

April 2015

# Important Note

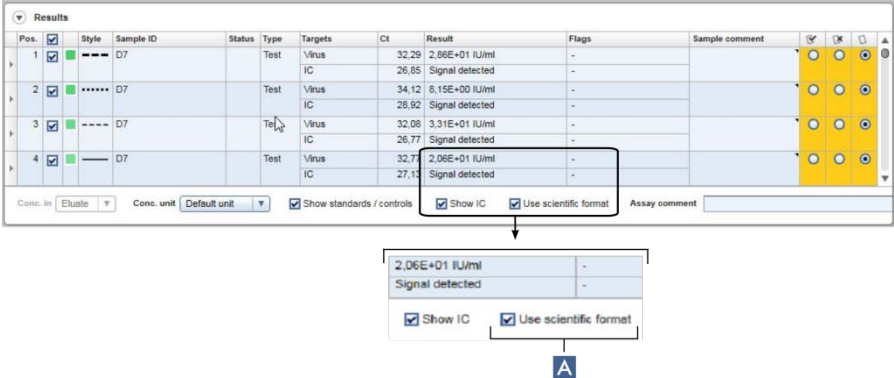
Dear Rotor-Gene AssayManager user,

We would like to inform you about new functionalities provided by **Rotor-Gene AssayManager® artus® Basic Plug-in software version 1.0.4**. These improvements are not yet described in the *Rotor-Gene AssayManager artus Basic Plug-in User Manual*. Therefore please read this Important Note carefully.

The improvements relate to Section 1.3.1.3 (*artus* Specific Tasks and Procedures → Approving Samples → Concept of Approval Buttons in *artus* Plug-in) in the *Rotor-Gene AssayManager artus Basic Plug-in User Manual*. The following changes have been made, described in greater detail below:

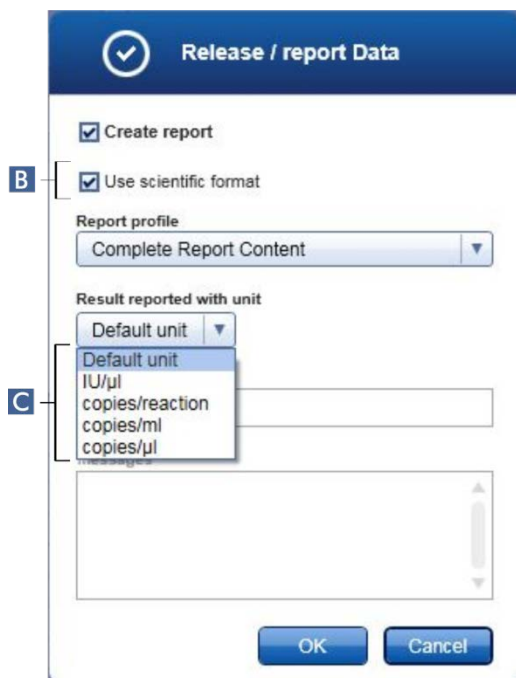
- Improvement to the “Results” area in the “Approval” environment
- Improvements to the “Release/ report Data” dialog
- Improvements to the “Test Results – Overview” report
- New error message

- **Improvement to the “Results” area in the “Approval” environment.** Option to display results in scientific format [A] has been added (Figure 1).



**Figure 1. Additional functionality in the “Results” area in the “Approval” environment.** [A] Additional option to display results in scientific format.

- **Improvements to the “Release/ report Data” dialog.** Options to display results in scientific format [B] and to choose the concentration unit in the report overview table [C] have been added (Figure 2).



**Figure 2. Improvements to the “Release/ report Data” dialog.** [B] Additional option to display results in the report in scientific format. [C] Additional option to select the concentration unit in the report overview table.

- **Improvements to the “Test Results – Overview” report.** A column that displays the approval status for each sample and control has been added [D], there are improved options for displaying results in terms of concentration unit and scientific format [E], and a column that indicates if flags are assigned to a target has been added [F] (Figure 3).

D

E

F

Id	Color	Approval	Target	Ct	Result	Flags
D7	<div></div>	✓	Virus	32,77	2,06E+01 IU/ml	
			IC	27,13	Signal detected	

All concentrations given in this table are concentrations in the eluate

! This target has flags

✓ Accepted

x Rejected

**Figure 3: Improvements to the “Test Results – Overview” report.** [D] New column displaying the approval status for each sample and control. [E] Improved options for displaying results in terms of concentration unit and scientific format. [F] New column indicating if flags are assigned to a target.

**Note on scientific format:** To display quantitative results, Rotor-Gene AssayManager *artus* Basic Plug-in software version 1.0.4 lets the user choose between scientific format and decimal format in the “Approval” environment and in the report. The approval screen contains a check box “Use scientific format” in the results area below the results table [A]. If the check box is activated, the concentrations in the result column [E] of the results report are displayed in scientific format (e.g., 222,732.63 IU/ml would be displayed as 2.23E+05 IU/ml).

You can also choose to display results in scientific format in the report. The “Release/ report Data” dialog contains an additional check box “Use scientific format” [B]. If that check box is activated, all concentrations in the report are displayed in scientific format [E].

**Note on configurable concentration unit in report overview table:** Rotor-Gene AssayManager *artus* Basic Plug-in software version 1.0.4 lets you select the concentration unit to be

displayed in the report. This functionality is provided if the assay is quantitative and if there is more than one concentration unit defined in the assay profile. The “Release/ report Data” dialog contains a dropdown box [C] from which you can choose the concentration unit for the selected assay. The selected unit will be used for concentrations in the overview table in the report [E].

**Note on approval status and flags information in reports:** Reports created with Rotor-Gene AssayManager *artus* Basic Plug-in software version 1.0.4 contain two additional columns in the overview table: a column that displays the approval status of each control or test sample (either accepted or rejected) [D], and a column that displays if there are any flags assigned to a target (indicated by the symbol “!”) [F].

- **New error message.** One new error message is now available in Rotor-Gene AssayManager *artus* Basic Plug-in software version 1.0.4, listed in the table below.

Error ID	Error Text
550272	The generated report could not be opened. Verify that you have installed a pdf viewer on your system.

If you have any questions, please contact your local QIAGEN Technical Services Department.

Best regards,

Your QIAGEN Rotor-Gene AssayManager Team

Trademarks: QIAGEN®, *artus*®, Rotor-Gene AssayManager® (QIAGEN Group).  
1094900 04/2015 © 2015 QIAGEN, all rights reserved.