

CERTIFICATE OF ANALYSIS

pNPP 1-Component AP Microwell Substrate (SUBP)			
Catalog#	Size	Lot#	Expiry
3328	Bulk	16N6	08/31/2018
6279	100 mL	16N7	08/31/2018

Specifications	Criteria	Lot Data	Results
pH	9.90 ± 0.30	9.87	Pass
Absorbance at 405 nm	≤ 0.150	0.068	Pass
Appearance	Colorless to very light yellow solution with charcoal pellets	Conforms	Pass
Storage	Refrigerated	Refrigerated	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications.

Signature: *Traug Murphy*
 Quality Assurance

Date: 06/03/2016
 06/03/16