

**COX-2 (HUMAN) ASSAY LAYOUT SHEET**  
for use with Enzo Life Sciences Catalog No. ADI-900-094

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

\* First add 500µL of DI water to the vial of Standard, let sit for 5 min., then mix gently.

<b>Std.</b>	<b>Assay Buffer Vol. (µL)</b>	<b>Vol. Added (µL)</b>	<b>human COX-II Conc. (ng/mL)</b>
1	230	230, Stock	70
2	230	230, Std. 1	35
3	230	230, Std. 2	17.5
4	230	230, Std. 3	8.75
5	230	230, Std. 4	4.38
6	230	230, Std. 5	2.19
7	230	230, Std. 6	1.09

• **ASSAY PROTOCOL FLOW CHART:**

<b>Well I.D.:</b>	<b>Blank A1, B1</b>	<b>S0 C1, D1</b>	<b>Stds. E1 - B3</b>	<b>Samples C3 - H12</b>
Assay Buffer	100µL	100µL	---	---
Std. and/or Sample	---	---	100µL	100µL
Tap plate gently & seal	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Incub. 1 hour @ 37°C	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 7 x 200µL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Labeled Antibody	---	100µL	100µL	100µL
Seal plate & Incub. 30 min. @ 4°C	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 9 x 200µL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Substrate	100µL	100µL	100µL	100µL
Incub. 30 min. @ RT, dark	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	100µL	100µL	100µL	100µL

• **COX-2 (HUMAN) PLATE LAYOUT:**

A1 Blank	A2 Std 3	A3 Std 7	A4	A5	A6	A7	A8	A9	A10	A11	A12
B1 Blank	B2 Std 3	B3 Std 7	B4	B5	B6	B7	B8	B9	B10	B11	B12
C1 S0	C2 Std 4	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1 S0	D2 Std 4	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E1 Std 1	E2 Std 5	E3	E4	E5	E6	E7	E8	E9	E10	E11	E12
F1 Std 1	F2 Std 5	F3	F4	F5	F6	F7	F8	F9	F10	F11	F12
G1 Std 2	G2 Std 6	G3	G4	G5	G6	G7	G8	G9	G10	G11	G12
H1 Std 2	H2 Std 6	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. \_\_\_\_\_  
 3rd Incub.: Start Time \_\_\_\_\_ Temp. \_\_\_\_\_  
 End Time \_\_\_\_\_ Temp. \_\_\_\_\_