

IgG₁ (mouse) ELISA kit, Assay Layout Sheet
for use with Enzo Life Sciences Catalog No. ADI-900-109A

• **TABLE FOR MAKING STANDARDS 1-6:**

Std.	Diluent Vol. (μL)	Vol. Added (μL)	mouse IgG₁ Conc. (ng/mL)
1	475	25 Stock	250
2	250	250 Std. 1	125
3	250	250 Std. 2	62.5
4	250	250 Std. 3	31.25
5	250	250 Std. 4	15.62
6	250	250 Std. 5	7.81

• **ASSAY PROTOCOL FLOW CHART:**

Well ID:	Blank A1, B1	S0 C1, D1	Standards E1-H2	Samples A3-H12
Standard Diluent	---	50μL	---	---
Std. and/or Sample	---	---	50μL	50μL
Conjugate	---	50μL	50μL	50μL
Incub. 1 hour @ RT, shaking, sealed	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒
Asp. & Wash 4 x 400μL	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒
Substrate	100μL	100μL	100μL	100μL
Incub. 30 min. @ RT, shaking	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒	⇒⇒⇒
Stop Solution 2	100μL	100μL	100μL	100μL

IgG₁ (mouse) ELISA, Plate Layout:

A1 Blank	A2 Std 3	A3	A4	A5	A6	A7	A8	A9	A10	A11	A12
B1 Blank	B2 Std 3	B3	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. _____