

High Sensitivity PGF_{2α} ASSAY LAYOUT SHEET
for use with Enzo Life Sciences Catalog Nos. ADI-930-069 and ADI-931-069

• **DILUTION TABLE FOR MAKING STANDARDS 1-6:**

Std.	Assay Buffer Vol. (μL)	Vol. Added (μL)	PGF_{2α} Conc. (pg/mL)
Tube A	900	100, Stock	50,000
Tube B	900	100, Tube A	5,000
1	600	400, Tube B	2,000
2	750	250, Std. 1	500
3	750	250, Std. 2	125
4	750	250, Std. 3	31.3
5	750	250, Std. 4	7.8
6	750	250, Std. 5	1.95

• **ASSAY PROTOCOL FLOW CHART:**

Well I.D.:	Blank A1, B1	TA C1, D1	NSB E1, F1	Bo G1, H1	Stds. A2 - D3	Samples E3 - H12
Standard Diluent	---	---	100 μL	100 μL	---	---
Assay Buffer	---	---	25 μL	---	---	---
Std. and/or Sample	---	---	---	---	100 μL	100 μL
Conjugate	---	---	25 μL	25 μL	25 μL	25 μL
Antibody	---	---	---	25 μL	25 μL	25 μL
Incub. 0/N @ 4°C	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x 400 μL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Conjugate	---	2.5μL	---	---	---	---
Substrate	200 μL	200 μL	200 μL	200 μL	200 μL	200 μL
Incub. 3 hr. @ 37°C	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	50 μL	50 μL	50 μL	50 μL	50 μL	50 μL

High Sensitivity PGF_{2α} PLATE LAYOUT:

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