

CERTIFICATE OF ANALYSIS

BCIP/NBT 1-Component AP Membrane Substrate (SUBB)			
Catalog#	Size	Lot#	Expiry
3330	Bulk	19R52	09/14/2021
6281	100 mL	19R53	09/14/2021

Specifications	Criteria	Lot Data	Results
pH	9.80 ± 0.20	9.93	Pass
Appearance	Colorless to very light yellow	Conforms	Pass
Storage	2-25°C	2-25°C	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications.
 This information is believed to be accurate and is offered in good faith.
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance: 
 Tracy Murphy

Date: 3/27/19
 03/27/2019