

SARS-COV-2 Nucleocapsid IgG Assay Layout Sheet
for use with Enzo Life Sciences Catalog No. ENZ-KIT193-0001

• **CALIBRATOR PREPARATION: Reconstitute each calibrator in 1mL Sample Diluent.**

Std.	Nucleocapsid IgG, ng (lyophilized)	Sample Diluent to add (μL)	Nucleocapsid IgG Conc. (ng/mL)
1	200	1000	200
2	50	1000	50
3	10	1000	10
4	0	1000	0

• **ASSAY PROTOCOL FLOW CHART:**

Well ID:	Blank A1, B1	C1 C1, D1	C2 E1, F1	C3 G1, H1	C4 A2-B2	Samples C2-H12
Standards or Samples	---	100 μL	100 μL	100 μL	100 μL	100 μL
Incub. 10 min. @ RT, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x Full well vol. (~350μL)	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Nucleocapsid Conjugate	---	100 μL	100 μL	100 μL	100 μL	100 μL
Incub. 10 min. @ RT, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x Full well vol. (~350μL)	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
TMB Substrate	100 μL	100 μL	100 μL	100 μL	100 μL	100 μL
Incub. 10 min. @ RT, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	100 μL	100 μL	100 μL	100 μL	100 μL	100 μL
Read OD at 450 nm	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

SARS-CoV-2 NUCLEOCAPSID IGG PLATE LAYOUT:

A1 Blank	A2 C4	A3	A4	A5	A6	A7	A8	A9	A10	A11	A12
B1 Blank	B2 C4	B3	B4	B5	B6	B7	B8	B9	B10	B11	B12
C1 C1	C2	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1 C1	D2	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E1 C2	E2	E3	E4	E5	E6	E7	E8	E9	E10	E11	E12
F1 C2	F2	F3	F4	F5	F6	F7	F8	F9	F10	F11	F12
G1 C3	G2	G3	G4	G5	G6	G7	G8	G9	G10	G11	G12
H1 C3	H2	H3	H4	H5	H6	H7	H8	H9	H10	H11	H12

Kit Lot No. _____ Exp. Date _____ Date _____ Tech. _____

1st Incub.: Start Time _____ Temp. _____ 3rd Incub.: Start Time _____ Temp. _____

End Time _____ Temp. _____ End Time _____ Temp. _____

2nd Incub.: Start Time _____ Temp. _____

End Time _____ Temp. _____