



Certificate of Analysis

DIY- Capture Antibody

Catalog No: 60127-1

Lot No: 6176

Expiration: February 1, 2018

Size: 180 ug/ml

Description: Mouse Anti-Human IL-29/IL-28B

Formulation: Lyophilized

Reconstitution: Reconstitute with 1 ml of Phosphate Buffered Saline (PBS) to a concentration of 180 ug/ml. Dilute to a working concentration of 1.0 ug/ml in PBS, without carrier protein.

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Lyophilized

Storage Conditions/Comments: Lyophilized products are stable for one year at 4°C until the expiration date listed above. Upon reconstitution, the product can be stored at 4°C for short term only (up to 60 days), or at -20°C to -80°C in aliquots for up to 6 months. Avoid repeated freeze-thaw cycles. Refreezing should be performed on dry ice or in a dry ice/alcohol bath. For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: December 3, 2014





pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: pbl assaysci.com

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.

Rev. 00

