



CERTIFICATE OF ANALYSIS

COA - 07577

Product Description: MSD[®] Total STAT3 Kit

Catalog #/Size: K150SND-1 / 1 Plate; K150SND-2 / 5 Plates; K150SND-3 / 20 Plates

Lot Number: K0013933

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: Z0055589

Plate Lot Catalog Number: N450SMA

Plate Lot Expiration Date: 30 SEP 2018

Antibody blend Lot Number: D0013535

Antibody blend Catalog Number: D20SN

Antibody blend Expiration Date: 31 AUG 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	Xiomara Talavera	Quality Control	<i>X Talavera</i>	02 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	12 DEC 2016