



CERTIFICATE OF ANALYSIS

COA - 06370

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

Catalog #/Size: K150IGD-1 / 1 Plate; K150IGD-2 / 5 Plates; K150IGD-3 / 20 Plates

Lot Number: K0013678

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: Z0055117

Plate Lot Catalog Number: N450IGA

Plate Lot Expiration Date: 31 JAN 2017

Antibody blend Lot Number: D0013251

Antibody blend Catalog Number: D20IG

Antibody blend Expiration Date: 31 JAN 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 |