

Acetyltransferase ASSAY LAYOUT SHEET

for use with Enzo Life Sciences Catalog No. ADI-907-026

AT End Point Assay Protocol Flow Chart:

Well ID:	Blank A1, A2	Enzyme Dilutions A3-A12
Mix all reagents gently	---	---
Transferase Assay Buffer	25 μ L	---
Acetyltransferase Dilution	---	25 μ L
Reaction Mix	25 μ L	25 μ L
Incubate 30 min. @ 25°C, shaking, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒
Isopropyl Alcohol	50 μ L	50 μ L
Detection Solution	100 μ L	100 μ L
Incubate 10 min. @ 25°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒
Read Fluorescence at 380 _{ex} /520 _{em}	⇒⇒⇒⇒	⇒⇒⇒⇒

AT Control Protocol Flow Chart:

Well ID:	Blank B1, B2	Positive Control B3, B4
Mix all reagents gently	---	---
Transferase Assay Buffer	50 μ L	---
AT Positive Control	---	50 μ L
Isopropyl Alcohol	50 μ L	50 μ L
Detection Solution	100 μ L	100 μ L
Incubate 10 min. @ 25°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒
Read Fluorescence at 380 _{ex} /520 _{em}	⇒⇒⇒⇒	⇒⇒⇒⇒

AT Inhibition Assay Protocol Flow Chart:

Well ID:	Blank C1, C2	Inhibition Samples C3-C10	D0 C11, C12
Mix all reagents gently	---	---	---
Transferase Assay Buffer	25 µL	---	10 µL
Inhibitor Dilution	---	10 µL	---
Acetyltransferase	---	15 µL	15 µL
Incubate 10 min. @ 25°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Reaction Mix	25 µL	25 µL	25 µL
Incubate 30 min. @ 25°C, shaking, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Isopropyl Alcohol	50 µL	50 µL	50 µL
Detection Solution	100 µL	100 µL	100 µL
Incubate 10 min. @ 25°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒
Read Fluorescence at 380 _{ex} /520 _{em}	⇒⇒⇒⇒	⇒⇒⇒⇒	⇒⇒⇒⇒

AT Kinetic Assay Protocol Flow Chart:

Well ID:	Blank D1, D2	Enzyme Dilutions E1-H12
Mix all reagents gently	---	---
Transferase Assay Buffer	25 µL	---
Acetyltransferase Dilution	---	25 µL
Reaction Mix	25 µL	25 µL
Incubate 30 min. @ 25°C, shaking, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒
Isopropyl Alcohol [Stop replicate dilutions at predetermined time intervals.]	50 µL	50 µL
Detection Solution	100 µL	100 µL
Incubate 10 min. @ 25°C, sealed	⇒⇒⇒⇒	⇒⇒⇒⇒
Read Fluorescence at 380 _{ex} /520 _{em}	⇒⇒⇒⇒	⇒⇒⇒⇒

Acetyltransferase PLATE LAYOUT:

A1 Blank	A2 Blank	A3 Dil.1	A4 Dil.1	A5 Dil.2	A6 Dil.2	A7 Dil.3	A8 Dil.3	A9 Dil.4	A10 Dil.4	A11 Dil.5	A12 Dil.5
B1 Blank	B2 Blank	B3 Pos. Control	B4 Pos. Control	B5	B6	B7	B8	B9	B10	B11	B12
C1 Blank	C2 Blank	C3 Inhib. Dil.1	C4 Inhib. Dil.1	C5 Inhib. Dil.2	C6 Inhib. Dil.2	C7 Inhib. Dil.3	C8 Inhib. Dil.3	C9 Inhib. Dil.4	C10 Inhib. Dil.4	C11 D0	C12 D0
D1 Kinetic Blank	D2 Kinetic Blank	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E1 Kinetic D1	E2 Kinetic D1	E3 Kinetic D1	E4 Kinetic D1	E5 Kinetic D1	E6 Kinetic D1	E7 Kinetic D1	E8 Kinetic D1	E9 Kinetic D1	E10 Kinetic D1	E11 Kinetic D1	E12 Kinetic D1
F1 Kinetic D2	F2 Kinetic D2	F3 Kinetic D2	F4 Kinetic D2	F5 Kinetic D2	F6 Kinetic D2	F7 Kinetic D2	F8 Kinetic D2	F9 Kinetic D2	F10 Kinetic D2	F11 Kinetic D2	F12 Kinetic D2
G1 Kinetic D3	G2 Kinetic D3	G3 Kinetic D3	G4 Kinetic D3	G5 Kinetic D3	G6 Kinetic D3	G7 Kinetic D3	G8 Kinetic D3	G9 Kinetic D3	G10 Kinetic D3	G11 Kinetic D3	G12 Kinetic D3
H1 Kinetic D4	H2 Kinetic D4	H3 Kinetic D4	H4 Kinetic D4	H5 Kinetic D4	H6 Kinetic D4	H7 Kinetic D4	H8 Kinetic D4	H9 Kinetic D4	H10 Kinetic D4	H11 Kinetic D4	H12 Kinetic D4