



CERTIFICATE OF ANALYSIS

Product Description: Phosphatase Inhibitor II

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

Lot Number: R0010591

Expiration Date: 30 APR 2021

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

Concentration/Specific Activity: 100X

Volume: 0.1 mL / 0.5 mL / 2.0 mL

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	23 JAN 2019