



# CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04356

**Product Description:** MSD® Total Akt Kit  
MSD® Total Akt Base Kit

**Catalog #/Size:** K150MOD-1 / 1 Plate; K150MOD-2 / 5 Plates; K150MOD-4 / 25 Plates  
K150MOA-4 / 25 Plates

**Lot Number:** K0013524

**Expiration Date:** 30 JUN 2016

### Acceptance Data:

Metric		Specification	Assay Total Akt
% Recovery Concentration	Calibrator	80% - 120%	Pass
	Control	70% - 130%	Pass
% Recovery CV	Calibrator	< 20%	Pass
	Control	< 20%	Pass

### Concentration of Materials Tested:

	Assay (pg/well) Total Akt
Calibrator Concentration (A0010342)	87 800
Control 1 Concentration (A00C0176)	485
Control 2 Concentration (A00C0177)	21.6

### Standard Curve Data from Kit QC:

Total Akt		
Conc., (pg/well)	Average Signal	%CV
1097	868 930	4.9
274	271 696	1.7
68.6	69 647	2.1
17.1	17 278	3.5
4.29	4413	4.4
1.07	1296	4.4
0.268	481	7.9
0	199	10.1



## CERTIFICATE OF ANALYSIS


Page 2 of 2

COA-04356

Total Akt Kit Components (K150MOD-1, K150MOD-2, K150MOD-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054865	2-8°C	30 JUN 2016
Total Akt Antibody	D0013084	2-8°C	30 SEP 2016
Akt Calibrator	A0010342	≤ -70°C	30 SEP 2016
Diluent 39	Not Kit Specific	2-8°C	N/A
Protease Inhibitor	Not Kit Specific	2-8°C	N/A
Phosphatase Inhibitor I	Not Kit Specific	2-8°C	N/A
Phosphatase Inhibitor II	Not Kit Specific	2-8°C	N/A
MSD Tris Wash Buffer (10X), 200mL	Not Kit Specific	2-8°C	N/A
MSD Blocker A	Not Kit Specific	Room Temperature	N/A
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Total Akt Base Kit Components (K150MOA-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054865	2-8°C	30 JUN 2016
Total Akt Antibody	D0013084	2-8°C	30 SEP 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

**Statement:** The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dana Short	Quality Control		18 APR 2014