

Understanding COVID-19: A Global Pandemic

Britany Tang, BS Microbiologist, ATCC

Credible Leads to Incredible™



About ATCC

- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's largest, most diverse biological materials and information resource for microbes – the "gold standard"
- Innovative R&D company featuring gene editing, microbiome, NGS, advanced models
- cGMP biorepository

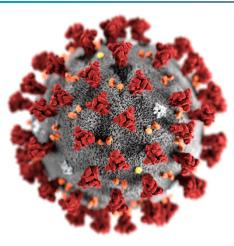
- Partner with government, industry, and academia
- Leading global supplier of authenticated cell lines, viral and microbial standards
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 450+ employees, over onethird with advanced degrees

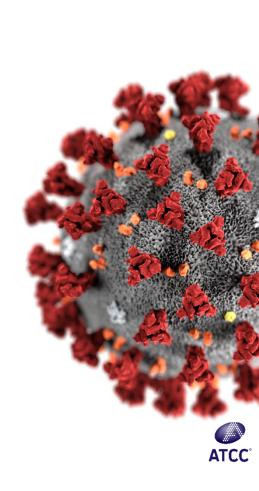


Agenda

- What is COVID-19 and SARS-CoV-2?
- What are coronaviruses?
- Diagnostics
- Vaccines
- Therapeutics
- ATCC solutions







Major areas in scientific research to combat the pandemic



Understanding the disease (COVID-19)

4



Understanding the infectious agent (SARS-CoV-2)



Diagnostics

Detection & Surveillance



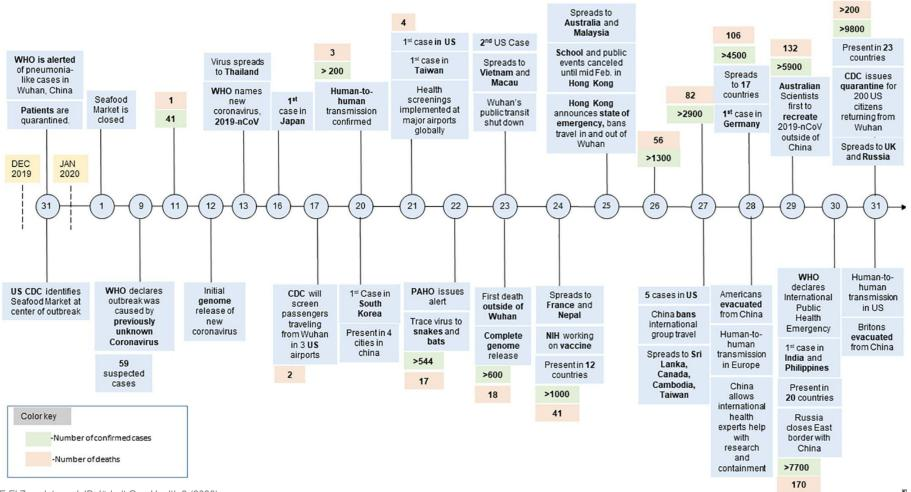
Development of Prophylactics Vaccines



Development of Therapeutics Antiviral drugs



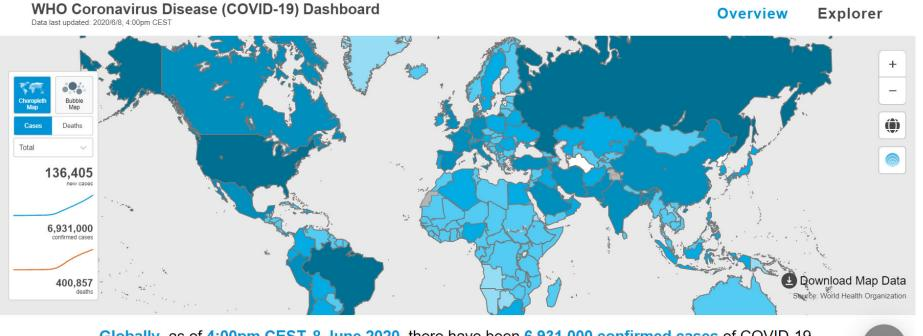
COVID-19 timeline



ATCC°

Epidemiology

- There are 215 countries, areas, or territories with cases
- More than 6.9 million cases of COVID-19 and 400,000 deaths have been reported to WHO

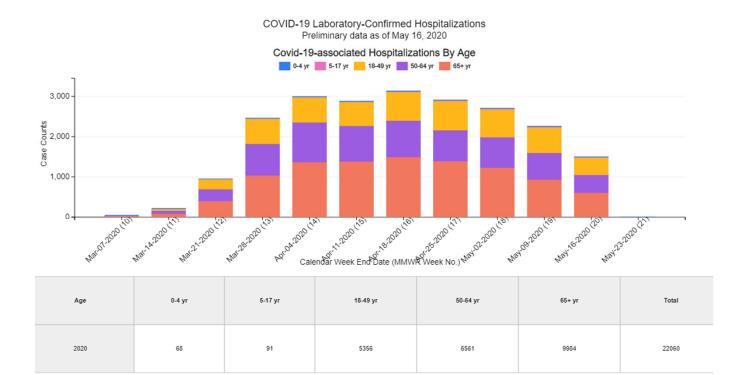


Globally, as of 4:00pm CEST, 8 June 2020, there have been 6,931,000 confirmed cases of COVID-19, including 400,857 deaths, reported to WHO.

ATCC[°]

Epidemiology

Risk factors: age groups



People of any age can be affected by COVID-19; however, **older adults** (65 years +) might be at higher risk for severe illness from COVID-19.

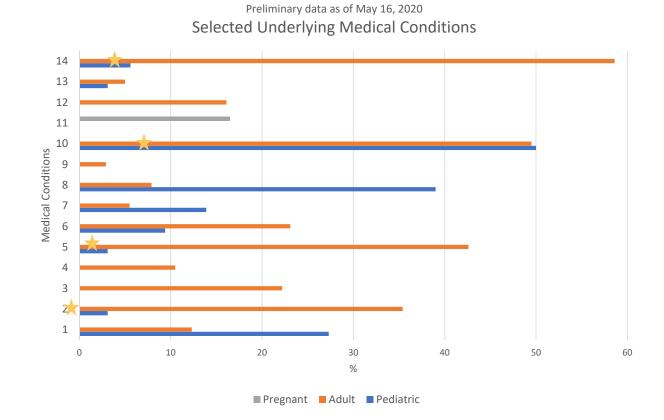
The Coronavirus Disease 2019 (COVID-19)-Associated Hospitalization Surveillance Network (COVID-NET) hospitalization data are preliminary and subject to change as more data become available. In particular, case counts and rates for recent hospital admissions are subject to lag. As data are received each week, prior case counts and rates are updated accordingly.



COVID-NET: COVID-19-Associated Hospitalization Surveillance Network, Centers for Disease Control and Prevention

Epidemiology

Risk factors: selected underlying medical conditions



COVID-19 Laboratory-Confirmed Hospitalizations

Based on the current information available and clinical expertise, **people of any age who have serious underlying medical conditions** like heart disease, lung disease, or diabetes might be at higher risk for severe illness from COVID-19.



COVID-19

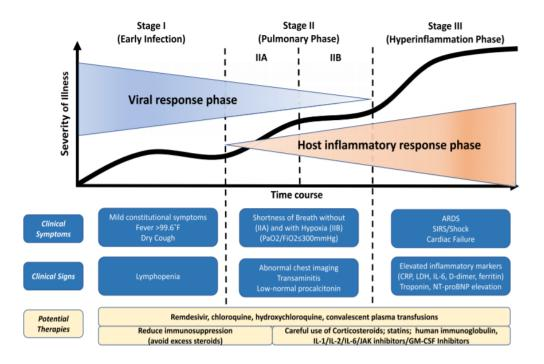
Pathogenesis & clinical manifestations

Symptoms appear 2 to 14 days after exposure

- Range from mild to severe illness with either of the following accompanying symptoms:
 - Flu or fever-like symptoms
 - Muscle pain
 - Shortness of breath or difficulty breathing
 - Temporary loss of taste or smell

Based on various risk factors, symptoms can progress to:

- Developing pneumonia in both lungs
- Extrapulmonary systemic hyperinflammation syndrome
- Acute respiratory distress syndrome (ARDS)



9

Major areas in scientific research to combat the pandemic



Understanding the disease (COVID-19)



Understanding the infectious agent (SARS-CoV-2)



Diagnostics

Detection & Surveillance



Development of Prophylactics Vaccines



Development of Therapeutics Antiviral drugs



What do we know about Coronaviruses?

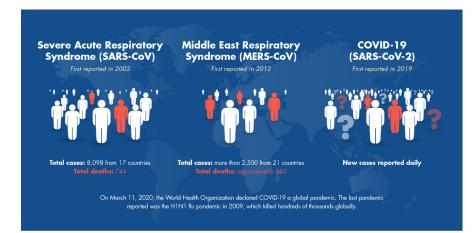


- Coronaviruses (CoVs) are large, enveloped RNA viruses of both medical and veterinary importance.
- These viruses cause a variety of diseases, including respiratory disease, enteric disease, neurological illness, and hepatitis.
- Zoonosis:
 - These viruses are described in various wildlife species such as swine, cattle, horses, camels, cats, and dogs. Many coronavirus infections are subclinical.
 - In humans, coronaviruses are included in the spectrum of viruses that cause the common cold.
 - Alphacoronaviruses (229E and NL63)
 - Betacoronaviruses (OC43 and HKU1)



Recent history of Coronavirus epidemics

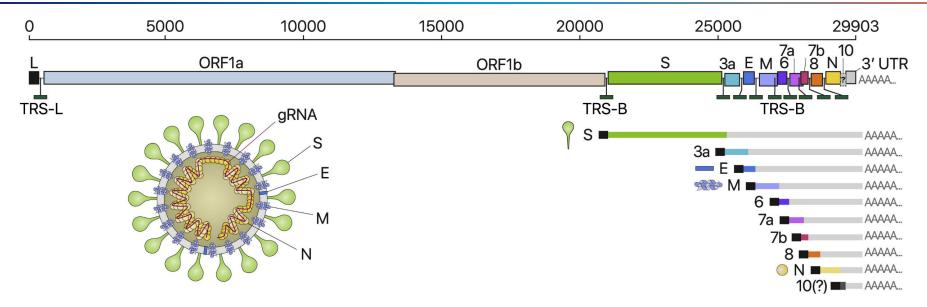
- During the past two decades, three zoonotic coronaviruses have been identified as the cause of large-scale disease outbreaks:
 - Severe Acute Respiratory Syndrome (SARS)
 - Middle East Respiratory Syndrome (MERS)
 - Swine Acute Diarrhea Syndrome (SADS)
- SARS and MERS emerged in 2003 and 2012, respectively, and caused worldwide pandemics that claimed thousands of human lives while SADS struck the swine industry in 2017.
- Common characteristics of these viruses:
 - Highly pathogenic to humans or livestock
 - They originated in bats
- It is highly likely that future SARS- or MERS-like coronavirus outbreaks will originate in bats.







Coronavirus genome biology



Coronaviruses are enveloped, positive-sense RNA viruses that are characterized by three main features:

Club-like spikes that project from their surface

An unusually large RNA genome

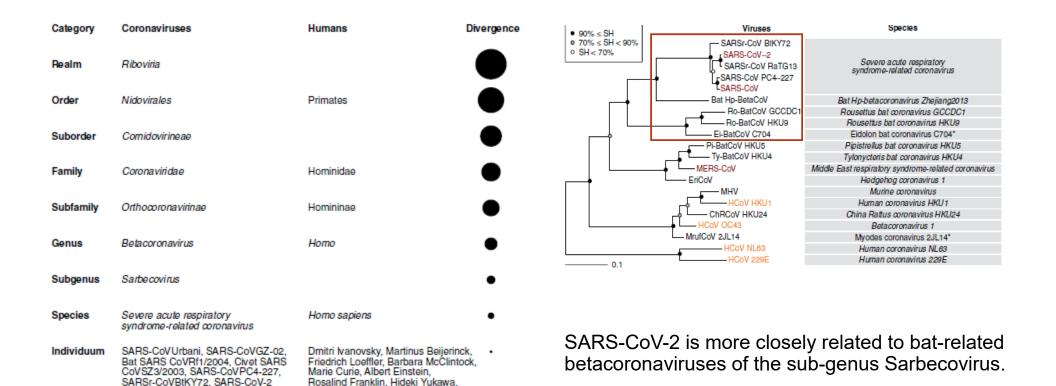
Unique replication strategy

AR Fehr & S Perlman Coronaviruses pp 1-23, 2015 D Kim, et al. Cell 181(4): 914-921, 2020



13

Phylogeny



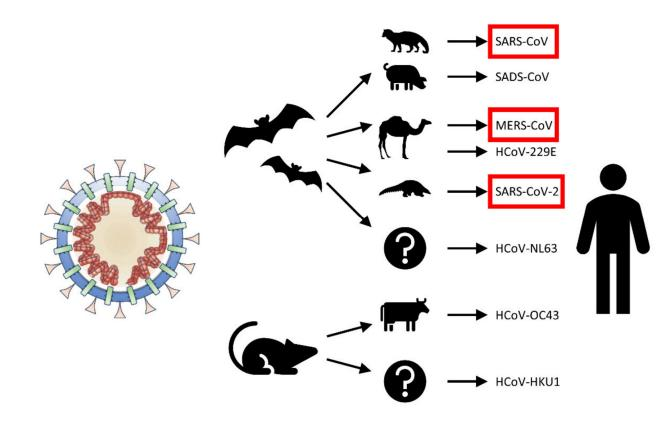


Wuhan-Hu-1, SARSr-CoVRatG13,

and so on.

and so on.

Host range

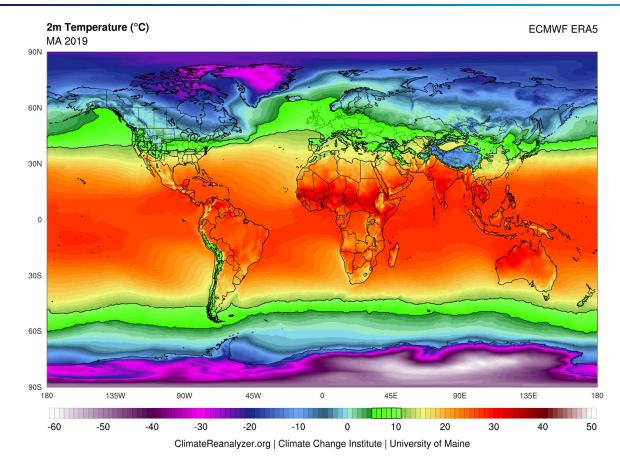


Proposed host reservoir: bats

Possibilities of intermediate hosts: Bamboo rats, pangolins, snakes, and others?



Seasonality



This map reflects the average temperature data from March-April 2019 and was developed to predict the areas that are at risk for community transmission of COVID-19. Zones at the highest risk are within the green bands.

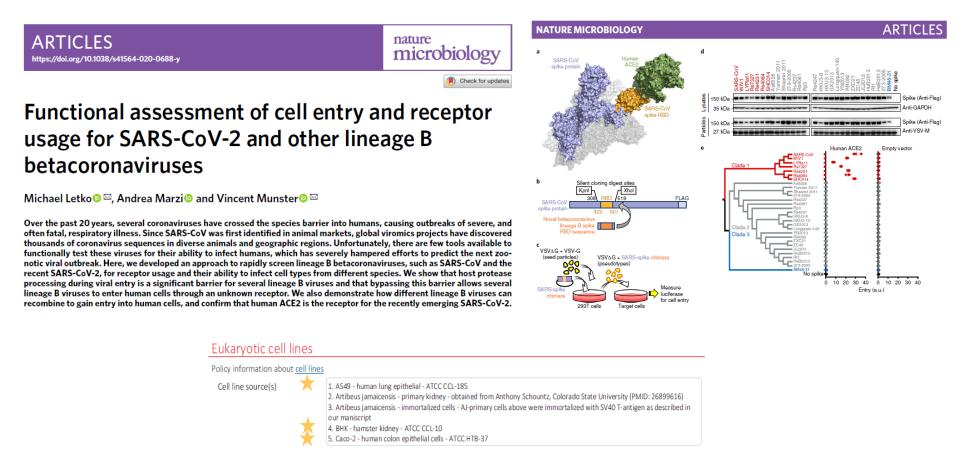
Given the consistent seasonal variation of the four endemic coronaviruses (229E, HKU1, NL63, and OC43), SARS-CoV-2 may be affected by the following factors:

- Climate
- Humidity
- Presence of UV light

Image from Climate Reanalyzer, Climate Change Institute, University of Maine, USA. Image manipulation by Cameron Gutierrez and Glenn Jameson. Z Sun, et al. International Journal of Environmental Research and Public Health 17(5): 1633, 2020 SM Kissler, et al. Science, 2020.



Discovering the molecular mechanisms of pathogenesis



ATCC

Major areas in scientific research to combat the pandemic



Understanding the disease (COVID-19)



Understanding the infectious agent (SARS-CoV-2)



Diagnostics

Detection & Surveillance



Development of Prophylactics Vaccines

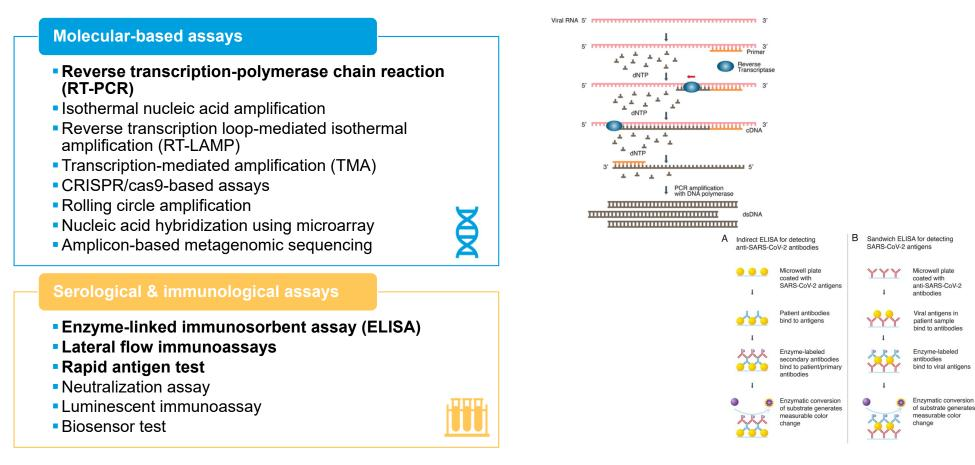


Development of Therapeutics Antiviral drugs



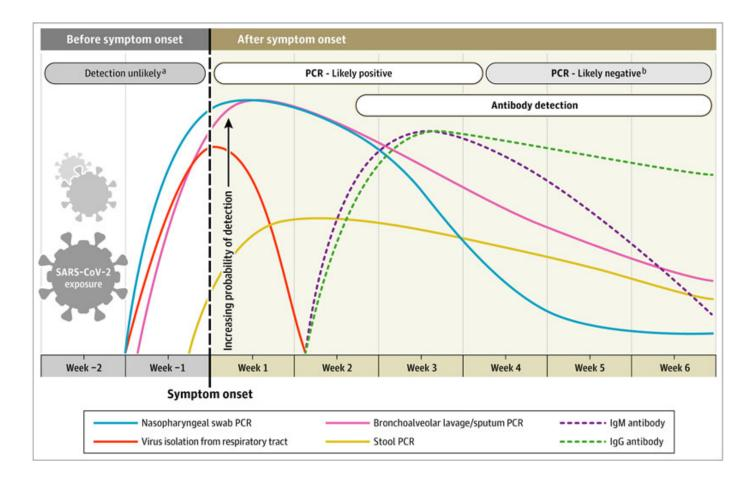
Current methods for detection & surveillance

Molecular and serological/immunological assays



ATCC°

Variation in diagnostic tests relative to symptom onset





Diagnostics

Molecular and immunological assays and targets

A majority of PCR-based assays target the ORF1ab, RNA-dependent RNA polymerase (RdRp), E, N, and spike regions. Currently authorized serological tests for SARS-CoV-2 measure IgM and/or IgG antibodies, total antibodies, or the spike protein.

• FDA:

https://www.fda.gov/medical-devices/emergencysituations-medical-devices/emergency-useauthorizations#covid19ivd

• CDC:

https://www.cdc.gov/coronavirus/2019-ncov/lab/rtpcr-panel-primer-probes.html

• FIND:

https://www.finddx.org/covid-19/pipeline/

• WHO:

https://www.who.int/who-documentsdetail/molecular-assays-to-diagnose-covid-19summary-table-of-available-protocols

| Institute | Gene targets |
|---|---|
| China CDC, China | ORF1ab and N |
| Institut Pasteur, Paris, France | Two targets in RdRP |
| US CDC, USA | Three targets in N gene |
| National Institute of Infectious Diseases, Japan | Pancorona and multiple targets, Spike protein |
| Charité, Germany | RdRP, E, N |
| HKU, Hong Kong SAR | ORF1b-nsp14, N |
| National Institute of Health, Thailand | N |

ΔΤϹϹ

Diagnostics

Validation methods & ATCC solutions





Cross-reactivity/Analytical specificity



SARS-CoV-2 Reference Materials SARS-CoV-2 Synthetic Molecular Standards Materials for Inclusivity Testing Materials for Exclusivity [Specificity] Testing



Introduction

Contains Nonbinding Recommendations

Policy for Coronavirus Disease-2019 Tests During the Public Health Emergency (Revised)

Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff

Document issued on the web on May 11, 2020.

This document supersedes "Policy for Diagnostic Tests for Coronavirus Disease-2019 during the Public Health Emergency: Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff" issued May 4, 2020.

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study



LOD/analytical sensitivity

Contains Nonbinding Recommendations

Policy for Coronavirus Disease-2019 Tests During the Public Health Emergency (Revised)

Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff

Document issued on the web on May 11, 2020.

This document supersedes "Policy for Diagnostic Tests for Coronavirus Disease-2019 during the Public Health Emergency: Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff" issued May 4, 2020.

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study

(1) Limit of Detection

FDA recommends that developers document the limit of detection (7) of their SARS-CoV-2 assay. FDA generally does not have concerns with spiking RNA or inactivated virus into artificial or real clinical matrix (e.g., Bronchoalveolar lavage [BAL] fluid, sputum, etc.) for LoD determination.

FDA recommends that developers test a dilution series of three replicates per concentration with inactivated virus on actual patient specimen, and then confirm the final concentration with 20 replicates. For this guidance, FDA defines LoD as the lowest concentration at which 19/20 replicates are positive. If multiple clinical matrices are intended for clinical testing, FDA recommends that developers submit in their EUA requests the results from the most challenging clinical matrix to FDA. For example, if testing respiratory specimens (e.g., sputum, BAL, nasopharyngeal (NP) swabs, etc.), laboratories should include only results from sputum in their EUA request.



Cross-reactivity/analytical specificity & microbial interference

Contains Nonbinding Recommendations

Policy for Coronavirus Disease-2019 Tests During the Public Health Emergency (Revised)

Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff

Document issued on the web on May 11, 2020.

This document supersedes "Policy for Diagnostic Tests for Coronavirus Disease-2019 during the Public Health Emergency: Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff" issued May 4, 2020.

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study

(3) Inclusivity

developers should document the results of an *in silico* analysis indicating the percent identity matches against publicly available SARS-CoV-2 sequences that can be detected by the proposed molecular assay. FDA anticipates that 100% of published SARS-CoV-2 sequences will be detectable with the selected primers and probes.

(4) Cross-reactivity

FDA recommends cross-reactivity wet testing on common respiratory flora and other viral pathogens at concentrations of 10^6 CFU/ml or higher for bacteria and 10^5 pfu/ml or higher for viruses, except for SARS-Coronavirus and MERS-Coronavirus, which can be accomplished by *in silico* analysis. As an alternative, FDA believes an *in silico* analysis of the assay primer and probes compared to common respiratory flora and other viral pathogens can be performed. For this guidance, FDA defines *in silico* cross-reactivity as greater than 80% homology between one of the primers/probes and any sequence present in the targeted microorganism. In addition, FDA recommends that developers follow recognized laboratory procedures in the context of the sample types intended for testing for any additional cross-reactivity testing.

Additional information for the validation of molecular diagnostics is included in the manufacturer and developers EUA templates available for download on our website.



Clinical agreement

Contains Nonbinding Recommendations

Policy for Coronavirus Disease-2019 Tests During the Public Health Emergency (Revised)

Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff

Document issued on the web on May 11, 2020.

This document supersedes "Policy for Diagnostic Tests for Coronavirus Disease-2019 during the Public Health Emergency: Immediately in Effect Guidance for Clinical Laboratories, Commercial Manufacturers, and Food and Drug Administration Staff" issued May 4, 2020.

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study

(2) Clinical Evaluation

The availability of positive samples has increased as the pandemic has progressed. As such, FDA now recommends that developers use positive clinical samples for clinical validation. Moreover, due to the increased availability of clinical samples, FDA recommends that developers confirm performance of their assay by testing a minimum of 30 positive specimens and 30 negative specimens as determined by an authorized assay. If you do not have access to clinical samples as determined by an authorized assay, contrived clinical specimens may be considered. Contrived reactive specimens can be created by spiking RNA or inactivated virus into leftover clinical specimens, of which the majority can be leftover upper respiratory specimens such as NP swabs, or lower respiratory tract specimens such as sputum, etc. If contrived samples are used, FDA recommends that twenty of the contrived clinical specimens be spiked at a concentration of 1x-2x LoD, with the remainder of specimens spanning the assay testing range. For this guidance, FDA defines the acceptance criteria for the performance as 95% agreement at 1x-2x LoD, and 100% agreement at all other concentrations and for negative specimens.

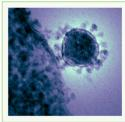


ပ္န ပြံ

ATCC portfolio

SARS-CoV-2 reference materials for inclusivity testing

| ATCC [®] No. | Product Description | Availability |
|-----------------------|---|--------------|
| VR-1986HK™ | Heat-inactivated SARS-CoV-2, Washington | Available |
| VR-1986D™ | Genomic SARS-CoV-2 RNA, Washington | Available |
| VR-1991D™ | Genomic SARS-CoV-2 RNA, Hong Kong | Available |
| VR-1992D™ | Genomic SARS-CoV-2 RNA, Italy | Available |
| VR-1994D™ | Genomic SARS-CoV-2 RNA, Germany | June 2020 |

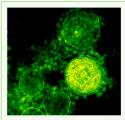


Heat-inactivated SARS-CoV-2

When developing and using a novel detection assay, researchers need access to clinically relevant positive controls to ensure the reliability and accuracy of their results. To meet this need, ATCC has developed a heat-inactivated preparation of the 2019-nCoV/USA-WA1/2020 strain (ATCC[®] VR-1986HK[™]).

- Confirmed to be inviable and non-replicative
- Quantitated by ddPCR[™]
- Useful for assays that include an extraction step

Order your preparation today at www.atcc.org/HKCoronavirus



Genomic RNA for SARS-CoV-2

Clinically relevant reference materials are an essential component of basic research and diagnostic development. That's why ATCC has made it a priority to provide heat-inactivated and genomic RNA preparations from a strain recently sourced from an infected patient in Washington state (2019-nCoV/USA-WA1/2020). This strain serves as the SARS-CoV-2 reference material for the United States.

- Fully sequenced (GenBank: MN985325.1)
- Prepared using methods known to inactivate viruses
- Suitable for RT-PCR or other RNA-based assays

Order yours today at www.atcc.org/CoronavirusRNA.

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study
- Downgraded from BSL-3 to BSL-2 (gRNA) and BSL-1 (heat-inactivated)*
- Applications:
 - Positive controls for RT-PCR or other RNAbased assays
 - Monitoring run-to-run variation within each step of the procedure, such as:
 - Nucleic acid extraction
 - Process verification
 - Amplification



www.atcc.org/coronavirus

<mark>မှု ပ</mark>ြ

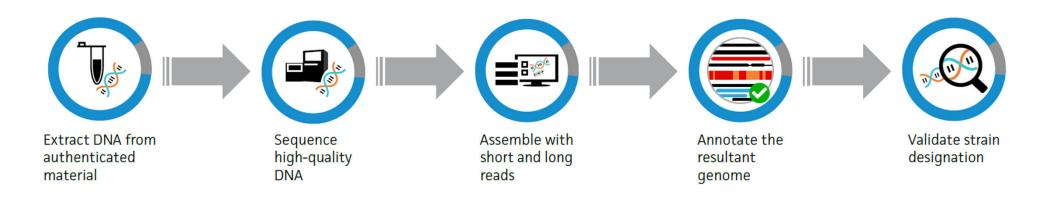
ATCC°

ATCC portfolio

SARS-CoV-2 reference materials for inclusivity testing

| ATCC [®] No. | Product Description | Availability |
|-----------------------|------------------------------------|--------------|
| VR-1986D™ | Genomic SARS-CoV-2 RNA, Washington | Available |
| VR-1991D™ | Genomic SARS-CoV-2 RNA, Hong Kong | Available |
| VR-1992D™ | Genomic SARS-CoV-2 RNA, Italy | Available |

- The genome of each strain is sequenced and assembled using our standardized workflow
- Genes are annotated and the species identity is confirmed
- Annotated genome sequences are provided on the ATCC Genome Portal.



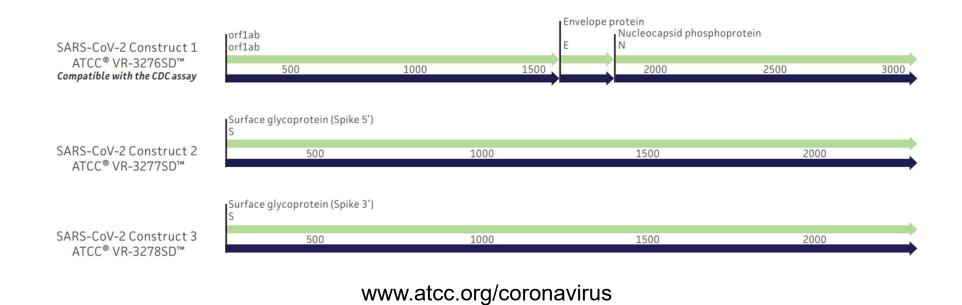
genomes.atcc.org

ATCC portfolio

Synthetic molecular standards

| ATCC [®] No. | Product Description | Availability |
|-----------------------|---|--------------|
| VR-3276SD™ | Quantitative Synthetic SARS-CoV-2 RNA: ORF, E, N | Available |
| VR-3277SD™ | Quantitative Synthetic SARS-CoV-2 RNA: Spike 5' | Available |
| VR-3278SD™ | Quantitative Synthetic SARS-CoV-2 RNA: Spike 3' | Available |
| VR-3279SD™ | Quantitative Synthetic SARS-CoV-2 RNA: nsp9, nsp12 (RdRp) | June 2020 |

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study

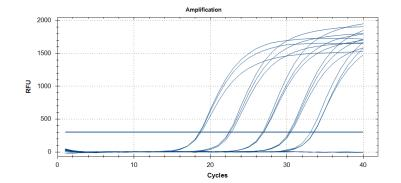


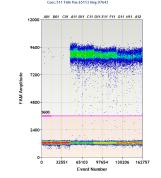


ATCC portfolio

Synthetic molecular standards

| ATCC [®] No. | Product Description | Compatible Assays |
|-----------------------|---|---|
| VR-3276SD™ | Quantitative Synthetic SARS-CoV-2 RNA: ORF, E, N | China CDC Primers and probes for detection 2019-nCoV (24 January 2020) Diagnostic detection of Wuhan coronavirus 2019 by real-time RT-PCR – Charité, Berlin Germany (17 January 2020) Detection of 2019 novel coronavirus (2019-nCoV) in suspected human cases by RT-PCR – Hong Kong University (23 January 2020) PCR and sequencing protocol for 2019-nCoV - Department of Medical Sciences, Ministry of Public Health, Thailand (Updated 28 January 2020) US CDC Real-Time RT-PCR Panel for Detection 2019-Novel Coronavirus (28 January 2020) US CDC panel primer and probes– U.S. CDC, USA (28 January 2020) |
| VR-3277SD™ | Quantitative Synthetic SARS-CoV-2 RNA: Spike 5' | Detection of WN-Human1 sequence from clinical specimen. – National Institute of Infectious Diseases Japan (17 January 2020) |
| VR-3278SD™ | Quantitative Synthetic SARS-CoV-2 RNA: Spike 3' | PCR and sequencing protocols for 2019-nCoV- National Institute of Infectious Diseases Japan (24 January 2020) |
| VR-3279SD™ | Quantitative Synthetic SARS-CoV-2 RNA: nsp9, nsp12 (RdRp) | Diagnostic detection of Wuhan coronavirus 2019 by real-time RT-PCR – Charité, Berlin Germany (17 January 2020) Real-time RT-PCR assays for the detection of SARS-CoV-2 - Institut Pasteur, Paris (02 March 2020) |





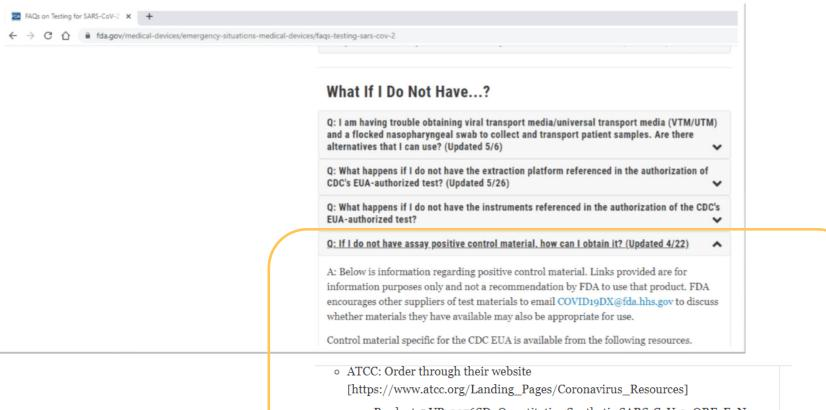
Applications:

- Positive controls for RT-PCR or other RNA-based assays
- Generation of a standard curve for quantitative RT-PCR to determine viral load
- Monitoring run-to-run, assay-to-assay, and lot-to-lot variation within each step of the procedure, such as:
 - Process verification
 - Amplification
- Assay development, verification, and validation
- To assign a genome copy number to secondary calibrators – for example, to establish a ratio of plaque or colony forming units to genome copies
- Can be used in BSL-1

Ĝ

ATCC°

ATCC Portfolio – Synthetic Molecular Standards



- Product # VR-3276SD: Quantitative Synthetic SARS-CoV-2: ORF, E, N
- Product # VR-3278SD: Quantitative Synthetic SARS-CoV-2 RNA: Spike 3'



ATCC portfolio

Exclusivity [specificity] testing materials

Related Viruses & Nucleic Acids

Human coronavirus 229E Human coronavirus OC43 Human coronavirus HKU1 Human coronavirus NL63 SARS-CoV MERS-CoV Quantitative genomic RNA from human coronavirus 229E Quantitative synthetic human coronavirus NL63 RNA Quantitative synthetic human coronavirus HKU1 RNA Quantitative synthetic MERS-CoV RNA Quantitative synthetic SARS-CoV [2003] RNA RNA from Betacoronavirus 1 OC43 RNA from human coronavirus 229E

Bacteria

Mycobacterium tuberculosis Streptococcus pyogenes Bordetella pertussis Mycoplasma pneumoniae Candida albicans Pseudomonas aeruginosa Staphylococcus epidermis Chlamydia pneumoniae Haemophilus influenzae Legionella pneumophila

Non-related Viruses

Adenovirus (e.g., C1 Ad. 71) Human metapneumovirus (hMPV) Parainfluenza virus 1-4 Influenza A & B Enterovirus (e.g., EV68) Respiratory syncytial virus Rhinovirus Measles Mumps Rubella virus Coxsackie virus Echovirus

- Limit of Detection/Analytical Sensitivity
- Cross-reactivity/Analytical Specificity
- Microbial Interference
- Clinical Agreement Study
- Having access to a variety of coronavirus strains is essential for establishing the inclusivity and exclusivity of a novel assay.
- To support this need, ATCC provides microbial strains that have a wide spectrum of temporal and geographical diversity.

www.atcc.org/coronavirus



Major areas in scientific research to combat the pandemic



Understanding the disease (COVID-19)



Understanding the infectious agent (SARS-CoV-2)



Diagnostics

Detection & Surveillance



Development of Prophylactics Vaccines

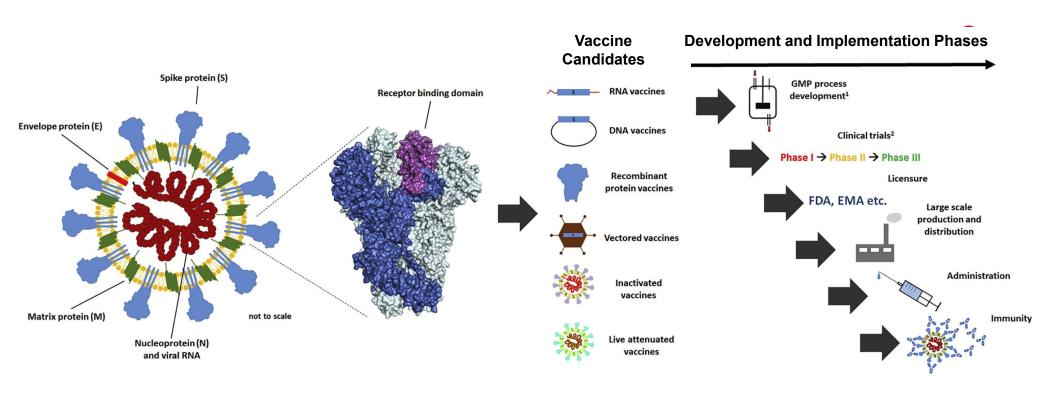


Development of Therapeutics Antiviral drugs



Vaccine development

Platforms



ATCC

Vaccine development

ATCC biological materials – Meeting the need

Cell Lines for SARS-CoV-2 Propagation

- Vero E6 cells (ATCC[®] CRL-1586[™])
- Vero CCL-81 (ATCC[®] CCL-81[™])
- MRC-5 (ATCC[®] CCL-171[™])
- HCT-8 (ATCC[®] CCL-244[™])
- Media and reagents to support cellular growth

Cell Lines for Enhanced Virus Production

- STAT1 knockout cell lines capable of producing high-titer viral stocks:
 - Vero.STAT1 KO (ATCC® CCL-81-VHG™)
 - MDCK.STAT1KO (ATCC® CCL-34-VHG™)
- Additional cell lines can be found on ATCC's website

Volume 26, Number 6—June 2020

Research

Severe Acute Respiratory Syndrome Coronavirus 2 from Patient with Coronavirus Disease, United States

Jennifer Harcourt¹, Azaibi Tamin¹, Xiaoyan Lu, Shifaq Kamili, Senthil K. Sakthivel, Janna Murray, Krista Queen, Ying Tao, Clinton R. Paden, Jing Zhang, Yan Li, Anna Uehara, Haibin Wang, Cynthia Goldsmith, Hannah A. Bullock, Lijuan Wang, Brett Whitaker, Brian Lynch, Rashi Gautam, Craig Schindewolf, Kumari G. Lokugamage, Dionna Scharton, Jessica A. Plante, Divya Mirchandani, Steven G. Widen, Krishna Narayanan, Shinji Makino, Thomas G. Ksiazek, Kenneth S. Plante, Scott C. Weaver, Stephen Lindstrom, Suxiang Tong, Vineet D. Menachery², and Natalie J. Thornburg³

Author affiliations: Centers for Disease Control and Prevention, Atlanta, Georgia, USA (J. Harcourt, A. Tamin, X. Lu, K. Queen, Y. Tao, C.R. Paden, Y. Li, C. Goldsmith, B. Whitaker, R. Gautam, S. Lindstrom, S. Tong, N.J. Thornburg): Eagle Medical Services, Atlanta (S. Kamili, S.K. Sakthivel, J. Murray, B. Lynch); IHRC, Atlanta (J. Zhang, H. Wang); Oak Ridge Institute for Science and Education, Oak Ridge, Tennessee, USA (A. Uehara); Synergy America, Inc., Atlanta (I.A. Bullock, L. Wang); University of Texas Medical Branch, Galveston, Texas, USA (C. Schindewolf, K.G. Lokugamage, D. Mirchandani, S. Widen, K. Narayanan, S. Makino, T.G. Ksiazek, S.C. Weaver, V.D. Menachery); World Reference Center for Emerging Viruses and Arboviruses, Galveston (D. Scharton, J.A. Plante, T.G. Ksiazek, K.S. Plante, S.C. Weaver, V.D. Menachery);

Cite This Article

Abstract

The etiologic agent of an outbreak of pneumonia in Wuhan, China, was identified as severe acute respiratory syndrome coronavirus 2 in January 2020. A patient in the United States was given a diagnosis of infection with this virus by the state of Washington and the US Centers for Disease Control and Prevention on January 20, 2020. We isolated virus from nasopharyngeal and oropharyngeal specimens from this patient and characterized the viral sequence, replication properties, and cell culture tropism. We found that the

virus replicates to high titer in Vero-CCL81 cells and Vero E6 cells in the absence of trypsin. We also deposited the virus into 2 virus repositories, making it broadly available to the public health and research communities. We hope that open access to this reagent will expedite development of medical countermeasures.

virus replicates to high titer in Vero-CCL81 cells and Vero E6 cells in the absence of trypsin





On This Page

Methods

Results

Discussion

Figures

Figure 1

Figure 2

Figure 3

Figure 4

Podcast

Downloads

Listen to audio/Podcast

Cite This Article

Vaccine development

Published research & news

Immunogenicity and protective efficacy in monkeys of purified inactivated Vero-cell SARS vaccine

Ede Qin^{a,*,1}, Huiying Shi^{b,1}, Lin Tang^{c,1}, Cuie Wang^{a,1}, Guohui Chang^a, Zhifen Ding^b, Kai Zhao^b, Jian Wang^c, Ze Chen^c, Man Yu^a, Bingyin Si^a, Jianyuan Liu^b, Donglai Wu^d, Xiaojie Cheng^c, Baoan Yang^a, Wenming Peng^a, Qingwen Meng^d, Bohua Liu^a, Weiguo Han^a, Xunnan Yin^d, Hongyuan Duan^a, Dawei Zhan^a, Long Tian^b, Shuangli Li^c, Jinsong Wu^a, Gang Tan^a, Yi Li^b, Yuchuan Li^a, Yonggang Liu^d, Hong Liu^a, Fushuang Lv^a, Yu Zhang^a, Xiangang Kong^d, Baochang Fan^a, Tao Jiang^a, Shuli Xu^c, Xiaomei M_{Research paper}

Xiaohong Wu^a, Yongqiang Deng^a, Min Zhao^b, Qingyu Microneedle array delivered recombinant coronavirus vaccines:

^a Institute of Microbiology and Epidemiology, Academy of Military Medical Sciences. No. 20 Dongda Street, Feng ^b National Vaccine and Serum Institute, No. 4 Nanti, Sanjianfang, Chaoyang District, Beijing ^c Beijing Genomics Institute (BGI), Chinese Academy of Sciences, I-Zone, Shunyi, Beijing I ^d Harbin Institute of Veterinary Medicine, The Chinese Academy of Agricultural Sciences, No. Eun Kim³, Geza Erdos^b, Shaohua Huang^a, Thomas W. Kenniston^a, Stephen C. Balmert^b,

Nangang District, Harbin 150001, PR China Received 2 September 2004; accepted 12 June 2005

Received 2 September 2004; accepted 12 June 2005 Available online 12 September 2005

Vero cells (ATCC[®] CCL-81[™])

<u>J Virol</u>. 2005 Feb; 79(3): 1635–1644. doi: <u>10.1128/JVI.79.3.1635-1644.2005</u>

Molecular and Biological Characterization of Human Monoclonal Antibodies Binding to the Spike and Nucleocapsid Proteins of Severe Acute Respiratory Syndrome Coronavirus

Edward N. van den Brink,¹ Jan ter Meulen,¹ Freek Cox,¹ Mandy A. C. Jongeneelen,¹ Alexandra Thijsse,¹ Mark Throsby,¹ Wilfred E. Marissen,¹ Pauline M. L. Rood,¹ Alexander B. H. Bakker,¹ Hans R. Gelderblom,² Byron E. Martina,³ Albert D. M. E. Osterhaus,³ Wolfgang Preiser,⁴ Hans Wilhelm Doerr,⁴ John de Kruif,¹ and Jaap Goudsmit^{1,*}

<u>Proc Natl Acad Sci U S A</u>. 2007 Jul 17; 104(29): 12123–12128. Published online 2007 Jul 9. doi: <u>10.1073/pnas.0701000104</u> Medical Sciences

Cara Donahue Carey^b, V. Stalin Raj^{e,1}, Michael W. Epperly^c, William B. Klimstra^d,

Department of Bioengineering, Swanson School of Engineering, University of Pittsburgh, Pittsburgh, PA 15231, USA

Chinical and Translational Science Institute, University of Pittsburgh, Pittsburgh, PA 15213, USA ¹⁰ The McGowan Institute for Regenerative Medicine, University of Pittsburgh, PA 15219, USA

ARTICLE INFO

Received 16 March 2020 Revised 18 March 2020

Accepted 18 March 2020

Available online xxx

Article History

Bart L. Haagmans^e, Emrullah Korkmaz^{b,f}, Louis D. Falo Jr.^{b,f,g,h,*}, Andrea Gambotto^{a,**}

ARSTRACT

¹ Department of Surgery, University of Pittsburgh School of Medicine, Pittsburgh, W1148 Biomedical Science Tower, 200 Lathrop St., Pennsylvania, PA 15213, USA ¹ Department of Demunologi, University of Pittsburgh School of Medicine, W1150 Biomedical Science Tower, 200 Lathrop St., Pittsburgh, PA 15213, USA ¹ Department of Demunologi, Diniversity of Pittsburgh, Pittsburgh, Pittsburgh, PA 15213, USA ² Center for Vaccine Research, Department of Immunology, University of Pittsburgh, Pittsburgh, PA 15213, USA ³ Department of Viscorience, Desman Medical Center Rotterdum, Rotterdum, the Ventherhands

Background: Coronaviruses pose a serio

drome (SARS), Middle East Respiratory

Coronavirus (MERS-CoV), and the no named SARS-CoV-2, are the causative a

Safe vaccines that rapidly induce potent

PMCID: PMC1924550 PMID: <u>17620608</u>

Potent cross-reactive neutralization of SARS coronavirus isolates by human monoclonal antibodies

<u>He</u>,[‡] <u>Anjeanette Roberts</u>,[§] <u>Tim Sheahan</u>,[¶] <u>Xiaodong Xiao</u>,^{*} <u>kx</u>,[¶] <u>Igor A. Sidorov</u>,^{*} <u>Davide Corti</u>,^{**} <u>Leatrice Vogel</u>,[§] <u>Yang Feng</u>,^{*} <u>itonio Lanzavecchia</u>,^{**} <u>Kristopher M. Curtis</u>,[¶] <u>Gary J. Nabel</u>,^{††} <u>)imitrov</u>^{*}§§

> Nat Med. 2004 Aug;10(8):871-5. doi: 10.1038/nm1080. Epub 2004 Jul 11.

An Efficient Method to Make Human Monoclonal Antibodies From Memory B Cells: Potent Neutralization of SARS Coronavirus

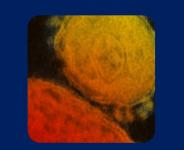
Elisabetta Traggiai ¹, Stephan Becker, Kanta Subbarao, Larissa Kolesnikova, Yasushi Uematsu, Maria Rita Gismondo, Brian R Murphy, Rino Rappuoli, Antonio Lanzavecchia

Affiliations + expand PMID: 15247913 PMCID: PMC7095806 DOI: 10.1038/nm1080

Major areas in scientific research to combat the pandemic



Understanding the disease (COVID-19)



Understanding the infectious agent (SARS-CoV-2)



Diagnostics

Detection & Surveillance



Development of Prophylactics Vaccines



Development of Therapeutics Antiviral drugs



Therapeutics

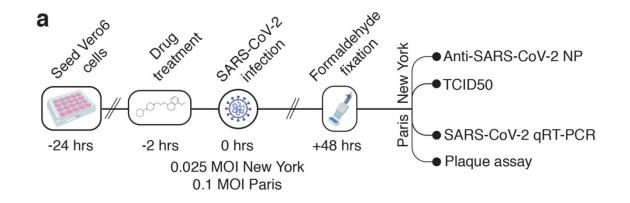
ATCC biological materials – Meeting the need

Cell Lines for SARS-CoV-2 Propagation

- Vero E6 cells (ATCC[®] CRL-1586[™])
- Vero CCL-81 (ATCC[®] CCL-81[™])
- MRC-5 (ATCC[®] CCL-171[™])
- HCT-8 (ATCC[®] CCL-244[™])
- Media and reagents to support cellular growth

Cell Lines for Enhanced Virus Production

- STAT1 knockout cell lines capable of producing high-titer viral stocks:
 - Vero.STAT1 KO (ATCC[®] CCL-81-VHG[™])
 - MDCK.STAT1KO (ATCC[®] CCL-34-VHG[™])



www.atcc.org/coronavirus



DE Gordon, et al. Nature https://doi.org/10.1038/s41586-020-2286-9 (2020)

Therapeutics

Published research & news

Nafamostat Shows High Antiviral Efficacy

Institut Pasteur Korea Finds COVID-19 Treatment Candidate 600 Time than Remdesivir

By Choi Moon-hee 🕴 🕑 May 15, 2020, 11:25

Accelerated Article Preview

A SARS-CoV-2 protein interaction map reveals targets for drug repurposing

| Received: 23 March 2020 | |
|---|--|
| Accepted: 22 April 2020 | |
| Accelerated Article Preview Published online 30 April 2020 | |

David E. Gordon, Gwendolyn M. Jang, Mehdi Bouhaddou, Jiewei Xu, Kirsten Obernier, Kris M White, Matthew J. O'Meara, Veronica V. Rezelj, Jeffrey Z. Guo, Danielle L. Swaney, Tia A. Tummino, Ruth Huettenhain, Robyn M. Kaake, Alicia L. Richards, Berlin Tutuncuoglu, Helene Foussard, Jyoti Batra, Kelsey Haas, Maya Modak, Minkyu Kim, Paige Haas, Benjamin J. Polacco, Hannes Braberg, Jacqueline M. Fabius, Manon Eckhardt, Margaret Soucheray, Verden M. Berlin M. Berline M. Fabius, Manon Eckhardt, Margaret Soucheray,

Vero E6 cells (ATCC[®] CCL-1586™)

Cell Research

www.nature.com/cr www.cell-research.com

LETTER TO THE EDITOR OPEN

Remdesivir and chloroquine effectively inhibit the recently emerged novel coronavirus (2019-nCoV) in vitro

Cell Research (2020) 30:269-271; https://doi.org/10.1038/s41422-020-0282-0



Antiviral Research 178 (2020) 104786

Contents lists available at ScienceDirect
Antiviral Research
journal homepage: www.elsevier.com/locate/antiviral

Short Communication

Remdesivir, lopinavir, emetine, and homoharringtonine inhibit SARS-CoV-2 replication in vitro



Ka-Tim Choy^a, Alvina Yin-Lam Wong^a, Prathanporn Kaewpreedee^a, Sin Fun Sia^a, Dongdong Chen^a, Kenrie Pui Yan Hui^a, Daniel Ka Wing Chu^a, Michael Chi Wai Chan^a, Peter Pak-Hang Cheung^b, Xuhui Huang^b, Malik Peiris^a, Hui-Ling Yen^{a,*}

^aSchool of Public Health, LKS Faculty of Medicine, The University of Hong Kong, Hong Kong SAR, China
^bDepartment of Chemistry, Hong Kong University of Science and Technology, Hong Kong SAR, China

Vero E6 cells (ATCC[®] CCL-1586[™])

Correspondence Open Access Published: 18 March 2020

Hydroxychloroquine, a less toxic derivative of chloroquine, is effective in inhibiting SARS-CoV-2 infection in vitro

Jia Liu, Ruiyuan Cao, Mingyue Xu, Xi Wang, Huanyu Zhang, Hengrui Hu, Yufeng Li, Zhihong Hu ⊠, Wu Zhong ⊠ & Manli Wang ⊠

Cell Discovery 6, Article number: 16 (2020) Cite this article

Vero E6 cells (ATCC[®] CCL-1586[™])

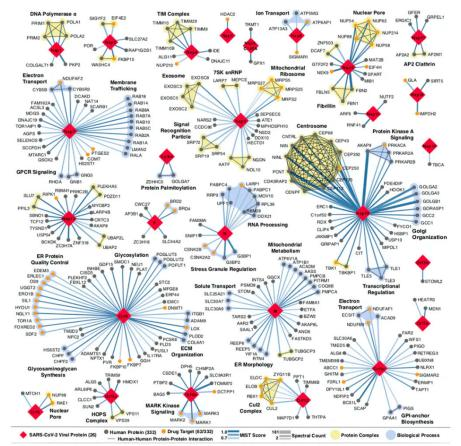


nature https://doi.org/10.1038/441586-020-22
Accelerated Article Preview

Therapeutics

A SARS-CoV-2 protein interaction map reveals targets for drug repurposing

Antiviral drug targets



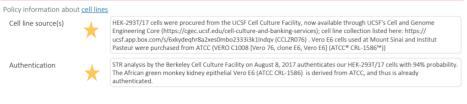
Two groups of drugs that affect the virus and the two different ways that it happens were identified from screening old drugs:

- Disrupting translation of the virus
 - Ternatin-4
 - Zotatifin
 - Plitidepsin

Sigma receptors

- Two antipsychotics: haloperidol and melperone
- Two potent antihistamines:
 - Clemastine
 - Cloperastine
- Compound PB28
- Female hormone progesterone

Eukaryotic cell lines

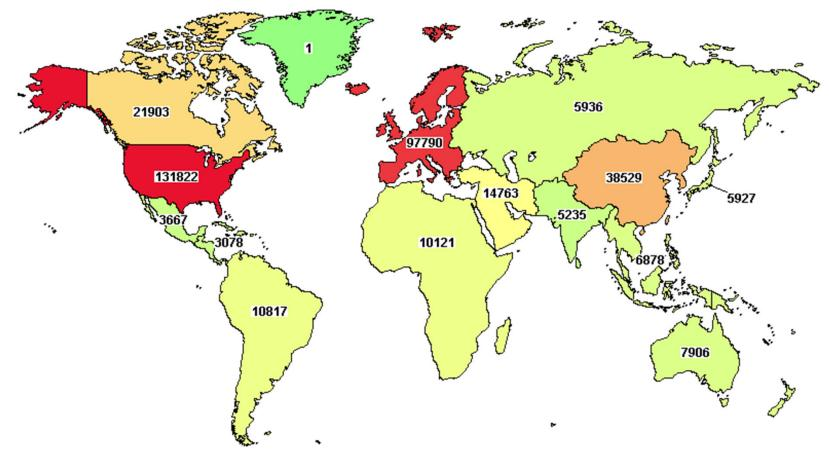


40

NIH U.S. National Library of Medicine ClinicalTrials.gov

Ongoing COVID-19 clinical trials

341,642 studies for COVID-19



ATCC

NIH. ClinicalTrials.gov, Accessed June 08, 2020.

41

Summary

- COVID-19 is a high global and public health threat.
- The development of safe and effective diagnostic methods, prophylactics, and therapeutics will depend on solid scientific research.
- ATCC has expedited scientific research by quickly providing a variety of research materials for the development of diagnostic assays, vaccines, and therapeutics.
- We must proactively protect ourselves and our community from COVID-19 infection.
- Everyone is a part of the solution.





Coming soon!



Viral Metagenomics and the Use of Standards: From Biology to Clinical Applications

Presented by Tasha M. Santiago-Rodriguez, Ph.D. 12 ET, June 25, 2020

© 2020 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise.

