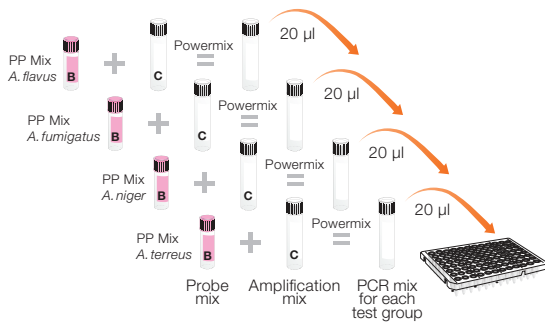




iQ-Design Assay

Aspergillus Identification Solution

Quick Guide



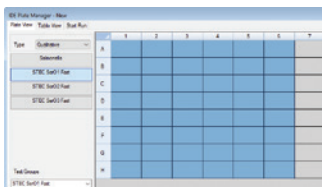
- Prepare the PCR mix for each Test Group: *A. flavus*, *A. fumigatus*, *A. niger*, and *A. terreus* (if applicable to local requirements) (see PCR Mix Calculation Guide on back)
- Use primer-probe vial (PP Mix) *A. flavus* for Test Group **iQ-DA *Aspergillus flavus***, PP Mix *A. fumigatus* for Test Group **iQ-DA *Aspergillus fumigatus***, PP Mix *A. niger* for Test Group **iQ-DA *Aspergillus niger***, and PP Mix *A. terreus* for Test Group **iQ-DA *Aspergillus terreus***
- Distribute 20 µl/well of each PCR mix in a microplate



- Add 5 µl each of positive and negative controls to designated wells
- Transfer 5 µl/well of extracted DNA (from iQ-Check™ *Aspergillus* spp. DNA extraction preparation) to the PCR microplate for each Test Group

Do not vortex before collecting the sample.
Check that there are no bubbles.

- Seal the PCR microplate

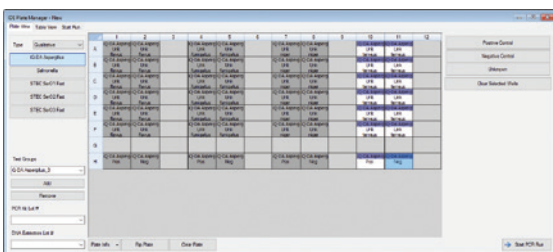


- Launch CFX Manager Industrial Diagnostic Edition (IDE) Software to get the Plate Manager window
- Select **iQ-DA *Aspergillus*** from the list of tests on the left side of the screen
- Select the required wells for the Test Group **iQ-DA *A. flavus*** (labeled as **iQ-DA *Aspergillus*** in the software)



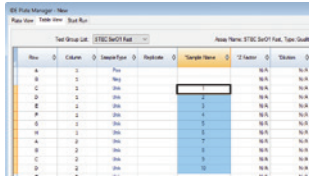
- Apply sample types to the wells by selecting **Positive Control**, **Negative Control**, or **Unknown**
- A separate positive control and negative control must be set up for each Test Group.**

- Enter the test group name (*flavus*, *fumigatus*, *niger*, or *terreus*), and sample name



- Select **Add** under the Test Group tab on the left-hand column of the window to repeat for Test Group **iQ-DA *A. fumigatus*** (labeled as **iQ-DA *Aspergillus_1***), Test Group **iQ-DA *A. niger*** (labeled as **iQ-DA *Aspergillus_2***), and Test Group **iQ-DA *A. terreus*** (labeled as **iQ-DA *Aspergillus_3***)

For detailed instructions, review the kit user guide.



- To duplicate the sample names for all Test Groups, select the Table View tab
- With **iQ-DA *Aspergillus*** selected in the Test Group dropdown, select the Sample Name column and then copy (Ctrl + C)
- In the Test Group List dropdown, move to **iQ-DA *Aspergillus*_1**, then select the Sample Name column and paste (Ctrl + V)
- Repeat for **iQ-DA *Aspergillus*_2** and **iQ-DA *Aspergillus*_3**
- Start the amplification by clicking **Start PCR Run**

PCR Mix Calculation Guide for the iQ-Design Assays

To find the correct volumes to use when preparing the PCR mix, add the total number of samples and controls to be analyzed and find the corresponding volumes of iQ-Design Assay Primer-Probe Mix and iQ Multiplex Powermix in the table.

| Total Number of Samples and Controls | iQ-Design Assay Primer-Probe Mix, µl | iQ Multiplex Powermix Amplification Mix, µl |
|--------------------------------------|--------------------------------------|---|
| 1 | 7.5 | 12.5 |
| 2 | 18 | 30 |
| 3 | 24 | 40 |
| 4 | 33 | 55 |
| 5 | 39 | 65 |
| 6 | 48 | 80 |
| 7 | 57 | 95 |
| 8 | 63 | 105 |
| 9 | 72 | 120 |
| 10 | 78 | 130 |
| 11 | 87 | 145 |
| 12 | 96 | 160 |
| 13 | 102 | 170 |
| 14 | 111 | 185 |
| 15 | 117 | 195 |
| 16 | 126 | 210 |

| Total Number of Samples and Controls | iQ-Design Assay Primer-Probe Mix, µl | iQ Multiplex Powermix Amplification Mix, µl |
|--------------------------------------|--------------------------------------|---|
| 17 | 135 | 225 |
| 18 | 141 | 235 |
| 19 | 150 | 250 |
| 20 | 159 | 265 |
| 21 | 165 | 275 |
| 22 | 174 | 290 |
| 23 | 180 | 300 |
| 24 | 189 | 315 |
| 25 | 198 | 390 |
| 26 | 204 | 340 |
| 27 | 213 | 355 |
| 28 | 222 | 370 |
| 29 | 228 | 380 |
| 30 | 237 | 395 |
| 31 | 246 | 410 |
| 32 | 252 | 420 |

Visit bio-rad.com/Cannabis for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. IQ-CHECK is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

