

For a detailed protocol, see the *Water-Glo™ System Technical Manual #TM547*.

Single Test Protocol



ATP Calculation Worksheet



Calculate ATP concentration using the following formula:



Or download the Water-Glo[™] ATP Calculator Tool at: www.promega.com/resources/tools/water-glo-atp-calculator-tool

Promega Quick Protocol

96-Well Batch Protocol



3. Calculate ATP Concentration

To calculate ATP concentration, copy exported data into the Water-Glo[™] ATP Calculator Tool, available at: www.promega.com/resources/tools/water-glo-atp-calculator-tool

Additional protocol information is in Technical Manual #TM547. For detailed instructions on instrument use, refer to the GloMax[®] System Operating Manuals #TM397, #TM433 or #TM470, available online at: **www.promega.com**