



I. Description

The Plexor™ qPCR and qRT-PCR Systems^(a,b) are compatible with a variety of real-time PCR instruments. Data from these instruments can be analyzed with one dedicated software program, the Plexor™ Analysis Software. This manual includes instructions and thermal cycling conditions specific for use of the Plexor™ qPCR System, Plexor™ Two-Step qRT-PCR System and Plexor™ One-Step qRT-PCR System with the Applied Biosystems 7900HT Real-Time PCR System. Instructions are included for instrument setup, data transfer from the instrument to the Plexor™ Analysis Software and data analysis.

II. Plate Preparation and Amplification

For detailed instructions describing assay setup, see the *Plexor™ qPCR System Technical Manual #TM262*, the *Plexor™ qRT-PCR One-Step System Technical Manual #TM263* and *Plexor™ qRT-PCR Two-Step System Technical Manual #TM264*. Additional information is available at www.promega.com/plexorresources/

When using the Plexor™ qPCR or qRT-PCR System for the first time, we recommend programming the thermal cycling conditions and checking that the instrument is compatible with the dyes used and is configured for those dyes before assembling the reactions, so that the reactions are not kept on ice for prolonged periods of time. Once you are familiar with the programming process, the instrument can be programmed after assembling the reactions.

Materials to Be Supplied By the User

- centrifuge capable of centrifuging a 96- or 384-well plate
 - optical adhesive covers and applicator
 - foam compression pad for 96-well plate
1. After the amplification reactions have been assembled, cover the reaction plate with an optical adhesive cover using the applicator. Centrifuge briefly to collect contents at the bottom of each well.
Note: Keep the plate on ice during reaction setup and programming of the thermal cycling conditions.
 2. Program the Applied Biosystems 7900HT Fast Real-Time PCR System. The proper thermal cycling conditions and instructions for programming the instrument are provided in Section III (qPCR and two-step qRT-PCR assays), Section IV (one-step qRT-PCR assays) and Section V (genotyping assays).