



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG Goat Anti-Rabbit Antibody
Catalog Number: R32AB-1
Lot Number: W0022942S
Expiration Date: 31 JUL 2029
Storage Temperature: 2-8°C

| Size | Total Volume | Concentration | L/P | Final Buffer |
|-------|--------------|---------------|-----|----------------------|
| 200µg | 400 µL | 500 µg/mL | 5.9 | 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 | | 17 APR 2026 |