



CERTIFICATE OF ANALYSIS

COA - 06629

Product Description: MSD[®] Insulin Signaling Panel (Phospho Protein) Kit

Catalog #/Size: K15151C-1 / 1 Plate; K15151C-2 / 5 Plates; K15151C-3 / 20 Plates

Lot Number: K0013868

Expiration Date: N/A (See expiration date on specific components below)

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Kit Contents:

Plate Lot Number: Z0055466

Plate Lot Catalog Number: N45151A

Plate Lot Expiration Date: 31 MAR 2018

Antibody blend Lot Number: D0013461

Antibody blend Catalog Number: D20HP

Antibody blend Expiration Date: 28 FEB 2018

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

No more than one plate with CV $> 18\%$

No plate with CV $> 20\%$

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		14 MAR 2016
Review/Approval	Dawn Stollar	Quality Assurance		14 MAR 2016