review</li> </ul>	UW and WC 900-2100 yr 1 400-1400 yr 2+
<b>Unity Desktop™ 1.0 UDT</b> <i>(PC-based)</i>	<ul style="list-style-type: none"> <li>• Single laboratory PC (some IT involvement)</li> <li>• Basic features: Rules, Graphs, Reports</li> </ul>	<ul style="list-style-type: none"> <li>• Minimal system requirements (Win 98 or higher)</li> <li>• Westgard Advisor™ is not available</li> </ul>	UDT and SC 1600-2500 yr 1 700-1200 yr 2+
<b>Unity Real Time™ 1.0 URT</b> <i>(PC or Network)</i>	<ul style="list-style-type: none"> <li>• Significant IT involvement with extensive setup &amp; configuration</li> <li>• “Gold standard” for QC management: Bench and Supervisor Review, Audits, Data Analysis Grid, Qual Data, Advanced</li> </ul>	LJ Options, Automatic Updates and Data Submission <ul style="list-style-type: none"> <li>• Can replace instrument or LIS QC packages</li> <li>• Westgard Advisor™ (optional purchase)</li> </ul>	URT and 1 DC 3200-5800 yr 1 1500-2600 yr 2+ <i>additional costs possible</i>
<b>Unity Real Time™ Online 1.0 URTO</b> <i>(Web-based)</i>	<ul style="list-style-type: none"> <li>• Comprehensive QC management on the internet (minimal IT involvement)</li> <li>• Alternative to URT for labs with limited IT resources or software installation restrictions</li> </ul>	<ul style="list-style-type: none"> <li>• Updates and support handled by Bio-Rad</li> <li>• Westgard Advisor™ Online (optional purchase)</li> </ul>	URTO and WC 2700-4800 yr 1 1800-3500 yr 2+

All solutions allow easy participation in the Unity Interlaboratory Program and facilitate regulatory compliance.

## Connectivity Options

Key Points		
<b>WebConnect™ 2.0 WC</b>	<ul style="list-style-type: none"> <li>• Downloads and imports a QC file to the internet</li> <li>• Saves time/effort by eliminating manual entry of results and transcription errors</li> </ul>	<ul style="list-style-type: none"> <li>• Simplified Vitros Slide Generation handling capabilities</li> <li>• Web-based solution that bypasses IT issues</li> </ul>
<b>SoftConnect™ SC</b>	<ul style="list-style-type: none"> <li>• Imports QC file to PC/network software</li> <li>• Saves time/effort by eliminating manual entry of results and transcription errors</li> </ul>	<ul style="list-style-type: none"> <li>• Some IT involvement</li> </ul>
<b>DirectConnect™ 1.0 DC</b>	<ul style="list-style-type: none"> <li>• Only used with URT</li> <li>• Wired and wireless options available</li> <li>• Requires significant IT involvement</li> </ul>	<ul style="list-style-type: none"> <li>• Directs QC data directly from instruments to URT in real time</li> </ul>
<b>Data Innovations DI</b> <i>(instrument manager purchased from DI)</i>	<ul style="list-style-type: none"> <li>• Integrated or older driver methods available</li> <li>• DI partnerships include Roche, Abbott, Ortho, McKesson and Dairyland</li> </ul>	<ul style="list-style-type: none"> <li>• QC from any instrument interfaced via DI to URT in real time</li> <li>• License or driver must be purchased from DI or one of its partners</li> </ul>

### Legacy Solutions

- Unity Post™, Unity PC™ and Unity Plus/Pro™ should be upgraded to a current solution in 2008
- QC OnCall™ is no longer available for new installations but support will be continued
- UnityWeb™ 1.0 has been replaced by UnityWeb™ 2.0
- WebConnect™ 1.0 will be replaced by WebConnect™ 2.0



**Bio-Rad  
Laboratories**

*For more information visit us on the web at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics). Call toll free at 1-800-2BIO-RAD (1-800-224-6723); outside the US, contact your local sales office.*