

## **A high-throughput neutralizing antibody assay for COVID-19 diagnosis and vaccine evaluation**

Antonio E. Muruato<sup>1,2#</sup>, Camila R. Fontes-Garfias<sup>1#</sup>, Ping Ren<sup>3#</sup>, Mariano A. Garcia-Blanco<sup>1,4,5</sup>,  
Vineet D. Menachery<sup>2,3,6</sup>, Xuping Xie<sup>1\*</sup>, Pei-Yong Shi<sup>1,6,7,8\*</sup>

<sup>1</sup>Department of Biochemistry and Molecular Biology, University of Texas Medical Branch, Galveston TX, USA

<sup>2</sup>Department of Microbiology and Immunology, University of Texas Medical Branch, Galveston TX, USA

<sup>3</sup>Department of Pathology, University of Texas Medical Branch, Galveston TX, USA

<sup>4</sup>Programme in Emerging Infectious Diseases, Duke-NUS Medical School, Singapore, Singapore

<sup>5</sup>Department of Internal Medicine, University of Texas Medical Branch, Galveston, TX, USA

<sup>6</sup>Institute for Human Infections and Immunity, University of Texas Medical Branch, Galveston, TX, USA

<sup>7</sup>Sealy Institute for Vaccine Sciences, University of Texas Medical Branch, Galveston, TX, USA

<sup>8</sup>Sealy Center for Structural Biology & Molecular Biophysics, University of Texas Medical Branch, Galveston, TX, USA

\*C.F.-G., A.E.M., and P.R. contributed equally to this study

\*Correspondence: X.X. ([xuxie@UTMB.edu](mailto:xuxie@UTMB.edu)) or P.-Y.S. ([peshi@UTMB.edu](mailto:peshi@UTMB.edu))

**Running title:** A high-throughput SARS-CoV-2 neutralizing assay

**Keywords:** SARS-CoV-2, COVID-19, SARS-CoV, diagnosis, vaccine, coronavirus

## 1 **Abstract**

2 Virus neutralization remains the gold standard for determining antibody efficacy.  
3 Therefore, a high-throughput assay to measure SARS-CoV-2 neutralizing antibodies is urgently  
4 needed for COVID-19 serodiagnosis, convalescent plasma therapy, and vaccine development.  
5 Here we report on a fluorescence-based SARS-CoV-2 neutralization assay that detects SARS-  
6 CoV-2 neutralizing antibodies in COVID-19 patient specimens and yields comparable results to  
7 plaque reduction neutralizing assay, the gold standard of serological testing. Our approach  
8 offers a rapid platform that can be scaled to screen people for antibody protection from COVID-  
9 19, a key parameter necessary to safely reopen local communities.

10

## 11 **Text**

12 The ongoing coronavirus disease 2019 (COVID-19) pandemic is caused by severe acute  
13 respiratory syndrome coronavirus 2 (SARS-CoV-2), first reported in Wuhan, China in late  
14 2019<sup>1,2</sup>. As of May 18, 2020, COVID-19 has caused 4.8 million confirmed infections and over  
15 318,028 deaths worldwide (<https://www.worldometers.info/coronavirus/>). Many areas of the  
16 world have been in lockdown mode to curb the viral transmission, but the reality is that COVID-  
17 19 is here to stay until a safe and efficacious vaccine becomes available. The pandemic's  
18 catastrophic economic impact is pushing governments to reopen their economies, and this  
19 creates a public health quandary. At this time, our only option is to minimize viral transmission  
20 through social distancing and contact tracing, which relies on the diagnosis of viral RNA through  
21 RT-PCR (<https://www.fda.gov/media/134922/download>). Proper public health policy would be  
22 greatly enhanced if we had a reliable and facile assay to measure the immune protection among  
23 COVID-19 recovered patients.

24 Coronavirus infections typically induce neutralizing antibody responses<sup>3</sup>. The  
25 seroconversion rates in COVID-19 patients are 50% and 100% on day 7 and 14 post symptom

26 onset, respectively<sup>4</sup>. Given the unknown scale of asymptomatic infections, there is a pressing  
27 need for serological diagnosis to determine the real number of infections. Such information is  
28 essential for defining the case-fatality rate and for making the policy on the scale and duration of  
29 social lockdowns. The serological assays are also required to identify donors with high-titers for  
30 convalescent plasma for therapy, and to define correlates of protection from SARS-CoV-2.  
31 While viral RNA-based testing for active infection is the current standard, surveying antibody  
32 protection is a necessary part of any return to social normality.

33 For serodiagnosis, several COVID-19 assay platforms have achieved FDA emergency  
34 use authorizations (EUA), including ELISA<sup>5</sup> (<https://www.fda.gov/media/137029/download>),  
35 lateral flow immunoassay (<https://www.fda.gov/media/136625/download>), and Microsphere  
36 Immunoassay (<https://www.fda.gov/media/137541/download>). These assays measure antibody  
37 binding to SARS-CoV-2 spike protein. Since not all spike-binding antibodies can block viral  
38 infection, these assay platforms do not functionally measure antibody inhibition of SARS-CoV-2  
39 infection. An ideal serological assay should measure neutralizing antibody levels, which should  
40 predict protection from reinfection. Conventionally, neutralizing antibodies are measured by  
41 plaque reduction neutralization test (PRNT). Although PRNT and ELISA results generally  
42 correlate with each other, the lack of complete fidelity of ELISA continues to make PRNT the  
43 gold-standard for determining immune protection<sup>6,7</sup>. However, due to its low throughput, PRNT  
44 is not practical for large scale serodiagnosis and vaccine evaluation. This is a major gap for  
45 COVID-19 surveillance and vaccine development.

46 To address the above gap, we developed a fluorescence-based assay that rapidly and  
47 reliably measures neutralization of a reporter SARS-CoV-2 by antibodies from patient  
48 specimens. The assay was built on a stable mNeonGreen SARS-CoV-2 where the  
49 mNeonGreen gene was engineered at the OFR7 of the viral genome<sup>8</sup>. Fig. 1a depicts the  
50 flowchart of the reporter neutralization assay in a 96-well format. Briefly, patient sera were

51 serially diluted and incubated with the reporter virus. After incubation at 37°C for 1 h, Vero E6  
52 cells (pre-seeded in a 96-well plate) were infected with the virus/serum mixtures at a multiplicity  
53 of infection (MOI) of 0.5. At 16 h post-infection, the mNeonGreen-positive cells were quantitated  
54 using a high-content imaging reader (Fig. 1a). Forty COVID-19 serum specimens from RT-PCR-  
55 confirmed patients and ten non-COVID-19 serum samples (archived before COVID-19  
56 emergence) were analyzed using the reporter virus. After reporter viral infection, the cells turned  
57 green in the absence of serum (Fig. 1b, bottom panel); in contrast, incubation of the reporter  
58 virus with COVID-19 patient serum decreased the number of fluorescent cells (top panel). A  
59 dose response curve was obtained between the number of fluorescent cells and the fold of  
60 serum dilution (Fig. 1c), which allowed for determination of the dilution fold that neutralized 50%  
61 of fluorescent cells (NT<sub>50</sub>). The reporter assay rapidly diagnosed fifty specimens in less than 20  
62 h: all forty COVID-19 sera (specimens 1-40) showed positive NT<sub>50</sub> of 80 to 5152, and all ten  
63 non-COVID-19 sera (specimens 41-50) showed negative NT<sub>50</sub> of <20 for (Fig. 1d).

64 To validate the reporter virus neutralization results, we performed the conventional  
65 PRNT on the same set of patient specimens. In agreement with the reporter virus results, the  
66 forty positive sera showed PRNT<sub>50</sub> of 40 to 3200, and the ten negative sera exhibited PRNT<sub>50</sub> of  
67 <20 (Fig. 1d). A strong correlation was observed between the reporter virus and PRNT results,  
68 with a correlation efficiency  $R^2$  of 0.9 (Fig. 1e). The results demonstrate that when diagnosing  
69 patient specimens, the reporter virus assay delivers neutralization results comparable to the  
70 PRNT assay, the gold standard of serological testing.

71 Next, we evaluated the specificity of reporter neutralization assay using potentially cross-  
72 reactive sera and interfering substances (Table 1). Two groups of specimens were tested for  
73 cross reactivity. Group I included 138 clinical sera from patients with antigens or antibodies  
74 against different viruses, bacteria, and parasites. Group II consisted of 19 samples with albumin,  
75 elevated bilirubin, cholesterol, rheumatoid factor, and autoimmune nuclear antibodies. None of

76 the specimens cross neutralized mNeonGreen SARS-CoV-2 (Table 1), including the four  
77 common cold coronaviruses (NL63, 229E, OC43, and HUK1). The latter result is consistent with  
78 the recent reports that sera from common cold coronavirus patients did not cross react with  
79 SARS-CoV-2<sup>5,9</sup>. However, more specimens are required to further validate the cross reactivity,  
80 particularly between SARS-CoV-2 and other human coronaviruses, including SARS-CoV-1 and  
81 MERS-CoV.

82 In this study, we developed a rapid fluorescence-based high-throughput assay for  
83 COVID-19 serodiagnosis. The reporter virus assay is superior to antigen/antibody binding  
84 assays because it measures functional SARS-CoV-2 neutralizing activity in the specimens.  
85 When diagnosing patient sera, the reporter virus assay generated NT<sub>50</sub> values comparable to  
86 the conventional PRNT assay. Compared with the PRNT assay, our reporter neutralization test  
87 has shortened the assay turnaround time by several days and increased the testing capacity to  
88 high throughput. Previously, lentiviruses or vesicular stomatitis virus (VSV) pseudotyped with  
89 SARS-CoV-2 spike protein have been reported for neutralization assays<sup>10</sup>. One weakness of the  
90 spike pseudotyped assay is that it lacks the same composition of an actual virion, including the  
91 SARS-CoV-2 M or E proteins. In addition, the spike protein conformation, either the trimer or  
92 monomer, may be different in the pseudotyped virus as compared with the authentic SARS-  
93 CoV-2 virion.

94 Since mNeonGreen SARS-CoV-2 is stable and replicates like wild-type virus, our  
95 reporter neutralization assay provides an ideal model for high-throughput serological testing. As  
96 the mNeonGreen SARS-CoV-2 grows to >10<sup>7</sup> PFU/ml in cell culture<sup>8</sup>, the reporter virus can be  
97 easily scaled up for testing large sample volumes. Besides mNeonGreen, we have begun to  
98 develop other reporter SARS-CoV-2 (e.g., luciferase or mCherry) that can also be used for such  
99 serological testing. Although the current study performed the assay in a 96-well format, the  
100 assay can be readily adapted to 384- and 1536-well formats. Despite the strengths of high

101 throughput and reliability, the current reporter neutralization assay must be performed in  
102 biosafety level 3 (BSL3) containment. Efforts are ongoing to engineer an attenuated version of  
103 SARS-CoV-2 so that the assay could be performed at a BSL2 facility. Nevertheless, the  
104 mNeonGreen reporter assay offers a rapid, high-throughput platform to test COVID-19 patient  
105 sera not previously available.

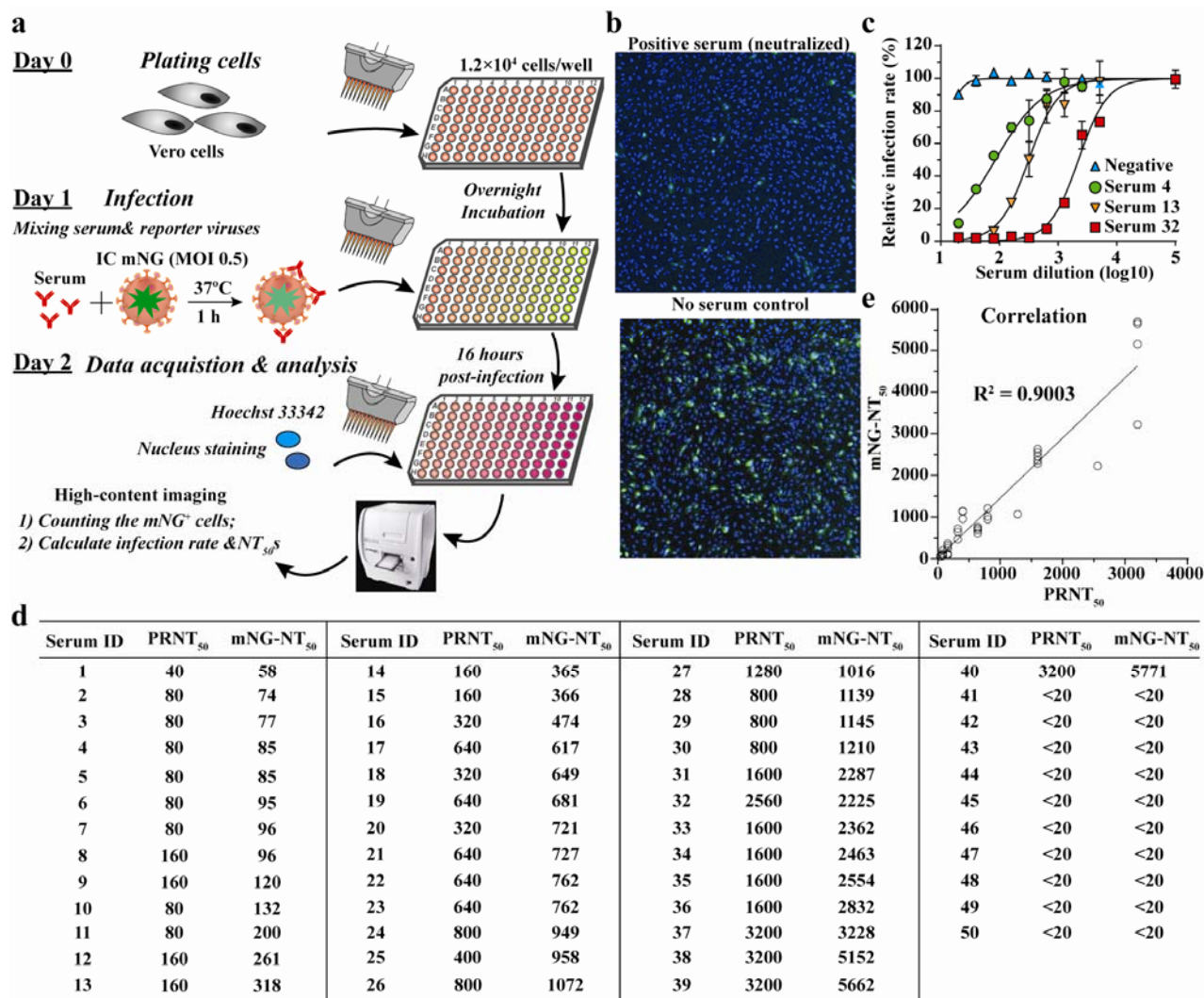
106 Because neutralizing titer is a key parameter to predict immunity, the reporter  
107 neutralization assay should be useful for high-throughput evaluation of COVID-19 vaccines and  
108 for identification of high neutralizing convalescent plasma for therapy. Indeed, treatment of  
109 severe COVID-19 patients with convalescent plasma shows clinical benefits<sup>11</sup>. For vaccine  
110 development, a standardized neutralizing assay will facilitate down selection of various  
111 candidates for clinical development. Furthermore, the reporter assay could be used over time to  
112 monitor the waning of protective neutralizing titers in COVID-19 patients and to study the  
113 correlates of protection from SARS-CoV-2. Thus, the ability to rapidly measure neutralizing  
114 antibody levels in populations is essential for guiding policymakers to reopen the economy and  
115 society, deploy healthcare workers, and prepare for SARS-CoV-2 reemergence.

116

## 117 **References**

- 118
- 119 1 Zhou, P. *et al.* A pneumonia outbreak associated with a new coronavirus of probable bat  
120 origin. *Nature* **579**, 270-273, doi:10.1038/s41586-020-2012-7 (2020).
  - 121 2 Zhu, N. *et al.* A Novel Coronavirus from Patients with Pneumonia in China, 2019. *N Engl*  
122 *J Med* **382**, 727-733, doi:10.1056/NEJMoa2001017 (2020).
  - 123 3 Huang, A. T. *et al.* A systematic review of antibody mediated immunity to coronaviruses:  
124 antibody kinetics, correlates of protection, and association of antibody responses with  
125 severity of disease. *medRxiv* (2020).
  - 126 4 Wolfel, R. *et al.* Virological assessment of hospitalized patients with COVID-2019.  
127 *Nature*, doi:10.1038/s41586-020-2196-x (2020).
  - 128 5 Amanat, F. *et al.* A serological assay to detect SARS-CoV-2 seroconversion in humans.  
129 *Nat Med*, doi:10.1038/s41591-020-0913-5 (2020).
  - 130 6 Perera, R. A. *et al.* Serological assays for severe acute respiratory syndrome  
131 coronavirus 2 (SARS-CoV-2), March 2020. *Euro Surveill* **25**, doi:10.2807/1560-  
132 7917.ES.2020.25.16.2000421 (2020).

- 133 7 Okba, N. M. A. *et al.* Severe Acute Respiratory Syndrome Coronavirus 2-Specific  
134 Antibody Responses in Coronavirus Disease 2019 Patients. *Emerg Infect Dis* **26**,  
135 doi:10.3201/eid2607.200841 (2020).
- 136 8 Xie, X. *et al.* An infectious cDNA clone of SARS-CoV-2. *Cell Host Microbe*,  
137 doi:10.1016/j.chom.2020.04.004 (2020).
- 138 9 Khan, S. *et al.* Analysis of Serologic Cross-Reactivity Between Common Human  
139 Coronaviruses and SARS-CoV-2 Using Coronavirus Antigen Microarray. *bioRxiv* (2020).
- 140 10 Nie, J. *et al.* Establishment and validation of a pseudovirus neutralization assay for  
141 SARS-CoV-2. *Emerg Microbes Infect* **9**, 680-686, doi:10.1080/22221751.2020.1743767  
142 (2020).
- 143 11 Shen, C. *et al.* Treatment of 5 Critically Ill Patients With COVID-19 With Convalescent  
144 Plasma. *JAMA*, doi:10.1001/jama.2020.4783 (2020).
- 145 12 Shan, C. *et al.* Evaluation of a Novel Reporter Virus Neutralization Test for Serological  
146 Diagnosis of Zika and Dengue Virus Infection. *J Clin Microbiol* **55**, 3028-3036,  
147 doi:10.1128/JCM.00975-17 (2017).
- 148 13 Shan, C. *et al.* A Rapid Zika Diagnostic Assay to Measure Neutralizing Antibodies in  
149 Patients. *EBioMedicine* **17**, 157-162, doi:10.1016/j.ebiom.2017.03.006 (2017).
- 150



151

152 **Figure 1. A high-throughput neutralizing antibody assay for COVID-19 diagnosis.** (a)

153 Assay flowchart. mNeonGreen SARS-CoV-2 was neutralized with COVID-19 patient sera. Vero

154 E6 cells were infected with the reporter virus/serum mixture with an MOI of 0.5. The

155 fluorescence of infected cells was quantified to estimate the NT<sub>50</sub> value for each serum. (b)

156 Representative images of reporter virus-infected Vero E6 cells. Images for a positive

157 neutralizing serum (top panel) and no serum control (bottom panel) are presented. (c)

158 Neutralization curves. Representative neutralization curves are presented for three positive sera

159 and one negative sera. (d) Summary of NT<sub>50</sub> values of fifty patient sera. The NT<sub>50</sub> values from

160 both reporter virus and conventional PRNT assays are presented. (e) Correlation analysis of



161 NT<sub>50</sub> values between the reporter virus and PRNT assays. The correlation efficiency  $R^2$  is  
162 indicated.

163 **Table 1. Cross reactivity of mNeonGreen SARS-CoV-2 neutralization assay**  
164

*Immune sera and #interfering substances	Sample number	Cross reactivity
Anti-Chikungunya virus	4	0
<i>Cryptococcus neoformans</i> antigen	2	0
Anti-Cytomegalovirus	8	0
Anti-Dengue virus	5	0
Anti-Epstein Barr Virus: capsid or nuclear antigen	8	0
Anti-Hepatitis A virus	5	0
Anti-Hepatitis B virus: surface antigen	14	0
Anti-Hepatitis C virus	3	0
Anti-Herpes simplex virus 1	7	0
Anti-Herpes simplex virus 2	5	0
Human coronavirus 229E	1	0
Human coronavirus HKU1	3	0
Human coronavirus NL63	1	0
Human coronavirus OC43	4	0
Anti-Human immunodeficiency virus 1	7	0
Human rhinovirus	3	0
Influenza B virus	2	0
Anti-Measles virus	7	0
Anti-Mumps virus	5	0
Parainfluenza virus 2	1	0
Parainfluenza virus 4	1	0
Anti-Parvovirus B19	4	0
Anti-Rubella virus	9	0
Anti-Syphilis	4	0
Anti-Toxoplasma	2	0
Anti-Typhus Fever	1	0
Varicella zoster virus	13	0
West Nile Virus	3	0
Anti-Yellow fever virus: vaccination	2	0
Anti-Zika virus	4	0
#Albumin (4.5 g/dL)	3	0
#Elevated bilirubin conjugated (>0.4 mg/dL)	3	0
#Elevated bilirubin unconjugated (>0.8 mg/dL)	3	0
#Elevated cholesterol (>200 mg/dL)	3	0
#Elevated rheumatoid factor (>100 IU/mL)	3	0
#Nuclear antibodies	4	0

165  
166 \*A total of 138 sera with antigens or antibodies against different infections (or immunizations)  
167 were tested against mNeonGreen SARS-CoV-2 neutralization assay. The immune sera are  
168 listed in alphabetical order.  
169 #Tested interfering substances and autoimmune disease nuclear antibodies.

170 **Methods**

171 **mNeonGreen SARS-CoV-2.** The virus stock of mNeonGreen SARS-CoV-2 was  
172 produced using an infectious cDNA clone of SARS-CoV-2 in which the ORF7 of the viral  
173 genome was replaced with reporter mNeonGreen gene<sup>8</sup>. After rescued from the genome-length  
174 viral RNA-electroporated cells, the viral stock was prepared by amplifying the mNeonGreen  
175 SARS-CoV-2 on Vero E6 cells for one or two rounds. The titer of the virus stock was determined  
176 by a standard plaque assay.

177 **Human sera and interfering substances.** All human serum specimens were obtained  
178 at the University of Texas Medical Branch (UTMB). All specimens were de-identified from  
179 patient information. A total of forty de-identified convalescent sera from COVID-19 patients  
180 (confirmed with viral RT-PCR positive) were tested in this study. Ten non-COVID-19 sera,  
181 collected before COVID-19 emergence<sup>12,13</sup>, were also tested in the reporter virus and PRNT  
182 assays. For testing cross reactivity, a total of 138 de-identified specimens from patients with  
183 antigens or antibodies against different viruses, bacteria, and parasites were tested in the  
184 mNeonGreen SARS-CoV-2 neutralization assay (Table 1). For testing interfering substances,  
185 nineteen de-identified serum specimens with albumin, elevated bilirubin, cholesterol, rheumatoid  
186 factor, and autoimmune nuclear antibodies were tested in the reporter neutralization assay. All  
187 human sera were heat-inactivated at 56°C for 30 min before testing.

188 **mNeonGreen SARS-CoV-2 reporter neutralization assay.** Vero E6 cells ( $1.2 \times 10^4$ ) in  
189 50  $\mu$ l of DMEM (Gibco) containing 2% FBS (Hyclone) and 100 U/ml Penicillium-Streptomycin  
190 (P/S; Gibco) were seeded in each well of black  $\mu$ CLEAR flat-bottom 96-well plate (Greiner Bio-  
191 one™). The cells were incubated overnight at 37°C with 5% CO<sub>2</sub>. On the following day, each  
192 serum was 2-fold serially diluted in 2% FBS and 100 U/ml P/S DMEM, and incubated with  
193 mNeonGreen SARS-CoV-2 at 37°C for 1 h. The virus-serum mixture was transferred to the  
194 Vero E6 cell plate with the final multiplicity of infection (MOI) of 0.5. For each serum, the starting  
195 dilution was 1/20 with nine 2-fold dilutions to the final dilution of 1/5120. After incubating the

196 infected cells at 37°C for 16 h, 25 µl of Hoechst 33342 Solution (400-fold diluted in Hank's  
197 Balanced Salt Solution; Gibco) were added to each well to stain cell nucleus. The plate was  
198 sealed with Breath-Easy sealing membrane (Diversified Biotech), incubated at 37°C for 20 min,  
199 and quantified for mNeonGreen fluorescence on Cytation™ 7 (BioTek). The raw images (2x2  
200 montage) were acquired using 4x objective, processed, and stitched using the default setting.  
201 The total cells (indicated by nucleus staining) and mNeonGreen-positive cells were quantified  
202 for each well. Infection rates were determined by dividing the mNeonGreen-positive cell number  
203 to total cell number. Relative infection rates were obtained by normalizing the infection rates of  
204 serum-treated groups to those of non-serum-treated controls. The curves of the relative  
205 infection rates versus the serum dilutions (log<sub>10</sub> values) were plotted using Prism 8 (GraphPad).  
206 A nonlinear regression method was used to determine the dilution fold that neutralized 50% of  
207 mNeonGreen fluorescence (NT<sub>50</sub>). Each serum was tested in duplicates. All mNeonGreen  
208 SARS-CoV-2 reporter neutralization assay was performed at the BSL-3 facility at UTMB.

209 **Plaque reduction neutralization test (PRNT).** Vero E6 cells (1.2×10<sup>6</sup> per well) were  
210 seeded to 6-well plates. On the following day, 100 PFU of infectious clone-derived wild-type  
211 SARS-CoV-2 was incubated with serially diluted serum (total volume of 200 µl) at 37°C for 1 h.  
212 The virus-serum mixture was added to the pre-seeded Vero E6 cells. After 1 h 37°C incubation,  
213 2 ml of 2% high gel temperature agar (SeaKem) in DMEM containing 5% FBS and 1% P/S was  
214 added to the infected cells. After 2 days of incubation, 2 ml neutral red (1 g/l in PBS; Sigma)  
215 was added to the agar-covered cells. After another 5-h incubation, neutral red was removed.  
216 Plaques were counted for NT<sub>50</sub> calculation. Each serum was tested in duplicates. The PRNT  
217 assay was performed at the BSL-3 facility at UTMB.

218 **Statistical analysis.** The correlation of the NT<sub>50</sub> values from mNeonGreen reporter  
219 SARS-CoV-2 assay and the PRNT<sub>50</sub> values from plaque neutralization assay was analyzed  
220 using a linear regression model in the software Prism 8 (GraphPad).

221

222 **Data availability**

223           The results presented in the study are available upon request from the corresponding  
224 authors. The mNeonGreen reporter SARS-CoV-2 has been deposited to the World Reference  
225 Center for Emerging Viruses and Arboviruses (<https://www.utmb.edu/wrceva>) at UTMB for  
226 distribution.

227

228 **Acknowledgements**

229           We thank colleagues at UTMB for helpful discussion during the course of this project.  
230 P.-Y.S. was supported by NIH grants AI142759, AI134907, AI145617, and UL1TR001439, and  
231 awards from the Sealy & Smith Foundation, Kleberg Foundation, John S. Dunn Foundation,  
232 Amon G. Carter Foundation, Gilson Longenbaugh Foundation, and Summerfield Robert  
233 Foundation. M.A.G.-B. was supported by NIH grant CA204806 and the Vacek Distinguished  
234 Chair. V.D.M. was supported by NIH grants U19AI100625, R00AG049092, R24AI120942, and  
235 STARs Award from the University of Texas System. A.E.M. is supported by a Clinical and  
236 Translational Science Award NRSA (TL1) Training Core (TL1TR001440) from NIH. C.R.F.-G. is  
237 supported by the predoctoral fellowship from the McLaughlin Fellowship Endowment at  
238 UTMB.

239

240 **Author contributions**

241           P.R., M.A.G.-B., V.D.M., X.X., and P.-Y.S conceived the study. A.E.M. and C.R.F.-G.  
242 performed the experiments and analyzed the results. P.R. prepared the serum specimens.  
243 M.A.G.-B., V.D.M., X.X., and P.-Y.S wrote the manuscript.

244

245 **Competing interests**

246 UTMB has filed a patent on the reverse genetic system and reporter SARS-CoV-2.