



CERTIFICATE OF ANALYSIS

COA - 05197

Product Description: Phosphatase Inhibitor II

Catalog #/Size: 3-3057-020009 / 2 mL

Lot Number: R0010498

Expiration Date: 31 MAY 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.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		10 FEB 2015
Review/Approval	Anastasia Belilovets	Quality Assurance		10 FEB 2015