



**CERTIFICATE OF ANALYSIS**

COA - 04874

**Product Description:** MSD<sup>®</sup> MAP Kinase Whole Cell Lysate Set

**Catalog #/Size:** C1101-1



**Lot Number:** CL0010019 / CL0010020

**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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		22 OCT 2014
<b>Review/Approval</b>	Dana Short	Quality Control		22 OCT 2014