





**CERTIFICATE OF ANALYSIS**

COA - 03505

**Product Description:** MSD<sup>®</sup> JAK-STAT Whole Cell Lysate Set  
**Catalog #/Size:** C10KQ-1  
**Lot Number:** RLL260262/261263  
**Expiration Date:** 31 MAR 2016  
**Statement:** This lysate set was evaluated by MSD assay and passed specifications.

Concentration: 2mg/mL

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

|                        | <b>Name</b>      | <b>Function</b>    | <b>Signature</b>   | <b>Date</b> |
|------------------------|------------------|--------------------|--|-------------|
| <b>Preparer</b>        | Rolanda Spratley | Biooigcs/Chemistry |   | 14 AUG 2014 |
| <b>Review/Approval</b> | Dana Short       | Quality Control    |  | 14 AUG 2014 |