



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-03302

Product Description: Phospho-Akt (Ser473) Kit

Kit Catalog Numbers: K150MND-1, K150MND-2, K150MND-4, K150MNA-4

Kit Lot Number: K0013050

Expiration Date: 24 Aug 2014

Acceptance Data:

Metric		Specifications	Assay Phospho Akt (Ser473)
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	75% - 125%	Pass
% Recovery CV	Calibrators	<20%	Pass
	Controls	<25%	Pass

Concentration of Materials Tested:

Units = U_{P473}/well	Assay Phospho Akt (Ser473)
Calibrator Concentration	906
Control 1 (A00C0143)	5.30
Control 2 (A00C0144)	0.200

Standard Curve Data from Kit QC:

Phospho Akt (Ser473)		
Conc., (U_{P473}/well)	Average Signal	%CV
45.3	359 340	6.1
11.3	121 421	4.8
2.83	24 315	5.2
0.707	4 824	7.7
0.177	1 181	7.4
0.0441	340	7.5
0.0111	130	7.7
0	70.1	12.0



CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-03302

Phospho-Akt (Ser473) Kit Components (K150MND-1, K150MND-2, K150MND-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054167	2-8° C	30 Jun 2015
SULFO-TAG Anti-Phospho-Akt (Ser473) Antibody	D0012597	2-8° C	30 Jun 2015
Phospho-Akt (Ser473) Calibrator *	A0010299	≤ -70° C	31 Aug 2014
Diluent 39	Not Kit Specific	2-8° C	N/A
Protease Inhibitor Solution	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
Tris Wash Buffer (10X)	Not Kit Specific	2-8° C	N/A
Blocker A	Not Kit Specific	Room Temperature	N/A
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

* The Phospho-Akt (Ser473) Calibrator and Akt Calibrator are equivalent products. They are both full length recombinant Akt1 protein expressed in baculovirus and phosphorylated in vitro.

Phospho-Akt (Ser473) Base Kit Components (K150MNA-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054167	2-8° C	30 Jun 2015
SULFO-TAG Anti-Phospho-Akt (Ser473) Antibody	D0012597	2-8° C	30 Jun 2015
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	Paula Eason	Quality Control		21 Mar 2013