



CERTIFICATE OF ANALYSIS

COA – 04221

Product Description: Akt Calibrator

Catalog Number: C00MN-2

Lot Number: A0010342

Expiration Date: 30 SEP 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	878	35 100

	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	87 800	3 510

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	Dana Short	Quality Control		04 APR 2014