



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** SULFO-TAG™ Streptavidin

**Catalog Number:** R32AD-Series

**Lot Number:** W0021617S

**Expiration Date:** 31 AUG 2026

**Storage Temperature:** 2-8°C

| Size   | Total Volume | Concentration | L/P | Final Buffer         |
|--------|--------------|---------------|-----|----------------------|
| 50 µg  | 100 µL       | 500 µg/mL     | 9.4 | Diluent 100 w/ Azide |
| 200 µg | 400 µL       | 500 µg/mL     | 9.4 | Diluent 100 w/ Azide |

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

|                 | Name          | Function | Signature                                                                          | Date        |
|-----------------|---------------|----------|------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 14 APR 2023 |