



CERTIFICATE OF ANALYSIS

COA – 04177

Product Description: Akt Calibrator

Catalog Number: C00MN-2

Lot Number: A0010341

Expiration Date: 31 JUL 2016


Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U _{P473} /well)	Phospho-Akt (Ser473) (U _{P473} /mL)
Calibrator Concentration	885	35 400

	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	88 500	3 540

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		21 FEB 2014