



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Certificate of Analysis and Data Sheet

Goat Anti Axl Receptor Tyrosine Kinase

With HRP-conjugated Secondary Antibody

Catalog No.
DS-PB-00216

Target Species
Mouse

Isotype
IgG1

Preparation

Immunogen: Peptide sequence RGSPAAPGQEDGA corresponding to the C-terminus of AXL (NP_068713).

Preparation: Antisera to AXL receptor tyrosine kinase were raised by repeated immunizations of goats with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Synonyms: ARK, UFO

Specificity

DS-PB-00216 recognizes AXL receptor tyrosine kinase, a type 1 transmembrane protein. AXL is a member of the receptor tyrosine kinase subfamily that binds growth factors like vitamin K-dependent protein growth-arrest-specific gene 6.

AXL has been reported to be involved in the stimulation of cell proliferation and has transforming potential in cases of chronic myelogenous leukaemia.

Species Crossreactivity: Based on sequence similarity, DS-PB-00216 is expected to react with human.

Formulation

Product Type: Polyclonal Antibody

Product Form: Purified IgG – liquid

Preservative Stabilizers: 0.02% Sodium Azide, 0.5% Bovine Serum Albumin

Buffer Solution: TRIS buffered saline pH7.3

Approx. Protein Concentrations: IgG concentration 0.5 mg/ml

Fusion Partners: Spleen cells from immunized BALB/c mice were fused with cells of the mouse X63-Ag.8.653 myeloma cell line.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Storage

Store at +4°C or at -20°C if preferred. This product should be stored undiluted.

Storage in frost-free freezers is not recommended. Avoid repeated freezing and thawing as this may denature the antibody. Should this product contain a precipitate we recommend microcentrifugation before use.

Shelf Life: 18 months from date of dispatch

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA			•	
Western Blotting (1)	•			0.5ug/ml - 1ug/ml
Immunohistology – frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunofluorescence staining			•	
Functional Assays			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

(1) Detects a band of approximately 90kDa in NIH3T3 cell lysates. An overnight incubation time of 12 hours is recommended for this antibody.

Secondary Antibody Applications

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Immunoassay (ELISA, Western blot)	•			1:5000-1:10000

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

References

1. Varnum, B.C. *et al.* (1995) Axl receptor tyrosine kinase stimulated by the vitamin K-dependent protein encoded by growth-arrest-specific gene 6 *Nature*. 373: 623-626.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**