

SOD ASSAY LAYOUT SHEET
for use with Enzo Life Sciences Catalog No. ADI-900-157

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

Std.	1X SOD Buffer Vol. (µL)	Vol. Added (µL)	SOD Conc. (units/25 µL)
1	620	5, Stock	10
2	100	100, Std. 1	5.0
3	150	100, Std. 2	2.0
4	100	100, Std. 3	1.0
5	100	100, Std. 4	0.5
6	150	100, Std. 5	0.2
7	100	100, Std. 6	0.1

• **ASSAY PROTOCOL FLOW CHART:**

Well I.D.:	Activity Control A1-A3	Stds. B1 - H3	Samples A4 - H12
1X SOD Buffer	25 µL	---	---
Std. and/or Sample	---	25 µL	25 µL
Master Mix	150 µL	150 µL	150 µL
1X Xanthine Solution	25 µL	25 µL	25 µL
Read at 450 nm once a min. for 10 min.	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

• SOD DISMUTASE PLATE LAYOUT:

A1 No SOD Activity Control	A2 No SOD Activity Control	A3 No SOD Activity Control	A4	A5	A6	A7	A8	A9	A10	A11	A12
B1 0.1 units SOD	B2 0.1 units SOD	B3 0.1 units SOD	B4	B5	B6	B7	B8	B9	B10	B11	B12
C1 0.2 units SOD	C2 0.2 units SOD	C3 0.2 units SOD	C4	C5	C6	C7	C8	C9	C10	C11	C12
D1 0.5 units SOD	D2 0.5 units SOD	D3 0.5 units SOD	D4	D5	D6	D7	D8	D9	D10	D11	D12
E1 1.0 units SOD	E2 1.0 units SOD	E3 1.0 units SOD	E4	E5	E6	E7	E8	E9	E10	E11	E12
F1 2.0 units SOD	F2 2.0 units SOD	F3 2.0 units SOD	F4	F5	F6	F7	F8	F9	F10	F11	F12
G1 5.0 units SOD	G2 5.0 units SOD	G3 5.0 units SOD	G4	G5	G6	G7	G8	G9	G10	G11	G12
H1 10 units SOD	H2 10 units SOD	H3 10 units SOD	H4	H5	H6	H7	H8	H9	H10	H11	H12

Kit Lot No. _____ Exp. Date _____ Date _____ Tech. _____

Notes: _____

