



Exact Diagnostics SARS-CoV-2, Flu, RSV

Whole organism run control for multiplexed respiratory PCR assays



Exact Diagnostics SARS-CoV-2, Flu, RSV Positive and Negative Run Control

Whole organism SARS-CoV-2, Influenza A, Influenza B, Respiratory Syncytial Virus (A) run control, with a human genomic DNA background

The SARS-CoV-2, Flu, RSV Positive Run Control is manufactured with whole, intact and inactivated virus for SARS-CoV-2, Influenza A, Influenza B and RSV (A). The product is formulated in a synthetic matrix and contains human genomic DNA. The SARS-CoV-2, Flu, RSV Negative Run Control is negative for SARS-CoV-2, Influenza A, Influenza B and RSV (A) and is formulated in a synthetic matrix and contains human genomic DNA.

- Calibrated with digital droplet PCR
- Human genomic DNA (gDNA) background
- High quality matrix to ensure lot-to-lot reproducibility
- 2 year shelf life from date of manufacture

Targets

Whole Organism:

Influenza A -H1N1
 Influenza B
 Respiratory Syncytial Virus A
 SARS-CoV-2

Ordering Information

Catalog No. Description

COVFLU	Positive	5 x 1.0 mL
COVFLUNEG	Negative	5 x 1.0 mL

Refer to exactdiagnostics.com or the package insert of currently available lots for specific analyte and stability claims.



**Bio-Rad
 Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/qualitycontrol

Clinical
 Diagnostics Group

Website www.bio-rad.com/qualitycontrol **U.S.** 1 800 224 6723 **Australia** +61 (2) 9914 2800 **Austria** +43 (0) 1 877 89 01 9 **Belgium** +32 (0) 3 710 53 00 **Brazil** +55 11 3065 7550
Canada +1 514 334 4372 **China** +86 21 6169 8500 **Czech Republic** +420 241 431 660 **Denmark** +45 44 52 10 00 **Finland** +358 9 804 22 00 **France** +33 (0) 1 47 95 60 00
Germany +49 (0) 89 31884 393 **Greece** +30 210 7774396 **Hong Kong** +85 2 2789 3300 **Hungary** +36 1 459 6190 **India** +91 124 4029300 **Israel** +972 03 963 6025
Italy +39 024 94 86 600 **Japan** +81 3 6361 7070 **Korea** +82 080 007 7373 **Mexico** +52 (55) 5488 7670 **The Netherlands** +31 (0) 318 540 666 **New Zealand** +64 (9) 415 2280
Norway +47 23 38 41 30 **Poland** +48 22 331 99 99 **Portugal** +351 21 47 27 700 **Russia** +7 495 721 1404 **Singapore** +65 6415 3170 **South Africa** +27 11 442 8508
Spain +34 91 490 6580 **Sweden** +46 844 98053 **Switzerland** +41 (0) 61 717 9555 **Taiwan** +886 (2) 2578 7189 **Thailand** (662) 651 8311 **United Kingdom** +44 (0) 1923 471301

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