

PGE₂ PLATE CLIA ASSAY LAYOUT SHEET
for use with Enzo Life Sciences Catalog No. ADI-910-001

• **DILUTION TABLE FOR MAKING STANDARDS 1-8:**

Std.	Diluent Vol. (μL)	Vol. Added (μL)	PGE ₂ Conc. (pg/mL)
1	980	20, Stock	1,000
2	500	500, Std. 1	500
3	500	500, Std. 2	250
4	500	500, Std. 3	125
5	500	500, Std. 4	62.5
6	500	500, Std. 5	31.25
7	500	500, Std. 6	15.63
8	500	500, Std. 7	7.81

• **ASSAY PROTOCOL FLOW CHART:**

Well I.D.:	Blank A1, A2	TA A3, A4	NSB A5, A6	Stds. A7 – B10	Bo B11, B12	Samples C1 - H12
Standard Diluent	---	---	100μL	---	100μL	---
Assay Buffer	---	---	50μL	---	---	---
Std. and/or Sample	---	---	---	100μL	---	100μL
Conjugate	---	---	50μL	50μL	50μL	50μL
Antibody	---	---	---	50μL	50μL	50μL
Incub. 2 hours @ RT, shaking	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x 400μL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Conjugate	---	5μL	---	---	---	---
Substrate	200μL	200μL	200μL	200μL	200μL	200μL
Incub. 60 min. @ RT,	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

• PGE₂ PLATE CLIA LAYOUT:

A1 Blank	A2 Blank	A3 TA	A4 TA	A5 NSB	A6 NSB	A7 Std 1	A8 Std 1	A9 Std 2	A10 Std 2	A11 Std 3	A12 Std 3
B1 Std 4	B2 Std 4	B3 Std 5	B4 Std 5	B5 Std 6	B6 Std 6	B7 Std 7	B8 Std 7	B9 Std 8	B10 Std 8	B11 Bo	B12 Bo
C1	C2	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1	D2	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E1	E2	E3	E4	E5	E6	E7	E8	E9	E10	E11	E12
F1	F2	F3	F4	F5	F6	F7	F8	F9	F10	F11	F12
G1	G2	G3	G4	G5	G6	G7	G8	G9	G10	G11	G12
H1	H2	H3	H4	H5	H6	H7	H8	H9	H10	H11	H12

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

1st Incub.: Start Time _____ Temp. _____ Notes: _____

End Time _____ Temp. _____

2nd Incub.: Start Time _____ Temp. _____

End Time _____ Temp. _____