

TECHNICAL DATA SHEET

WR304 MOUSE MONOCLONAL ANTIBODY (IGM), BIOTINYLATED ANTI-(PIP/PIP₂)

Background:

The WR304 antibody was produced by the Walter Reed Army Institute of Research (WRAIR) using an advanced, targeted Lipid-A adjuvant protocolⁱ. The antigen was a lipid, porcine brain PI(4)Pⁱ. WR304 is a monoclonal antibody that binds specifically to PIP, and PIP₂ⁱⁱ. It neutralizes infectious HIV-1 virusⁱⁱⁱ.

Antibody Information:

Antigens: Brain PI(4)P.

Ig Class: Mouse IgM (kappa).

Specificity: WR304 recognizes PIP and PIP₂.

Antibody Source: Monoclonal antibody from BALB/c-derived hybridoma WR304.

Production: In vitro cell culture.

Purification: Ultra filtration through 100 KDa cut-off filters.

Purity: ≥ 95%.

Formulation: WR304 Biotin is provided as a sterile-filtered solution in Tris buffered saline (TBS). WR304 binding to porcine brain PI(4)P is inhibited by high concentration of Ca²⁺ (10 mM)^{iv} and addition of 1 mM EDTA stimulates this binding activity. PBS and other phosphate-containing buffers should not be used as they inhibit binding to the phospholipid head groupⁱⁱⁱ.

Mass and Concentration: Refer to Product Label.

Recommended Applications: ELISA, IHC, Dot Blot, Flow cytometry and HIV-1 infectivity.

Storage conditions: Store undiluted at either -20°C or -80°C.

Hazardous/Non-hazardous Components: This product contains no substances that, at their given concentration, are known to be hazardous to health. Therefore, there is no MSDS for this product.

Product use:

The WR304 antibody has been used for the quantitation of PIP in direct ELISA and Dot Blot analysis. The WR304 antibody specifically binds to PIPⁱ and inhibits the infectivity of HIV-1 to peripheral blood

Avanti No.	Description	No. of Assays
330021S	WR304 monoclonal	100
330022S	WR304 monoclonal biotinylated	100

mononuclear cells (PBMC)ⁱⁱⁱ. A biotinylated form of WR304 is recommended for use in enzyme-linked immunosorbent assay (ELISA) determination of PIP in serum or plasma samples.

Note:

After thawing, centrifuge this product at > 1,000 g for 5 minutes to collect any antibody solution that may be retained in the cap.

The recommended long-term storage for WR304 is at -80°C. After initial thawing remaining product should be refrozen at -80°C.

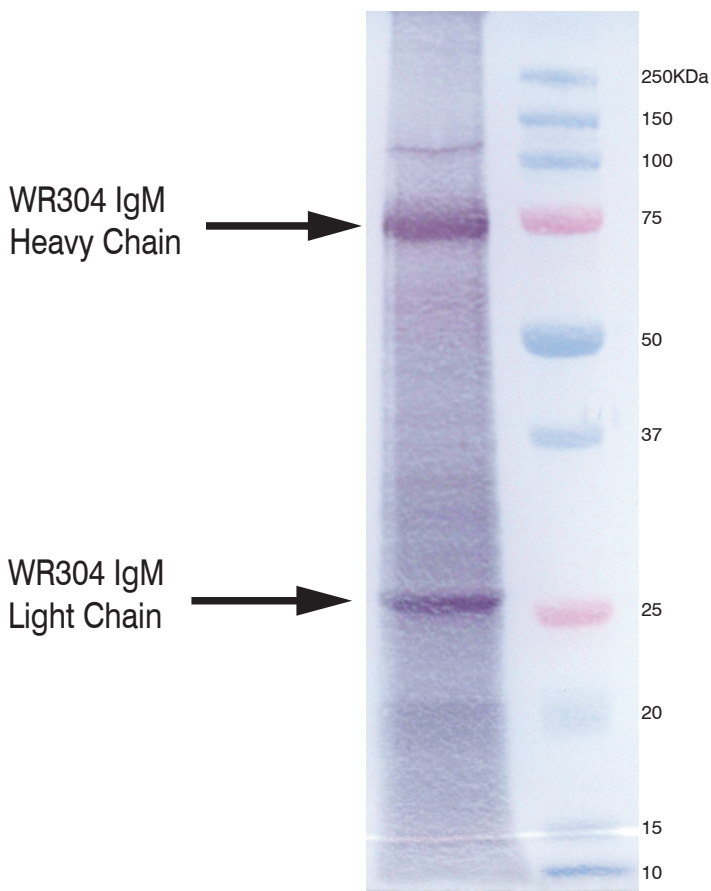


Fig. 1: Western blot of a SDS-PAGE gel of purified WR304-Biotin antibody. To detect biotin-labeled IgM protein subunits, the blot was probed with streptavidin-HRP using TMB as the substrate.

APPLICATIONS:

1. Direct ELISA Protocol

96 well “U” bottom Immulon 2HB plates (Thermo 3655) were used for lipid ELISA. Stock solutions of brain PI(4)P in CHCl_3 :MeOH:H₂O (20:9:1) were diluted in methanol to 10 μM . 1 nmol in 100 μL was added to each well and the solvent was allowed to evaporate O/Nⁱⁱ. The plates were blocked with 0.3% Gelatin in TBS (0.8% NaCl, 20 mM Tris-HCl pH 7.4, Mediatech) containing 1 mM EDTA for 1h at room temperature^{iv}. The wells were washed four times with TBS containing 1 mM EDTA. 10 μL aliquots of a serial dilution the WR304 antibody (400-1.6 ng/100 μL) in TBS containing 0.3% Gelatin, 1 mM EDTA were added to each well and incubated for 1h at RT. The amount of IgM bound to each well was quantified with streptavidin-HRP. After 1h incubation at room temperature, unbound secondary was washed

off with TBS. ABTS (KPL, Inc.) was added as substrate and absorbance was read at 405 nm with a correction for light scattering at 650 nm. Alternatively, using 96 well “U” bottom Microfluor 1 white plates (Thermo 6905) with the above protocol, the amount of biotinylated IgM bound to each well was quantified with streptavidin-AP and a LumiPhos 530 substrate after 1hr incubation at room temperature (Fig 2).

FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.

References:

ⁱWassef NM, Roerdink F, Swartz GM Jr, Lyon JA, Berson BJ, Alving CR. Phosphate-binding specificities of monoclonal antibodies against phosphoinositides in liposomes. *Mol. Immunol.*(1984) 21: 863-868.

ⁱⁱMatyas GR, Beck Z, Karasavvas N, Alving CR. Lipid binding properties of 4E10, 2F5, and WR304 monoclonal antibodies that neutralize HIV-1. *Biochim. Biophys. Acta* (2009) 1788: 660-665.

ⁱⁱⁱBrown BK, Karasavvas N, Beck Z, Matyas GR, Birx DL, Polonis VR, Alving CR. Monoclonal antibodies to phosphatidylinositol phosphate neutralize human immunodeficiency virus type 1: role of phosphate-binding subsites. *J Virol.* (2007) 81:2087-2091

^{iv}Beck Z, Karasavvas N, Tong J, Matyas GR, Rao M, Alving CR. Calcium modulation of monoclonal antibody binding to phosphatidylinositol phosphate. *Biochem. Biophys. Res. Commun.* (2007) 354: 747-751.

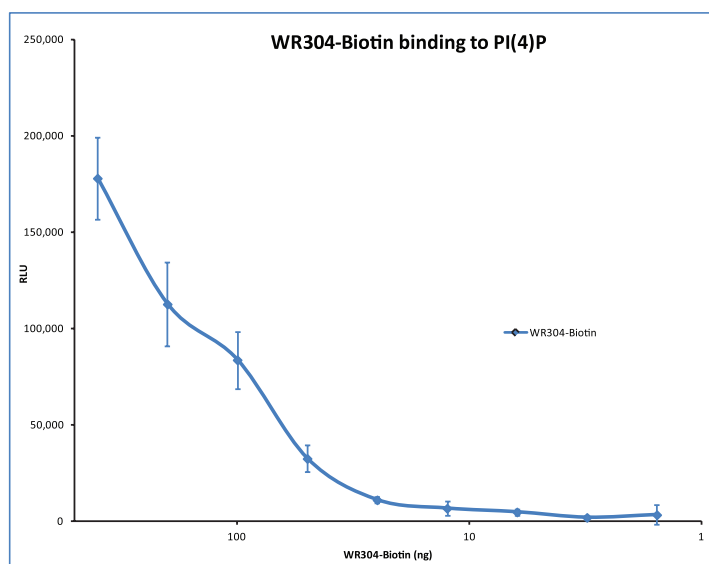
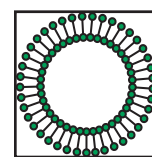


Fig. 2. Sample Direct ELISA

Immobilized antigen, porcine brain PI(4)P probed with WR304 biotin antibody followed by a streptavidin-AP. Bound WR304 biotin was measured as AP-dependent luminescence using Lumi-Phos 530 as substrate.



For more details: www.avantilipids.com
Phone: 800-227-0651
Email: lipidomics@avantilipids.com



Avanti[®]
POLAR LIPIDS, INC.

Revised 061714