



CERTIFICATE OF ANALYSIS

COA - 07124

Product Description: MSD[®] Phospho-STAT3 Whole Cell Lysate Kit

Catalog #/Size: K150DID-1 / 1 Plate; K150DID-2 / 5 Plates; K150DID-3 / 20 Plates

Lot Number: K0013734

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: Z0055273

Plate Lot Catalog Number: N450DIA

Plate Lot Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013320

Antibody blend Catalog Number: D21DI

Antibody blend Expiration Date: 30 APR 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		15 JUL 2016
Review/Approval	Vanessa Castleberry	Quality Assurance		15 JUL 2016