



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Phosphatase Inhibitor I

Catalog Number: 3-3057-020001-C / 3-3057-020006-B / 3-3057-020007-B

Lot Number: R0010624

Expiration Date: 31 MAR 2020

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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	09 JUL 2019