



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD Free TAG 300 000 ECL

**Catalog #/Size:** 3-3026-420003-B / 250 mL

**Lot Number:** Y0250026

**Expiration Date:** 29 FEB 2020

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

pH:  $7.8 \pm 0.1$

Signal test

ECL (tolerance range): 270 000 – 400 000

CV%:  $\leq 5\%$

**The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 JAN 2019