



CERTIFICATE OF ANALYSIS

COA – 3545

Product Description: Akt Calibrator
Catalog #/Size: C00MN-2
Lot Number: A0010306
Expiration Date: 31 Oct 2014
Storage Temperature: $\leq -70^{\circ}$ C

Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U _{P473} /well)	Phospho-Akt (Ser473) (U _{P473} /mL)
Calibrator Concentration	862	34 500

	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	86 200	3 450

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

	Name	Function	Signature	Date
Review/Approval	Paula Eason	Quality Control		20 Nov 2013