

Anti-Human ISG15, MAb

Mouse Monoclonal Antibody against Human ISG15, Clone 2.1
Catalog No. 21900-1

- Detects conjugated and free ISG15
- Tested applications include Western Blot
- Greater than 95% purity
- No cross-reactivity observed with ubiquitin

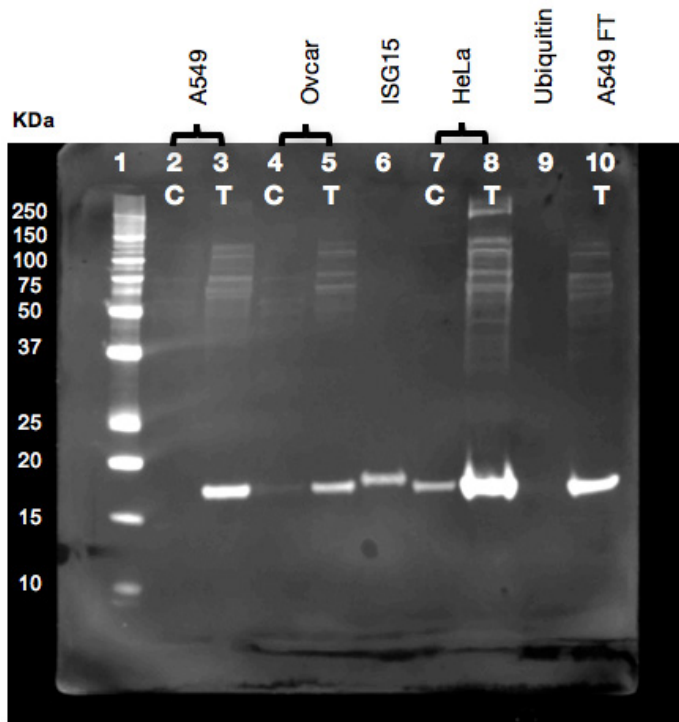


Figure 1. Western Blot. Control (C) and Treated (T) 40 µg samples from A549, Ovarian carcinoma (Ovcar) and HeLa cell lysates were run alongside 50 ng recombinant ISG15 and 100 µg ubiquitin under reducing conditions in a 12% SDS-PAGE gel. Cells were incubated for 48 hours with or without 1000 U/ml of Human IFN beta 1a (PBL Cat No. 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.

Many published reports indicate that ISG15 (G1P2) is commonly upregulated in response to Type I interferon stimulation. Though a wide range of cellular activities are influenced by ISG15 expression including Type I interferon signaling, translation, chromatin remodeling, cell motility, protein trafficking, and protein conjugation (ISGylation), the complete spectrum of ISG15-dependent biological sequelae remains to be fully elucidated. Further definition of new ISG15 target proteins and the role of free ISG15 will offer new insights into mechanistic and immunotherapeutic approaches to human diseases.

Research Citations and General 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 β 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.