



CERTIFICATE OF ANALYSIS

COA - 06371

Product Description: MSD[®] Total STAT5a,b Whole Cell Lysate Kit

Catalog #/Size: K150IHD-1 / 1 Plate; K150IHD-2 / 5 Plates; K150IHD-3 / 20 Plates

Lot Number: K0013813

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: Z0055326

Plate Lot Catalog Number: N450IHA

Plate Lot Expiration Date: 30 SEP 2017

Antibody blend Lot Number: D0013399

Antibody blend Catalog Number: D20IH

Antibody blend Expiration Date: 30 SEP 2017

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		06 JAN 2016
Review/Approval	Dawn Stollar	Quality Assurance		06 JAN 2016