

Certificate of Analysis

Anti-Human ISG15 (MAb)

Catalog No: 21900-1

Lot No:

Size: 100 µg/vial

Description: Mouse Monoclonal Antibody against Human ISG15

Clone: 2.1

Volume: ml

Concentration: mg/ml

Buffer: 50mM Tris HCl, 75mM NaCl, 50% Glycerol pH 8

Antigen: Human ISG15

Isotype: Mouse IgG₁ kappa

Endotoxin: < 1 EU/µg

Purity: > 95%

Purification Method: Protein G purified

Tested Applications: Western blot

Suggested concentration: In our hands good visualization of ISGylated proteins was obtained using an antibody concentration of 1 µg/ml; however, 0.1 µg/ml or lower can be used to give reasonable visualization of the free ISG band. *Optimal dilutions should be determined by each laboratory for each application.*

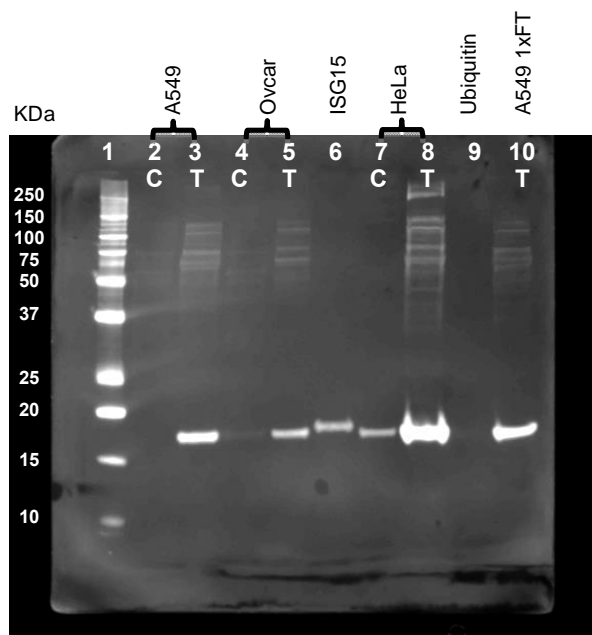


Figure 1. Western Blot. Control (C) and Treated (T) 40 µg samples from A549, Ovarcar and HeLa cell lysates were run alongside 50 ng recombinant ISG15 and 100 µg ubiquitin under reducing conditions in a 12% SDS-PAGE gel. Cells were incubated for 48 hours with or without 1000 U/ml of Human IFN beta 1a (PBL cat# 11410-2). No cross-reactivity was observed with ubiquitin at tested concentrations up to 100 µg. The antibody was additionally tested on Daudi and U937 cell lysates. Please contact us for further information.



References:

1. N Arnaud, S Dabo, D Akazawa, M Fukasawa, F Shinkai-Ouchi, J Hugon, T Wakita, and EF Meurs. **Hepatitis C Virus Reveals a Novel Early Control in Acute Immune Response,***+PLoS Pathog*, 2011, 7(10): e1002289.
2. C Dieterich and DA Relman. **Modulation of the Host Interferon Response and ISGylation Pathway by B. pertussis Filamentous Hemagglutinin,***+PLoS ONE*, 2011, 6(11): e27535.
3. YB Zhang, YL Wang, and JF Gui. **Identification and characterization of two homologues of interferon-stimulated gene ISG15 in crucian carp,***+Fish and Shellfish Immunology*, 2007, 23(1): 52-61.
4. T Takeuchi, T Kobayashi, S Tamura, and H Yokosawa. **Negative regulation of protein phosphatase 2C by ISG15 conjugation,***+FEBS Letters*, 2006, 580(18): 4521-4526.
5. MP Malakhov, K Il Kim, OA Malakhova, BS Jacobs, EC Borden, and D Zhang. **High-throughput Immunoblotting: Ubiquitin-like Protein ISG15 Modifies Key Regulators of Signal Transduction,***+JBC*, 2003, 278, 16608-16613.
6. J D Cunha, S Ramanujam, RJ Wagner, PL Witt, E Knight, Jr and EC Borden. **In vitro and in vivo secretion of human ISG15, an IFN-induced immunomodulatory cytokine,***+Journal of Immunology*, 1996, 157:4100.
7. J D Cunha, E Knight, Jr, AL Hass, RL Trutt and EC Borden. **Immunoregulatory properties of ISG15, an interferon-induced cytokine,***+PNAS*, 1996, 93:211.

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: After receipt, this product should be kept at -20°C for retention of full activity. Do not store below -20°C. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.