



## CERTIFICATE OF ANALYSIS

COA - 07121

**Product Description:** Phosphatase Inhibitor II

**Catalog #/Size:** 3-3057-020009 / 2.0 mL; 3-3057-020008 / 0.5 mL; 3-3057-020002 / 0.1 mL

**Lot Number:** R0010559

**Expiration Date:** 31 OCT 2019



**Statement:** This product was manufactured according to approved MSD documents.

Concentration/Specific Activity: 100X

Storage buffer: N/A

Storage conditions: 2-8°C

**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		15 JUL 2016
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance		15 JUL 2016