

## 6-keto-PGF<sub>1α</sub> ASSAY LAYOUT SHEET

for use with Enzo Life Sciences Catalog Nos. ADI-900-004 and ADI-901-004  
(formerly 90004 and 90104)

• **DILUTION TABLE FOR MAKING STANDARDS 1-7:**

Std.	Std. Diluent Vol. (μL)	Vol. Added (μL)	6-keto PGF <sub>1α</sub> Conc. (pg/mL)
1	900	100, Stock	50,000
2	800	200, Std. 1	10,000
3	800	200, Std. 2	2,000
4	800	200, Std. 3	400
5	800	200, Std. 4	80
6	800	200, Std. 5	16
7	800	200, Std. 6	3.2

• **ASSAY PROTOCOL FLOW CHART:**

Well I.D.:	Blank A1, B1	TA C1, D1	NSB E1, F1	Bo G1, H1	Stds. A2 - F3	Samples G3 - H12
Standard Diluent (AB or FGF 5)	---	---	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	---	<b>5μL*</b>	---	---	---	---
Substrate	200μL	200μL	200μL	200μL	200μL	200μL
Incub. 45 min. @ RT	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	50μL	50μL	50μL	50μL	50μL	50μL

\* Conjugate for TA must be diluted 1:10 into Assay Buffer. Add 500 μL of Assay Buffer to a clean test tube. From this volume, remove 50 μL of Assay Buffer and, using the same measuring device, add 50 μL of 6-keto-PGF<sub>1α</sub> Conjugate.

• 6-keto PGF<sub>1α</sub> 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 Std 7	E4	E5	E6	E7	E8	E9	E10	E11	E12
F1 NSB	F2 Std 3	F3 Std 7	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. \_\_\_\_\_