



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07218

Product Description: U-PLEX[®] 5-Assay, 96-Well SECTOR[®] Plate

Catalog Number: N05230A-1

Lot Number: Z00U0023D

Expiration Date: 31 JAN 2019

Storage Condition: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	$\leq 10.0\%$	Pass
Average Interplate CV	$\leq 13.0\%$	Pass
Signal Range at Binding Capacity	$\leq 30.0\%$	Pass

Statement: The above plate lot was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	12Aug2016