

## DIRECT cAMP ASSAY LAYOUT SHEET

for use with Enzo Life Sciences Catalog Nos. ADI-900-066A and ADI-901-066A

• **DILUTION TABLES FOR MAKING STANDARDS 1-5:**

**Non-Acetylated Version:**

Std.	0.1M HCl Vol. (μL)	Vol. Added (μL)	cAMP Conc. (pmol/mL)
1	900	100, Stock	200
2	750	250, Std. 1	50
3	750	250, Std. 2	12.5
4	750	250, Std. 3	3.12
5	750	250, Std. 4	0.78

**Acetylated Version:**

Std.	0.1M HCl Vol. (μL)	Vol. Added (μL)	cAMP Conc. (pmol/mL)
1	990	10, Stock	20
2	750	250, Std. 1	5
3	750	250, Std. 2	1.25
4	750	250, Std. 3	0.312
5	750	250, Std. 4	0.078

• **ACETYLATION PROCEDURE (Optional):**

- 1) Prepare Acetylating Reagent by mixing 0.5 mL of Acetic Anhydride with 1 mL of Triethylamine.
- 2) Add 10 μL of the Acetylating Reagent to 200 μL of standard (including 0 and NSB) or sample. Vortex immediately.
- 3) Assay the acetylated standards and samples within 30 minutes.

• **ASSAY PROTOCOL FLOW CHART:**

	Blank	TA	NSB	Zero Std. (Bo)	Stds.	Samples
Well I.D.:	A1, B1	C1, D1	E1, F1	G1, H1	A2 - B3	C3 - H12
Neutralizing Reagent	---	---	50 μL	50 μL	50 μL	50 μL
0.1M HCl	---	---	150 μL	100 μL	---	---
Std. and/or Sample	---	---	---	---	100 μL	100 μL
Conjugate	---	---	50 μL	50 μL	50 μL	50 μL
Antibody	---	---	---	50 μL	50 μL	50 μL
Incub. 2 hours @ RT, shaking	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x 200μL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Conjugate	---	5 μL	---	---	---	---
Substrate	200 μL	200 μL	200 μL	200 μL	200 μL	200 μL
Incub. 1.5 hours @ RT, shaking	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	50 μL	50 μL	50 μL	50 μL	50 μL	50 μL

• **DIRECT cAMP 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	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1 TA	D2 Std 2	D3	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. \_\_\_\_\_