

CERTIFICATE OF ANALYSIS

Antigen-Down Assay Diluent (AD4)

Catalog#	Size	Lot#	Expiry at 2-8°C
523	Bulk	18B37	11/2020
629	100 mL	NA	NA
630	500 mL	NA	NA
631	1 L	18B38	11/2020
674	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.2-7.6	7.40	Pass
Appearance	Clear to slightly yellow liquid	Conforms	Pass
Conductivity μ MHOs	30,000-55,000	33,600	Pass
A ₂₈₀ of 1:20 dilution	0.28-0.42	0.365	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control


Kristi Strandberg

Date

11/15/18
11/15/2018