

15(S)-HETE ASSAY LAYOUT SHEET

for use with Enzo Life Sciences Catalog Nos. ADI-900-051 and ADI-901-051

• **TABLE FOR MAKING STANDARDS 1-5:**

Std.	Assay Buffer Volume (μL)	Vol. added (μL)	15(S)-HETE Conc. (pg/mL)
1	900	100 μL Stock	20,000
2	750	250 μL Std. 1	5,000
3	750	250 μL Std. 2	1250
4	750	250 μL Std. 3	312.5
5	750	250 μL Std. 4	78.12

• **ASSAY PROTOCOL FLOW CHART:**

Well ID:	Blank A1, B1	TA C1, D1	NSB E1, F1	Bo G1, H1	Standards A2-B3	Samples C3-H12
Standard Diluent	---	---	100μL	100μL	---	---
Assay Buffer	---	---	50μL	---	---	---
Std and/or Sample	---	---	---	---	100μL	100μL
Conjugate	---	---	50μL	50μL	50μL	50μL
Antibody	---	---	---	50μL	50μL	50μL
Tap plate gently to mix	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Incub. 18-24 hours @ 4°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Wash 3x w/400 μL wash buffer	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Conjugate	---	5μL	---	---	---	---
Substrate	200μL	200μL	200μL	200μL	200μL	200μL
Tap plate gently to mix	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Incubate 1 hour @ 37°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Stop Solution	50μL	50μL	50μL	50μL	50μL	50μL
Read @ 405 nm	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

15(S)-HETE 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. _____