

1311 Orleans Drive  
Sunnyvale, CA  
94089-1136  
Ph +1-800-635-5577  
Fax +1-408-548-6461

## GxP Admin Software Version 2.1

### Primary Administrator password change

The changing of the primary GxP administrator's password by a secondary administrator was not recognized by the GxP Admin 2.0 application. The updated password would have to be re-verified and changed through the SoftMax Pro GxP application..

**Fix** – An update to the GxP Admin component (v2.1) addresses this issue.

**Note:** GxP Admin 2.1 may be used with SoftMax Pro GxP v5.2. It is recommended for users upgrading to copy the mdclisen.ini and mdcadmgr.ini files from the GxP Admin 2.0 program folder to that of GxP Admin 2.1 to preserve license codes and user permission templates.

## GxP Admin Software Version 2.1.1 Software Release Notes

Introduction The GxP Admin Software version 2.1.1 is a minor update release. The following are the change incorporated in this update as compared to version 2.1, the last general release of the GxP Admin Software.

### New in GxP Admin v2.1.1

**Software Rebranding** The software splash screen and artworks are updated to reflect Molecular Devices Inc. All references to MDS Analytical Technologies are removed.

**Impact:** This fix has no impact on current workflow or data.

**Improving the path to the User Account File** In order to enhance the usability, the user interface has been updated to reflect the full path to User Account file. Previously only the file name was visible with no indications to the path.

**Impact:** This fix has no impact on current workflow or data.