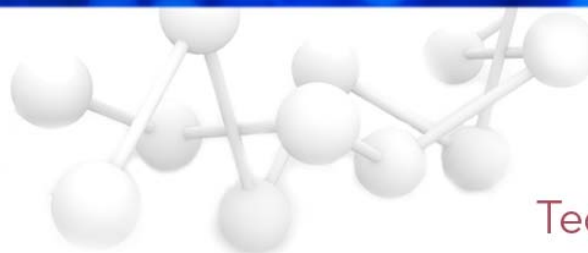


Errata Notice

This document contains references to BioTek.
Please note that BioTek is now Agilent.
For more information, go to
www.agilent.com/lifesciences/biotek



Gen5™ File Export

Creating a File to Export to LIMs/LIS

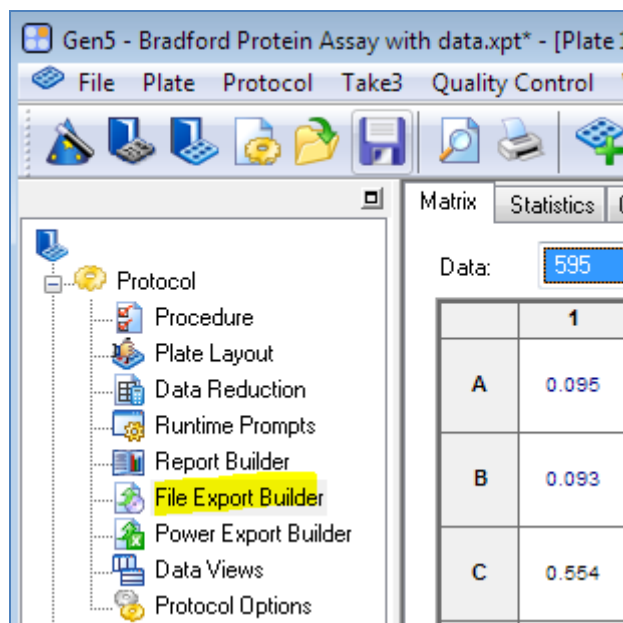
Introduction

Many commercially available LIMs/LIS applications can easily import text files (*.txt) from instrument control and analysis software programs like Gen5. This Tech Note describes the steps required to create a custom *.txt containing the required data, using the File Export Builder feature in the Gen5 protocol. The File Export Builder allows a user to create a template containing only the data that needs to be exported in the file, in either a column (table) or matrix (grid) format. See the example screens in the Appendix.

This Tech Note also describes the Gen5 protocol options available to define the exported file path, manual or automatic exporting and information about exporting each plate's data to its own file or appending multiple plates to a single file.

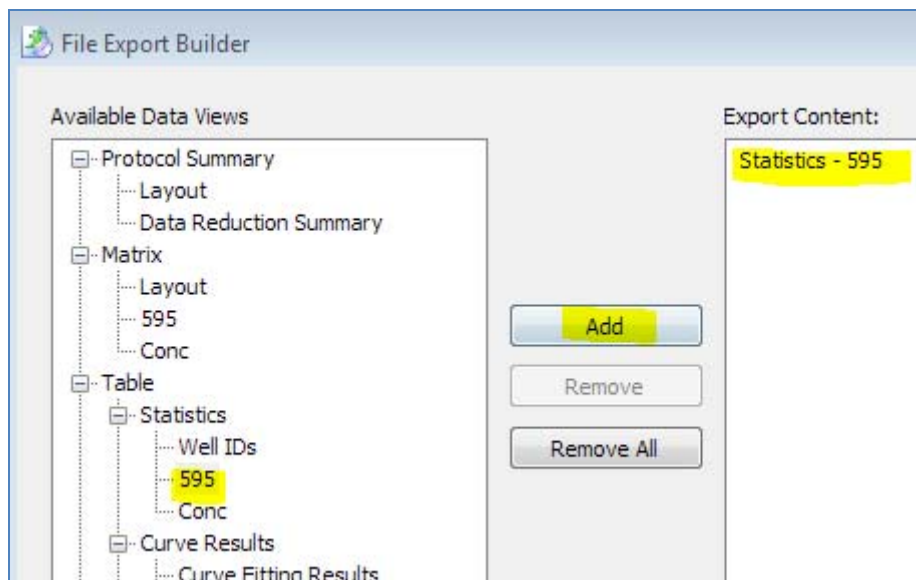
Step 1: Select the data to be exported using the **File Export Builder**

- Open the Gen5 protocol file
- Double-click the File Export Builder in the protocol menu:

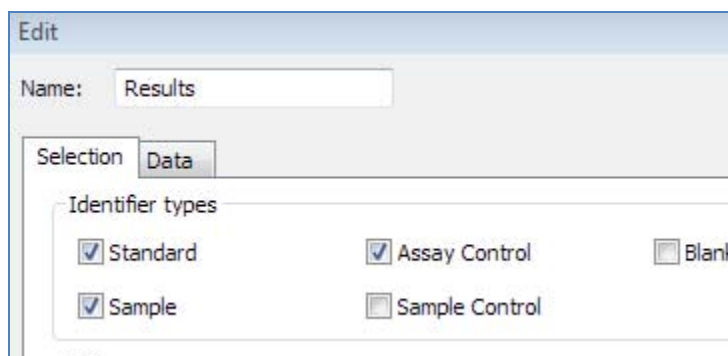


- c) In the File Export Builder, select data to be included in the export file from the **Available Data Views** list. “Matrix” data views will export data in a microplate matrix or grid format. “Table” views will export data in a column format. Table format is more commonly used for data that will be sent to a LIMS/LIS.

In the screen below, we’ve selected 595 nm data, then clicked Add to add this to the **Export Content**:



- d) Double-click on the item in the Export Content window to edit the contents. In this dialog, you may:
1. Change the name of the contents
 2. Choose the well types to be included in the file (in the Selection tab, check all desired types)



3. In the Data tab, select the type of data to be included. Several pieces of data are included by default. Any can be removed by right-mouse clicking and choosing Delete. Others may be added by clicking in an open row under the Data header:

Edit

Name: Results

Selection Data

	Data	Stat. Op.	Title
1	Well ID		<Well ID
2	Name		<Name
3	Well		<Well
4	595	Repl.	<595
5	Conc	Repl.	<Conc
6			

Step 2: Define the File Export Settings

- a) The **Protocol Options|File Export Settings** screen is shown below. In this screen, several default selections are made, but all can be modified.

Protocol Options

Protocol Type
Calculation Options
Report Options
Export Options
File Export Settings
Power Export Settings
Quick Export Settings
Clipboard Settings
Save Options
Auto-Output Options

File Export Settings

File Naming Convention

File Name: Export Extension: txt

Example: Export.txt

File Location

Last folder used
 Folder: S:\TempOCT

Example: S:\TempOCT

When Exporting, prompt before saving file:

Always
 Only if the file already exists
 Never, if the file exists...
 Append Overwrite

Include:

Headings
 Matrix column & row labels
 Statistic column labels

Separator

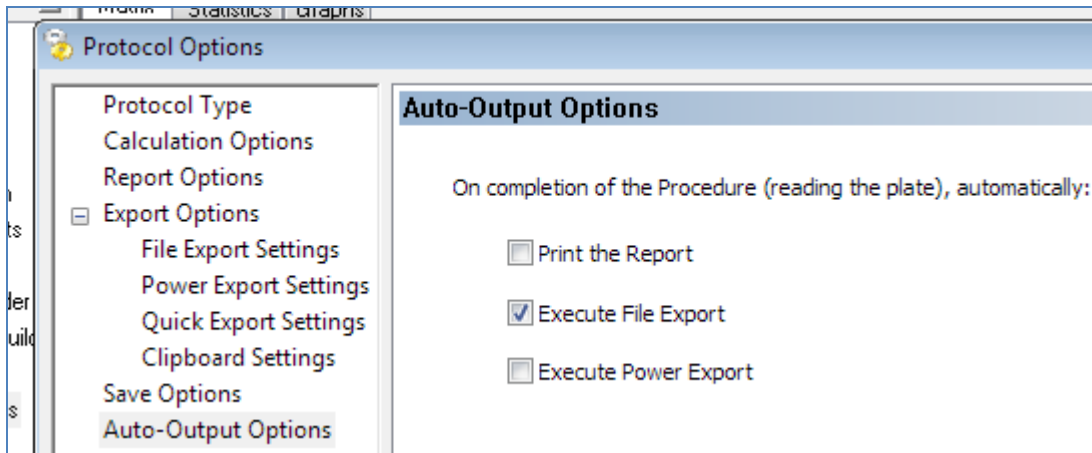
TAB
 ;
 ,
 Other:

1. The File Naming Convention should use the Extension "txt". By default, the Experiment name is selected as the File Name. To change this, click the right arrow at the end of the File Name field, and choose the desired File Name parameter from the list.
 2. Select the appropriate path for the File Location
 3. It is best to de-select all of the "Include" options; Separator should be selected by the LIMs/LIS requirement.
 4. In the "When Exporting ..." box, select the desired prompt. *Note that if you wish to send multiple plates to a single Export file, you must choose the "Never, ..." and Append option.*
- b) To Export the data, highlight the **Plate** containing the data to be exported - right-mouse click to see the menu. Choose "File Export". The data will be sent to the path specified in Step 2 above.

Step 3: Export the Data

Data can be exported from Gen5 either manually or automatically after each plate is read.

- a) To Export the data automatically after each plate is read, go to Protocol|Auto Output Options and select Execute File Export:

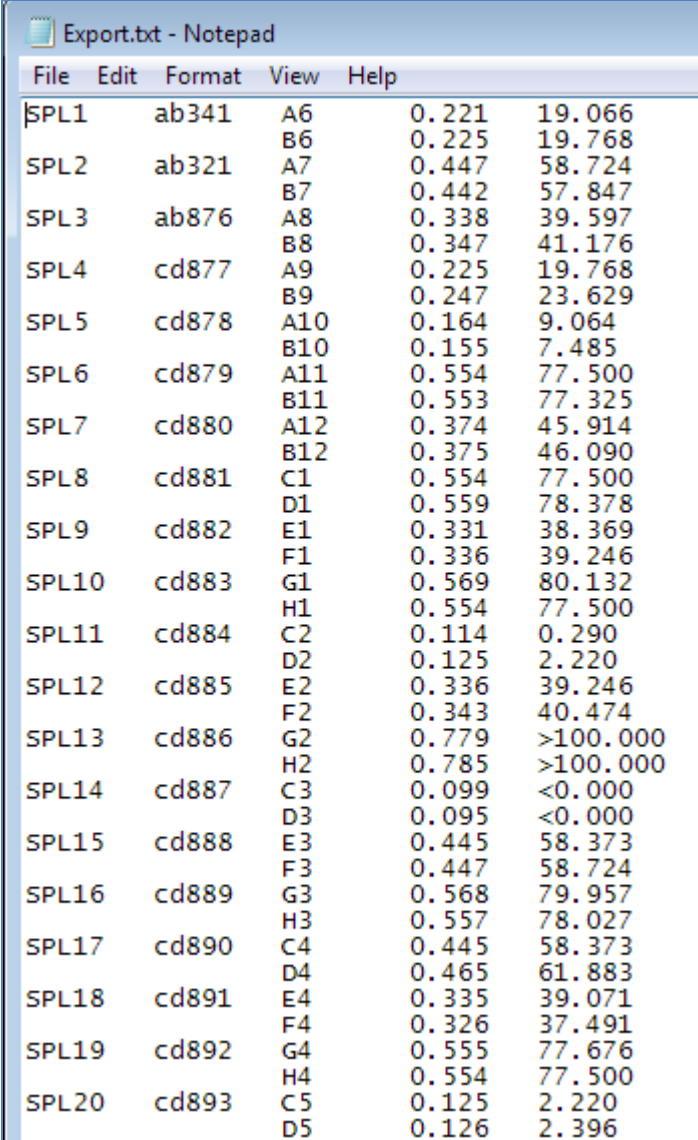


Each time a plate is read in the experiment file, the data will be automatically exported to the path defined in Step 2.

- b) To Export data manually from Gen5 to a *.txt file, first Read the plate. When reading is complete, right mouse click on the Plate in the menu tree and choose File Export. The data will be exported to the path defined in Step 2. If more than one plate exists in the Experiment, all plates can be highlighted and then exported.

APPENDIX

Example of data exported in the **Table** format:



The image shows a Notepad window titled "Export.txt - Notepad" with a menu bar containing "File", "Edit", "Format", "View", and "Help". The main text area contains a table with 20 rows of data. Each row represents a sample (SPL1 to SPL20) with associated codes and numerical values.

SPL1	ab341	A6	0.221	19.066
		B6	0.225	19.768
SPL2	ab321	A7	0.447	58.724
		B7	0.442	57.847
SPL3	ab876	A8	0.338	39.597
		B8	0.347	41.176
SPL4	cd877	A9	0.225	19.768
		B9	0.247	23.629
SPL5	cd878	A10	0.164	9.064
		B10	0.155	7.485
SPL6	cd879	A11	0.554	77.500
		B11	0.553	77.325
SPL7	cd880	A12	0.374	45.914
		B12	0.375	46.090
SPL8	cd881	C1	0.554	77.500
		D1	0.559	78.378
SPL9	cd882	E1	0.331	38.369
		F1	0.336	39.246
SPL10	cd883	G1	0.569	80.132
		H1	0.554	77.500
SPL11	cd884	C2	0.114	0.290
		D2	0.125	2.220
SPL12	cd885	E2	0.336	39.246
		F2	0.343	40.474
SPL13	cd886	G2	0.779	>100.000
		H2	0.785	>100.000
SPL14	cd887	C3	0.099	<0.000
		D3	0.095	<0.000
SPL15	cd888	E3	0.445	58.373
		F3	0.447	58.724
SPL16	cd889	G3	0.568	79.957
		H3	0.557	78.027
SPL17	cd890	C4	0.445	58.373
		D4	0.465	61.883
SPL18	cd891	E4	0.335	39.071
		F4	0.326	37.491
SPL19	cd892	G4	0.555	77.676
		H4	0.554	77.500
SPL20	cd893	C5	0.125	2.220
		D5	0.126	2.396

Example of data exported in the **Matrix** format:

Export2.txt - Notepad											
File	Edit	Format	View	Help							
STD1	STD2	STD3	STD4	STD5	SPL1	SPL2	SPL3	SPL4	SPL5	SPL6	SPL7
0.095	0.173	0.271	0.405	0.675	ab341	ab321	ab876	cd877	cd878	cd879	cd880
0.095	0.173	0.271	0.405	0.675	0.221	0.447	0.338	0.225	0.164	0.554	0.374
STD1	STD2	STD3	STD4	STD5	SPL1	SPL2	SPL3	SPL4	SPL5	SPL6	SPL7
0.093	0.167	0.269	0.415	0.669	ab341	ab321	ab876	cd877	cd878	cd879	cd880
0.093	0.167	0.269	0.415	0.669	0.225	0.442	0.347	0.247	0.155	0.553	0.375
SPL8	SPL11	SPL14	SPL17	SPL20	SPL23	SPL26	SPL29	SPL32	SPL35	SPL38	SPL41
cd881	cd884	cd887	cd890	cd893	cd896	cd899	cd902	cd905	cd908	cd911	cd914
0.554	0.114	0.099	0.445	0.125	0.442	0.664	0.442	0.354	0.114	0.339	0.487
0.554	0.114	0.099	0.445	0.125	0.442	0.664	0.442	0.354	0.114	0.339	0.487
SPL8	SPL11	SPL14	SPL17	SPL20	SPL23	SPL26	SPL29	SPL32	SPL35	SPL38	SPL41
cd881	cd884	cd887	cd890	cd893	cd896	cd899	cd902	cd905	cd908	cd911	cd914
0.559	0.125	0.095	0.465	0.126	0.457	0.655	0.451	0.345	0.116	0.334	0.468
0.559	0.125	0.095	0.465	0.126	0.457	0.655	0.451	0.345	0.116	0.334	0.468
SPL9	SPL12	SPL15	SPL18	SPL21	SPL24	SPL27	SPL30	SPL33	SPL36	SPL39	SPL42
cd882	cd885	cd888	cd891	cd894	cd897	cd900	cd903	cd906	cd909	cd912	cd915
0.331	0.336	0.445	0.335	0.198	0.337	0.472	0.231	0.332	0.098	0.246	0.214
0.331	0.336	0.445	0.335	0.198	0.337	0.472	0.231	0.332	0.098	0.246	0.214
SPL9	SPL12	SPL15	SPL18	SPL21	SPL24	SPL27	SPL30	SPL33	SPL36	SPL39	SPL42
cd882	cd885	cd888	cd891	cd894	cd897	cd900	cd903	cd906	cd909	cd912	cd915
0.336	0.343	0.447	0.326	0.187	0.329	0.493	0.216	0.390	0.102	0.258	0.199
0.336	0.343	0.447	0.326	0.187	0.329	0.493	0.216	0.390	0.102	0.258	0.199
SPL10	SPL13	SPL16	SPL19	SPL22	SPL25	SPL28	SPL31	SPL34	SPL37	SPL40	SPL43
cd883	cd886	cd889	cd892	cd895	cd898	cd901	cd904	cd907	cd910	cd913	cd916
0.569	0.779	0.568	0.555	0.201	0.998	0.553	0.222	0.559	0.241	0.557	0.661
0.569	0.779	0.568	0.555	0.201	0.998	0.553	0.222	0.559	0.241	0.557	0.661
SPL10	SPL13	SPL16	SPL19	SPL22	SPL25	SPL28	SPL31	SPL34	SPL37	SPL40	SPL43
cd883	cd886	cd889	cd892	cd895	cd898	cd901	cd904	cd907	cd910	cd913	cd916
0.554	0.785	0.557	0.554	0.195	1.025	0.549	0.221	0.567	0.251	0.568	0.641
0.554	0.785	0.557	0.554	0.195	1.025	0.549	0.221	0.567	0.251	0.568	0.641

TN102110_06

DE44335.0151157407

5994-3431EN
August 1, 2021