

# CERTIFICATE of ANALYSIS



## Nitrotyrosine (Peptide) Polyclonal Antibody

Item No. 10006778 • Batch No. 0562893 • Expires 04DEC2026

### Overview

|                            |                                                                                                               |
|----------------------------|---------------------------------------------------------------------------------------------------------------|
| <b>Formulation:</b>        | 500µl peptide affinity purified polyclonal antibody in PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide |
| <b>Concentration:</b>      | 0.2 mg/ml                                                                                                     |
| <b>Storage/Stability:</b>  | -20°C/1yr                                                                                                     |
| <b>Applications:</b>       | WB                                                                                                            |
| <b>Species reactivity:</b> | (Independent)Species                                                                                          |

Reviewed and approved by: Paul Domanski

Approved date: 05JUN2019

A handwritten signature in black ink, appearing to read "Paul Domanski".

**WARNING**  
THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2018

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM