



## CERTIFICATE OF ANALYSIS

COA – 3544

**Product Description:** Phospho-Akt (Ser473) Calibrator

**Catalog Number:** C00MN-2

**Lot Number:** A0010299

**Expiration Date:** 31 Aug 2014

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

### Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U <sub>P473</sub> /well)	Phospho-Akt (Ser473) (U <sub>P473</sub> /mL)
Calibrator Concentration	906	36 200

	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	90 600	3 620

\* 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.

**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