

LVV Hemorphin 7 ASSAY LAYOUT SHEET
for use with Enzo Life Sciences Catalog No. ADI-900-205

• **TABLE FOR MAKING STANDARDS 1-7:**

Reconstitute lyophilized standard with 1mL Assay Buffer, vortex gently to dissolve. Then proceed to making standards.

Std	Assay Buffer Vol. (µL)	Standard Vol. Added (µL)	LVV Hemorphin 7 Conc. (pg/mL)
1	750	250, Stock	10000
2	750	250, Std. 1	2500
3	750	250, Std. 2	625
4	750	250, Std. 3	156
5	750	250, Std. 4	39.1
6	750	250, Std. 5	9.8

• **ASSAY PROTOCOL FLOW CHART Vol. Added (µL):**

Well I.D.:	Blank A1, B1	NSB C1, D1	Bo A3, B3	Stds E1-B3	Samples C3-H12
Assay Buffer	---	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 hrs. @ RT, shaking (sealed)	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 4 x 400 µL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
SA-HRP	---	200 µL	200 µL	200 µL	200 µL
Incub. 30 mins. @ RT, shaking (sealed)	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Asp. & Wash 4 x 400 µL	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
TMB Substrate	200 µL	200 µL	200 µL	200 µL	200 µL
Incub. 30min @ RT	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	50 µL	50 µL	50 µL	50 µL	50 µL
read @ 450 nm	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

LVV Hemorphin 7 Plate Layout:

A1 Blank	A2 Std 3	A3 S0	A4	A5	A6	A7	A8	A9	A10	A11	A12
B1 Blank	B2 Std 3	B3 S0	B4	B5	B6	B7	B8	B9	B10	B11	B12
C1 NSB	C2 Std 4	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1 NSB	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. _____
 End Time _____ Temp. _____

2nd Incub.: Start Time _____ Temp. _____
 End Time _____ Temp. _____

3rd Incub.: Start Time _____ Temp. _____
 End Time _____ Temp. _____