



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-03301

Product Description: Total Akt Kit

Kit Catalog Numbers: K150MOD-1, K150MOD-2, K150MOD-4, K150MOA-4

Kit Lot Number: K0013054

Expiration Date: 19 Oct 2014

Acceptance Data:

Metric		Specifications	Assay
			Total Akt
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	75% - 125%	Pass
% Recovery CV	Calibrators	<20%	Pass
	Controls	<25%	Pass

Concentration of Materials Tested:

Units = pg/well	Assay
	Total Akt
Calibrator Concentration	86 200
Control 1 (A00C0143)	420
Control 2 (A00C0144)	25

Standard Curve Data from Kit QC:

Total Akt		
Conc., (pg/well)	Average Signal	%CV
1080	801 678	3.9
269	247 009	5.1
67.3	59 181	6.0
16.8	13 854	7.8
4.21	3 490	9.2
1.05	939	9.9
0.263	297	11.0
0	84.6	9.5



CERTIFICATE OF ANALYSIS


Page 2 of 2

COA-03301

Total Akt Kit Components (K150MOD-1, K150MOD-2, K150MOD-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054173	2-8° C	30 Jun 2015
SULFO-TAG Anti-Total Akt Antibody	D0012601	2-8° C	30 Jun 2015
Akt Calibrator	A0010306	≤ -70° C	31 Oct 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

Total Akt Base Kit Components (K150MOA-4)			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0054173	2-8° C	30 Jun 2015
SULFO-TAG Anti-Total Akt Antibody	D0012601	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