

Council on Dairy Cattle Breeding (CDCB) Export Tool

Pub. No. MAN1002001 Rev. A

The CDCB Export Tool is a companion application to Axiom™ Analysis Suite. It formats the genotype export and sample sheet to enable direct uploading to the Council on Dairy Cattle Breeding website.

Install the Council on Dairy Cattle Breeding (CDCB) Export Tool

IMPORTANT! Axiom™ Analysis Suite must be installed on your computer *before* installing the CDCB Export Tool.

1. After receiving/downloading the CDCB Export Tool zip package, save it to a local folder on your system.
2. Unzip the file, then double-click **CDCBExportToolSetup.exe** to install.
3. Follow the installer instructions.

Run the CDCB Export Tool

IMPORTANT! You need to have a previously run analysis in Axiom™ Analysis Suite to import into the CDCB Export Tool. For instructions about creating and running an analysis, see the *Axiom™ Analysis Suite User Guide* (Pub. No. [MAN0027928](#)).

1. In the Windows™ taskbar **Search** field, enter **CDCB**, then open the **CDCB Export Tool** from the results.

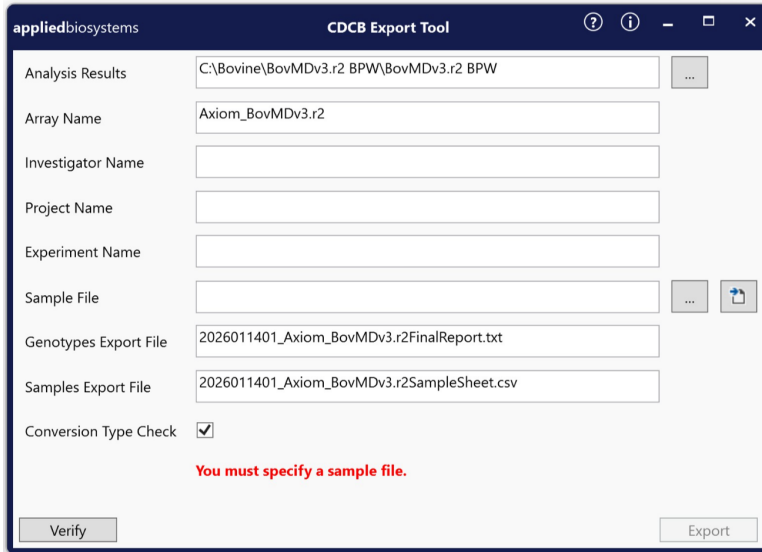
The screenshot shows the CDCB Export Tool application window. The title bar reads 'appliedbiosystems CDCB Export Tool'. The window contains the following elements:

- Analysis Results:** A text input field with a folder selection button (three dots) to its right.
- Array Name:** A text input field.
- Investigator Name:** A text input field.
- Project Name:** A text input field.
- Experiment Name:** A text input field.
- Sample File:** A text input field with folder selection and refresh buttons to its right.
- Genotypes Export File:** A text input field.
- Samples Export File:** A text input field.
- Conversion Type Check:** A checked checkbox.
- Error Message:** A red text message: "You must specify an analysis results folder."
- Buttons:** "Verify" and "Export" buttons at the bottom.

2. Click **...** (**open**) to the right of the **Analysis Results** field.

3. Navigate to your Analysis Results folder, click to highlight it, then click **Select Folder**.

The **Analysis Results**, **Array Name**, **Genotypes Export File**, and **Samples Export File** fields populate.



4. Optional: If needed, change the auto-generated **Genotypes Export File** and **Samples Export File** file names.
 - a. Click the text field.
 - b. Edit the file name, then click outside the text field to save.
5. If a Sample File already exists, go to step 7.

Note: The Sample File must be a tab-delimited TXT file with 8 columns.

6. Click  (**Create Sample File Template**) to the right of the **Sample File** field.

An 8-column template file is generated in the analysis folder with the CEL files, plate barcode, and well position columns automatically populated based on the selected analysis.

- a. Navigate to the Analysis Results folder, find the **CDCBExportSampleFileTemplate.txt** file, then open it using an Excel™ spreadsheet or a text editor.

	A	B	C	D	E	F	G	H
1	cel_files	sample_id	sample_name	sample_source	plate_barcode	well_position	sentrix_barcode	sentrix_position_a
2	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
3	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
4	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
5	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
6	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
7	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
8	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
9	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
10	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
11	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
12	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
13	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
14	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
15	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
16	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
17	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
18	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
19	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
20	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
21	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>
22	a550627-4	<sample id>	<sample name>	<sample source>	<plate barcode>	<well position>	<sentrix barcode	<sentrix position a>

Figure 1 CDCBExportSampleFileTemplate.txt example.

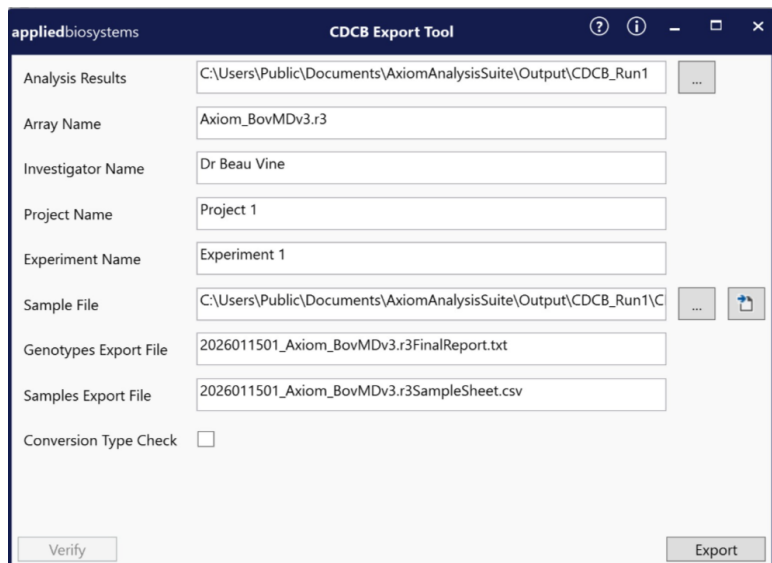
b. Complete the remaining/required fields.

Each column heading must use the same naming convention and be in the same left to right order as found in the template file.

c. Rename the template, then save it as a tab-delimited TXT file.

7. Click **...** (**open**) to the right of the **Sample File** field, navigate to the sample file, then click **Open**.

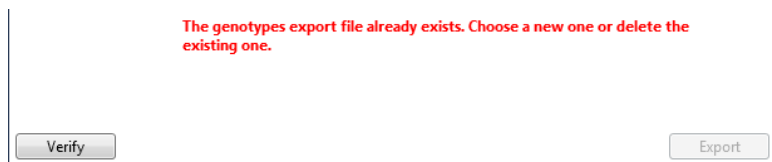
8. Optional: Enter an **Investigator Name**, **Project Name**, and **Experiment Name**.



The screenshot shows the 'CDCB Export Tool' window with the following fields and controls:

- Analysis Results:** C:\Users\Public\Documents\AxiomAnalysisSuite\Output\CDCB_Run1 (with a browse button)
- Array Name:** Axiom_BovMDv3.r3
- Investigator Name:** Dr Beau Vine
- Project Name:** Project 1
- Experiment Name:** Experiment 1
- Sample File:** C:\Users\Public\Documents\AxiomAnalysisSuite\Output\CDCB_Run1\C (with browse and refresh buttons)
- Genotypes Export File:** 2026011501_Axiom_BovMDv3.r3FinalReport.txt
- Samples Export File:** 2026011501_Axiom_BovMDv3.r3SampleSheet.csv
- Conversion Type Check:**
- Buttons:** Verify and Export

At any point, the software may detect an error or a duplicate file problem. If so, an error message appears.



The error message dialog box contains the following text:

The genotypes export file already exists. Choose a new one or delete the existing one.

Buttons: Verify and Export

a. Correct the reported problem, then click **Verify** to recheck/clear the displayed message.

9. Ensure that the **Conversion Type Check** option is *unchecked*.

Selecting this option converts all probesets classified as “Non Recommended” to 100% NoCall.

IMPORTANT! It is recommended to deselect this option as enabling it can significantly decrease call rates.

10. Click **Export**.

A progress bar appears, then after a few moments an **Export Complete** message appears.

11. Click the **X** button in the upper right corner to close the **CDCB Export Tool** window.

The generated Final Report (.txt) (genotype idata) and Samples Sheet (.csv) files reside in your Analysis Results folder and are ready for use.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Affymetrix, Inc. | 6055 Sunol Blvd | Pleasanton, California 94566 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN1002001 A

Revision	Date	Description
A	6 May 2026	Initial release.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF THE DOCUMENT OR THE PRODUCT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Microsoft, Windows, and Excel are either registered trademarks or trademarks of Microsoft Corporation in the United States and/or other countries.