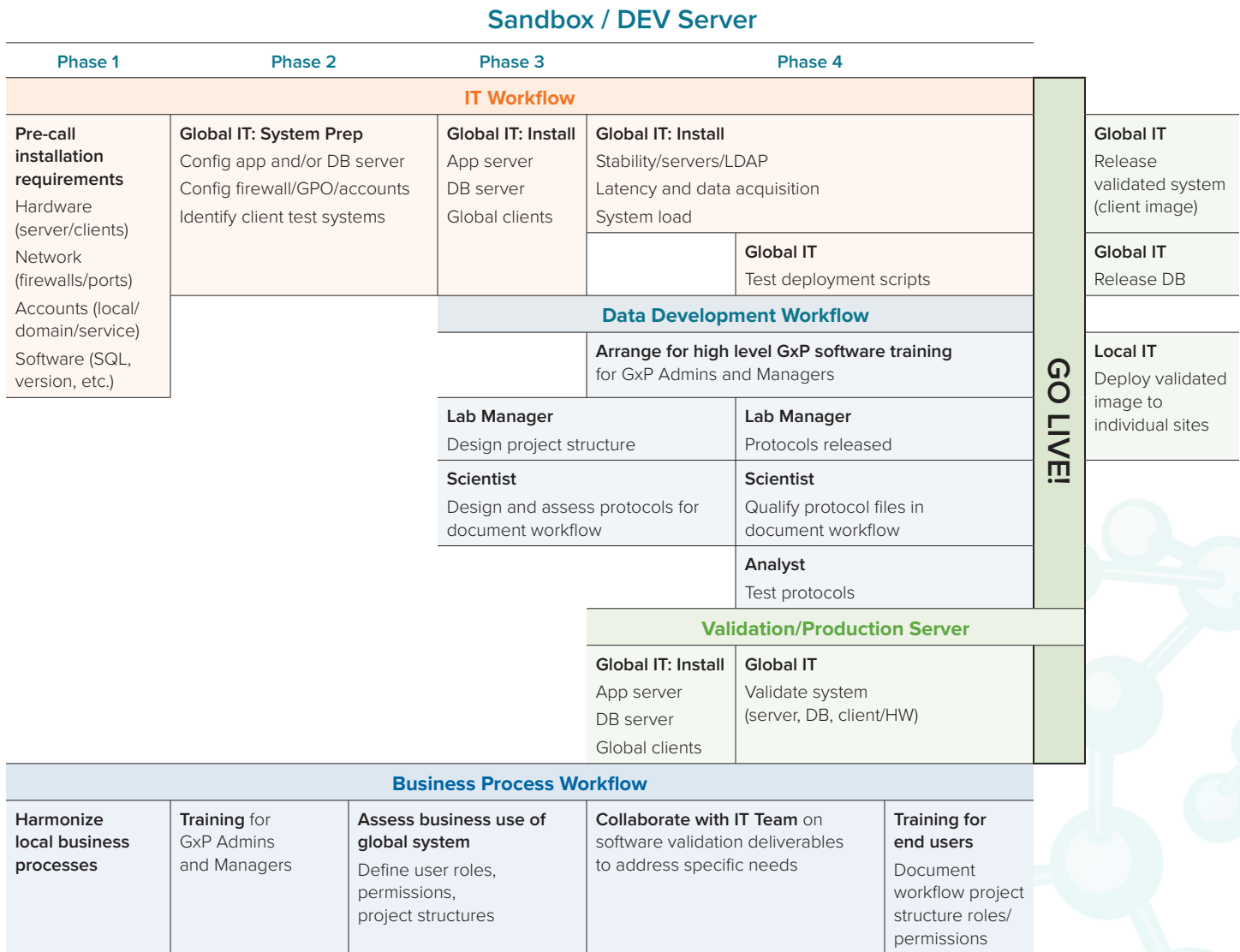


SoftMax Pro 7.1.2 GxP Software Enterprise Rollout

SoftMax® Pro GxP Software scales from a single computer installation to an enterprise solution with multiple client computers connected to a database hosted on a central server. For enterprise rollouts, our professional services team is experienced installing our software in a range of IT environments and can assist in planning and scaling your implementation according to your requirements.

Representation of a SoftMax Pro GxP Software enterprise rollout is shown below.



Contact Us

Phone: +1.800.635.5577
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom