

# Amersham ECL Anti-mouse IgG, peroxidase-linked whole antibody (from sheep) affinity purified general purpose reagent

## Product Specification Sheet

### Introduction

#### Important

Read these instructions carefully before using the products.

#### Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

#### Safety

All chemicals should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.

#### Storage

Store at 2–8°C. Do not freeze. Avoid repeated freeze-thaw cycles. Under these conditions the product is stable for six months from date of despatch.

#### Expiry

See outer packaging.

### Component

Horseradish Peroxidase conjugated antibody is supplied in Phosphate Buffered Saline (Sodium Phosphate 0.1 M, NaCl 0.1 M) pH7.5, containing 1%(w/v) Bovine Serum Albumin and an anti-microbial agent.

### Description

#### Purification to ensure species-specificity

The antibody is prepared by hyper-immunizing sheep with purified immunoglobulin fractions from normal mouse serum to produce high affinity antibodies.

The pooled antiserum is used to produce an immunoglobulin preparation which is then purified on affinity columns of mouse IgG. After washing to remove any remaining non-specific serum components and low affinity antibodies, the specific antibodies are eluted using carefully selected, mild conditions which minimize aggregation and preserve immunological activity, yet which elute high affinity antibodies.

#### Preparation of labelled antibody

The enzyme Horseradish Peroxidase is attached to the immunoglobulin molecules using an adaptation of the periodate oxidation technique (1). This method has been found not to affect the effective binding of the antibody to the antigen or the activity of the enzyme.

#### Quality control

For every batch of enzyme-linked antibody that is produced the antibody titre is determined in an ELISA. The substrate used for the peroxidase is 2,2'-Azinobis[3-Ethylbenzothiazoline Sulphonate, diammonium salt], ABTS™.

Every batch is also QC tested in a Western blotting system. This is performed using Hybond™ ECL™ membrane containing tubulin protein and immunodetected with: primary antibody, Monoclonal anti-tubulin; and secondary antibody NXA931, anti-mouse IgG, HRP GPR. Blots are detected using ECL and ECL Plus detection systems.

### Applications

#### Protein blotting

##### 1. Detection with ECL(2) Western blotting reagents

This reagent has been shown to be suitable for use in ECL Western blotting applications. The control system used was the detection of monoclonal anti-tubulin.

We have found in our laboratories that dilutions of: 1:2000 for monoclonal anti-tubulin; and 1:5000 for anti-mouse IgG, HRP GPR are suitable for the detection of 3 ng of tubulin on Hybond ECL membrane, exposed to Hyperfilm™ ECL for 5 minutes.

To achieve the same sensitivity level on Hybond-P PVDF, concentrations would typically be: anti-tubulin - 1:2000; and NXA931 - 1:10 000.

##### 2. Detection with ECL Plus(3,4) Western blotting reagents

ECL Plus Western blotting reagent is highly sensitive, giving an increase, for this antibody, of 4-20 fold over ECL detection.

This property can be utilized in 2 ways:

- Preservation of antibodies that are rare or costly
- Increase in detectable sensitivity levels

The control system used was the same as for ECL.

The suitable antibody dilutions, to detect 3 ng of tubulin on Hybond ECL membrane are: anti-tubulin - 1:5000; and NXA931 - 1:25 000.

For Hybond-P PVDF antibody dilutions are typically: anti-tubulin - 1:10 000; and NXA931 - 1:50 000.

### 3. Colorimetric detection

A dilution of 1:300 is recommended.

## ELISA

If this reagent is to be used to detect mouse immunoglobulins, we have found in our laboratories that a dilution of 1:6000 is suitable for the detection of 1 µg of IgG. For greater sensitivity (for example down to 300 pg) the reagent should be diluted rather less (for example 1:500). Thus 1.0 ml of stock reagent will be sufficient for up to 60000 wells at the higher dilution if used at 0.1 ml per well in standard microtitre plates. A suitable diluent is Phosphate-Buffered Saline containing 0.05%(v/v) Tween™ 20.

## Immunocytochemistry

Whilst potentially applicable to immunocytochemistry this product is not primarily intended for this use. It has not been adsorbed against immunoglobulins of other species to remove cross-reactivities. Therefore, it is likely that if the reagent is used for immunocytochemistry there may be some 'background' staining. More suitable reagents for immunocytochemistry are peroxidase labelled anti-mouse IgG, whole antibody (NA931) or peroxidase labelled anti-mouse IgG, F(ab')<sub>2</sub> fragments (NA9310), both of which are species-specific.

## Protocol recommendations

### Membranes

Nitrocellulose and PVDF membranes are suitable for use with both detection systems. PVDF membrane is highly recommended for use with ECL Plus detection reagents.

For high quality results the following guidelines should be followed:

**Blocking:** Use enough blocking agent to block all non-specific sites. A typical block is 5% non-fat dried milk in PBS Tween or TBS Tween. See 'Tech-Tips' No. 136 available from Cytiva, for further details.

**Washing:** The volume of wash buffer (eg PBS-T or TBS-T) must be sufficient to cover the membrane completely.

## Optimization of primary and secondary antibodies

### ECL detection

ECL Western blotting is a very sensitive technique. As such it is essential to optimize the system under study for high signal and low background for both primary and secondary antibodies.

Dot blots are a quick and effective method of determining the optimum dilutions required for primary and secondary antibodies. Optimization details are set out in the RPN2106/2108/2109/2209/2134 booklets and 'Tech-Tips' No. 129 available from Cytiva.

## ECL Plus detection

Due to the improved sensitivity of ECL Plus compared to ECL, optimization details as set out in the RPN2132/2133 booklets and 'Tech-Tips' No. 169 available from Cytiva are recommended.

## Typical anti-mouse secondary antibody dilution ranges:

ECL for nitrocellulose membrane	1:1000 to 1:5000
ECL Plus for nitrocellulose membrane	1:2000 to 1:10000

For PVDF membrane the use of higher dilutions may be necessary.

The exact concentration of the secondary antibody will always be dependent upon the primary antibody used and the sensitivity and exposure times required.

Detection: Ensure any excess ECL or ECL Plus detection reagents are sufficiently drained prior to exposure.

### Exposure times:

ECL - exposure times of 1 to 15 minutes are suggested.

ECL Plus - initial exposure times of 1 to 5 minutes are suggested.

Signal can still be obtained up to 24 hours after the application of ECL Plus reagents, and for this exposure times of 1 to 2 hours may be required.

## Related Products

ECL Western blotting detection reagents	RPN2106/2108/2109/2209/2134
ECL Plus Western blotting detection system	RPN2132/2133
Hybond ECL membrane	RPN2020D
Hybond-P PVDF membrane	PN2020F
Hyperfilm ECL	RPN2103/2104/1681/1674
ECL protein molecular weight markers	RPN2107NA9320

## References

1. NAKANE, P.K. and KAWAOI, A., *Journal of Histochemistry and Cytochemistry*, **22**, pp.1084-1091, 1974.
2. WHITEHEAD, T.P. *et al.*, *Clin. Chem.*, **25**, pp.1531-1546, 1979.
3. AKHAVEN-TAFTI, H. *et al.*, *Clin. Chem.*, **41**, pp.1368-1369, 1995.
4. AKHAVEN-TAFTI, H. *et al.*, *Biolum. And Chemilum. Fundamentals and Applied Aspects*, pp.199-202, Chichester, 1994.

## cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Amersham™, ECL, Hybond, and Hyperfilm are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

ABTS is a trademark of Boehringer Mannheim Corp.

Tween is a trademark of Croda Group of Companies.

All other third-party trademarks are the property of their respective owners.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

NXA-931PS AE V:4 03/2021

