



CERTIFICATE OF ANALYSIS

COA - 04528

Product Description: MSD[®] JAK-STAT Whole Cell Lysate Set

Catalog #/Size: C10KQ-1



Lot Number: CL0010003 / CL0010004

Expiration Date: 30 NOV 2017

Statement: This lysate set was evaluated by MSD assay and has passed specifications.

Concentration: 2mg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biooigcs/Chemistry		14 AUG 2014
Review/Approval	Dana Short	Quality Control		14 AUG 2014