

**ARG<sup>8</sup>-VASOPRESSIN ASSAY LAYOUT SHEET**  
**for use with Enzo Life Sciences Catalog Nos. ADI-900-017A and ADI-901-017A**

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

<b>Std.</b>	<b>Assay Buffer Vol. (μL)</b>	<b>Vol. Added (μL)</b>	<b>Vasopressin Conc. (pg/mL)</b>
1	900	100, Stock	1,000
2	600	400, Std. 1	400
3	600	400, Std. 2	160
4	600	400, Std. 3	64
5	600	400, Std. 4	25.6
6	600	400, Std. 5	10.24
7	600	400, Std. 6	4.10

• **ASSAY PROTOCOL FLOW CHART:**

<b>Well I.D.:</b>	<b>Blank A1, B1</b>	<b>NSB C1, D1</b>	<b>Bo E1, F1</b>	<b>Stds. G1 - D3</b>	<b>Samples E3 - H12</b>
Assay Buffer	---	150μL	100μL	---	---
Std. and/or Sample	---	---	---	100μL	100μL
Conjugate	---	50μL	50μL	50μL	50μL
Antibody	---	---	50μL	50μL	50μL
Incub. 18-24 hours @ 4°C	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x 400μL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
SA-HRP	---	200μL	200μL	200μL	200μL
Incub. 30 min. @ RT, shaking	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 3 x 400μL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
SA-HRP	---	---	---	---	---
Substrate	200μL	200μL	200μL	200μL	200μL
Incub. 30 min. @ RT, shaking	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	100μL	100μL	100μL	100μL	100μL
Read at 450nm	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

• **VASOPRESSIN PLATE LAYOUT:**

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

2nd Incub.: Start Time \_\_\_\_\_ Temp. \_\_\_\_\_ Notes: \_\_\_\_\_

End Time \_\_\_\_\_ Temp. \_\_\_\_\_ Notes: \_\_\_\_\_

3rd Incub.: Start Time \_\_\_\_\_ Temp. \_\_\_\_\_ Notes: \_\_\_\_\_

End Time \_\_\_\_\_ Temp. \_\_\_\_\_ Notes: \_\_\_\_\_