



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA- 04731

Product Description: Phospho-Akt (Ser473) Kit

Phospho-Akt (Ser473) Base Kit

Catalog #/Size: K150MND-1 / 1 Plate; K150MND-2 / 5 Plates; K150MND-4 / 25 Plates

K150MNA-4 / 25 Plates

Kit Lot Number: K0013627

Expiration Date: 31 OCT 2016

Acceptance Data:

Metric		Specification	Assay
			Phospho-Akt (Ser473)
% Recovery Concentration	Calibrator	80% - 120%	Pass
	Control	70% - 130%	Pass
% Recovery CV	Calibrator	< 20%	Pass
	Control	< 20%	Pass

Concentration of Materials Tested:

	Assay (U _{P473} /well)
	Phospho-Akt (Ser473)
20X Calibrator Concentration (A0010357)	872
Control 1 Concentration (A00C0176)	1.80
Control 2 Concentration (A00C0177)	0.112

Standard Curve Data from Kit QC:

Phospho-Akt (Ser473)		
Conc., (U _{P473} /well)	Average Signal	%CV
43.6	474 619	3.9
10.9	160 678	2.7
2.73	31 109	2.1
0.681	5732	1.8
0.170	1354	2.6
0.0426	400	4.0
0.0106	180	3.6
0	106	7.9



CERTIFICATE OF ANALYSIS


Page 2 of 2

COA- 04731

Phospho-Akt (Ser473) Kit Components (K150MND-1, K150MND-2, K150MND-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0055041	2-8° C	30 NOV 2016
Phospho-Akt (Ser473) Antibody	D0013197	2-8° C	31 JAN 2017
Akt Calibrator	A0010357	≤ -70° C	31 OCT 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)	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

Phospho-Akt (Ser473) Base Kit Components (K150MNA-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0055041	2-8° C	30 NOV 2016
Phospho-Akt (Ser473) Antibody	D0013197	2-8° C	31 JAN 2017
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		22 SEP 2014