

Anti-HA-Biotin, High Affinity (3F10)

Monoclonal antibody for the highly sensitive detection of HA-tagged recombinant proteins, F_{ab} fragments conjugated with biotin

Cat. No. 12 158 167 001 50 µg

Version 06
Content version: July 2019

Store at +2 to +8°C

1. Product characteristics

| | |
|-----------------------------------|--|
| Antibody type | Clone BMG-3F10, rat IgG ₁ , F _{ab} fragments |
| Specificity | Anti-HA-Biotin, High Affinity (3F10) recognizes the HA peptide sequence [YPYDVPDYA] derived from the human influenza hemagglutinin protein (1). The antibody recognizes its antigenic determinant even when the HA peptide epitope is introduced into unrelated recombinant proteins by a technique known as "epitope tagging". |
| Formulation | White lyophilizate, lyophilized in the presence of proteinous stabilizers. |
| Storage and stability | The lyophilized Anti-HA-Biotin, High Affinity (3F10) is stable for 24 months or through the expiration date printed on the label when stored at 2–8°C. After reconstitution, the conjugate is stable for 2 month at +2 to +8°C. Alternatively, it can be stored in aliquots at –15 to –25°C. The reconstituted conjugate is stable for 6 month at –15 to –25°C. Repeated freezing and thawing must be avoided. The Anti-HA-Biotin, High Affinity (3F10) is shipped at RT. |
| Reconstitution and storage | <ul style="list-style-type: none">• Reconstitute the lyophilizate in 1.0 ml double distilled water for 10 minutes at 15 to 25°C (RT), and mix thoroughly but do not vortex.• This results in a final concentration of 50 µg/ml.• The reconstituted antibody is stable for –2 months when stored at 2–8°C, or for–6 months when stored in aliquots at –15 to –25°C. Repeated freeze/thaw cycles must be avoided ! |

Application The following table lists the possible applications and recommended working concentrations:

| Application | Working concentration |
|------------------------|-----------------------|
| Western blot, Dot blot | 100 ng/ml |
| ELISA | 100 ng/ml |

Quality control The Anti-HA-Biotin, High Affinity (3F10) antibody is function tested by Western blot analysis of an HA-tagged fusion protein.

Advantages

| Benefits | Features |
|-------------|---|
| Sensitivity | Anti-HA-Biotin, High Affinity (3F10) provides superior detection of HA-tagged proteins in the picogram range. |
| Specificity | No cross reactivity compared to other Anti-HA antibodies. |
| Flexibility | The biotin conjugated antibody allows assay establishing using the universal biotin-streptavidin platform. |

2. Background information

Epitope tagging Epitope tagging, the fusion of a short stretch of amino acids to a protein of interest by recombinant techniques, is a widely used method that allows the surveillance of the fusion protein with tag-specific monoclonal antibodies. The epitope tagging approach offers the ability to determine

- the size, cellular localization, and abundance of proteins produced by newly discovered genes
- post-translational modifications of proteins
- the movement of proteins within cells
- the identity of proteins within functional protein complexes
- the function of proteins that are unstable, difficult to purify, or share epitopes with a number of other proteins and
- eliminates need to generate specific antibodies recognizing the protein of interest (1-6).

Anti-HA antibodies in epitope tagging Among the different epitope-tags described in the literature the most commonly used tag is derived from the hemagglutinin of the influenza virus (HA1; 2). Several antibodies have been described that react with this epitope tag, the most prominent one is Anti-HA (2; 12CA5). However these antibodies are restricted by requiring additional amino acids adjacent to the HA tag or by recognizing HA-tagged proteins with only moderate affinity, as demonstrated by cross-reacting bands that have been reported in certain Western blot experiments (7).

Anti-HA, High Affinity The Anti-HA High Affinity antibody (clone 3F10) recognizes the same epitope as clone 12CA5. It is a monoclonal antibody whose high affinity and low working concentrations result in less cross reactivity than with other antibodies to the HA-epitope. Anti-HA-Biotin, High Affinity (3F10) is a biotin conjugate of this clone which is especially useful in Western blotting, ELISA applications and assays using the universal biotin-streptavidin platform by allowing specific and highly sensitive detection of HA-tagged proteins.

3. Applications

3.1 Procedure for Western blotting

Introduction The following procedure describes the detection of a HA-tagged protein by enzyme-mediated chemiluminescence.
If using other detection systems *e.g.*, colorimetric detection, the conditions may vary and should be adapted.

Before you begin For separation by gel electrophoresis and blotting, please refer to reference 8.

Additional material required

The following table lists additional products from Roche Diagnostics necessary to perform the Western blotting procedure.

| Product | Cat. No. |
|---|--|
| PVDF Western Blotting Membranes | 11 722 034 001 11 722 026 001 |
| Tween 20 | 11 332 465 001 |
| BM Chemiluminescence Blotting Substrate (POD) | 11 500 708 001 11 500 694 001 |
| Blocking Reagent | 11 096 176 001 |
| Lumi-Film Chemiluminescent Detection Film | 11 666 916 001 11 666 657 001 11 666 711 001 |
| Na ₂ HPO ₄ , analysis grade | |
| NaH ₂ PO ₄ , analysis grade | |

Preparation of working solutions

| Working solution | Composition or preparation | Storage and stability | Use |
|---------------------------------------|--|--|---|
| Phosphate buffered saline (PBS), 10 × | 100 mM phosphate, 1.5 M NaCl, pH 7.2 | stable for • 1 week at +2 to +8°C, or • at least 2 years at -15 to -25°C | Preparation of 1 × PBS |
| PBS, 1 × | Dilute 10 ml 10× PBS with double dist. water to make 100 ml | stable for • 1 week at +2 to +8°C, or • at least 2 years at -15 to -25°C | • Preparation of blocking solution • Preparation of Washing solution |
| Washing solution | PBS, 1 ×, containing 0.1% Tween 20 (v/v) | stable for 1 week at +2 to +8°C | Washing |
| Blocking solution | PBS, 1 ×, containing 1% Blocking Reagent (w/v) | stable for • 1 week at +2 to +8°C, or • at least 2 years at -15 to -25°C | • Blocking • Preparation of Anti-HA-Biotin Working solution |
| Anti-HA-Biotin solution | Dilute the reconstituted antibody to 100 ng/ml using the Blocking solution | unstable, prepare shortly before use | Detection |

Important notes regarding sample preparation

Prepare protein extracts containing the HA tagged protein of interest using a variety of standard methods (8). The following lysis buffers have performed well:

- Bacterial extracts: 20 mM Tris, pH 8.0; 100 mM NaCl, Complete¹ Protease Inhibitor. (followed by sonication/freeze-thaw).
- Mammalian extracts: 50 mM Tris, pH 7.5; 150 mM NaCl, 1% Nonidet 40, 0.05% Deoxycholate, Complete Protease Inhibitor.

Other cell lysis buffers may be more appropriate for individual applications.

- Include protease inhibitors to reduce proteolytic activity. Use Complete tablets* for most applications.
- Limit detergents to the lowest concentration levels necessary to obtain adequate cell lysis.

Procedure for immunodetection

| Step | Action |
|------|--|
| 1 | After electrophoresis and transfer of the proteins to a PVDF membrane, block the membrane with blocking solution for 1 h at +37 °C or for 3 h at +15 to +25°C. |
| 2 | Incubate the blot with 100 ng/ml Anti-HA-Biotin working solution for 1 h at +15 to +25°C. |
| 3 | Wash 3×, 5 min each, with washing solution. |
| 4 | Incubate the blot with 15 mU/ml Anti-Biotin-Peroxidase* or 15 mU/ml Streptavidin-Peroxidase* diluted in Blocking solution for 1 h at +15 to +25°C. |
| 5 | Wash 3×, 5 min each, with washing solution. |
| 6 | Detect bound immunocomplexes with a chemiluminescence substrate as described in the package insert of the BM Chemiluminescence Blotting Substrate (POD). |

Typical result

The following picture shows a typical result regarding the specificity and sensitivity of the detection of HA-tagged proteins by Western blotting.

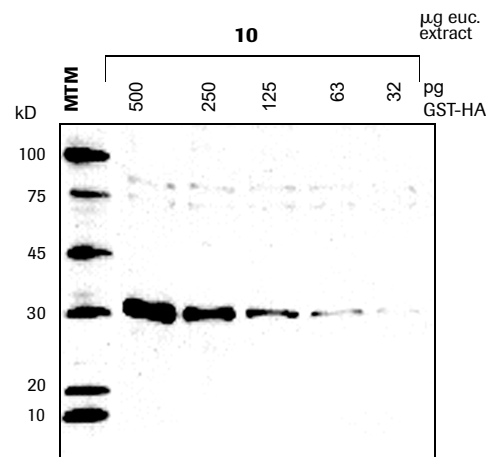


Fig. 1: Western blot analysis of HA-tagged Glutathion-S-transferase (GST-HA) detected with Anti-HA-Biotin, High Affinity (3F10).

Purified GST-HA was serially diluted to the indicated amounts in 10 µg of protein from eucaryotic cell extract. HA-tagged proteins were detected with 100 ng/ml Anti-HA-Biotin; High affinity (3F10), 20 mU/ml Anti-Biotin-Peroxidase, and BM Chemiluminescence Blotting substrate (POD), used according to the substrate's package insert (3 min exposure). The observed background activity is derived from non-specific binding of the secondary detection antibody (data not shown).

MTM: Multi-Tag-Marker

3.2 Procedure for ELISA

Before you begin

For detailed information, please refer to reference No. 8.

Additional equipment required

- Microtiter plates (e.g., Nunc Maxisorp)
- Microtiter plate washer (optional)
- Microtiter plate reader

Additional reagents required

The following table lists additional products required for the ELISA procedure

| Product | Cat. No. |
|---------------------------------------|----------------|
| Tween 20 | 11 332 465 001 |
| Blocking Reagent | 11 096 176 001 |
| BM Blue POD Substrate, soluble | 11 484 281 001 |
| Sodium carbonate, analysis grade | |
| Sulfuric acid, 95–97%, analysis grade | |

Preparation of working solutions

| Working solution | Composition or preparation | Storage and stability | Use |
|---------------------------------------|---|--|---|
| Sodium carbonate solution | 50 mM, pH 9.6 | prepare shortly before use | Coating |
| Phosphate buffered saline (PBS), 10 × | 100 mM phosphate, 1.5 M NaCl, pH 7.2 | stable for <ul style="list-style-type: none"> • 1 week at 2–8°C, or • at least 2 years at –15 to –25°C | Preparation of 1× PBS |
| PBS, 1× | Dilute 10 ml 10 × PBS with double dist. water to make 100 ml | stable for <ul style="list-style-type: none"> • 1 week at 2–8°C, or • at least 2 years at –15 to –25°C | <ul style="list-style-type: none"> • Preparation of blocking solution • Preparation of Washing solution |
| Washing solution | 1× PBS containing 0.1 % Tween 20 (v/v) | stable for 1 week at 2–8°C | Washing |
| Blocking solution | 1× PBS containing 1% Blocking Reagent (w/v) | stable for <ul style="list-style-type: none"> • 1 week at 2–8°C, or • at least 2 years at –15 to –25°C | <ul style="list-style-type: none"> • Blocking • Preparation of Anti-HA-Biotin Working solution |
| Coating solution | Dilute 1–10 µg of the appropriate protein in 1 ml sodium carbonate solution | prepare shortly before use | Coating |
| Anti-HA-Biotin working solution | Dilute the reconstituted antibody to 100 ng/ml using the Blocking solution | unstable; prepare shortly before use | Detection |

Procedure for ELISA

Cover the plate either with microtiter plate covers or adhesive sealing film during all incubation steps in order to avoid evaporation of the solutions.

| Step | Action |
|------|--|
| 1 | Coat the wells with 100 µl/well coating solution for 1–2 h at 37°C or over night at +2 to +8°C. |
| 2 | Wash 5 × with washing solution and remove residual washing solution. |
| 3 | Add 300 µl blocking solution per well and incubate for 1–2 h at 37 °C or over night at +2 to +8°C. |
| 4 | Wash 5 × with washing solution and remove residual washing solution. |
| 5 | Add 100 µl Anti-HA-Biotin working solution per well, and incubate for 1 h at +15 to +25°C. |
| 6 | Wash 5 × with washing solution and remove residual washing solution. |
| 7 | Add 100 µl Anti-Biotin-Peroxidase (15 mU/ml in blocking solution) per well, and incubate for 10 min at +15 to +25°C. |
| 8 | Wash 5× with washing solution and remove residual washing solution. |
| 9 | Add 100 µl/well RMB Blue POD Substrate, soluble, prewarmed to +15 to +25°C, and incubate at +15 to +25°C and under constant shaking until the color development is sufficient. |
| 10 | To stop the color development, add 100 µl/well 2 N sulphuric acid. |
| 11 | Read the absorbance at 450 nm (reference wavelength: 690 nm) within 30 min after stopping the reaction. |

4. Appendix

4.1 Trouble shooting

| Problem | Possible Cause | Recommendation |
|--|---|---|
| Nonspecific reactivity especially with high total protein loading. | <ul style="list-style-type: none"> • Nonspecific binding of secondary antibody. • Inadequate buffer conditions. • High Anti-HA-Biotin antibody concentration | <ul style="list-style-type: none"> • Optimise assay conditions by reducing the concentration of the secondary antibody. • Prolong time for blocking the membrane. • Reduce amount of total protein loaded • Use PBS containing Blocking reagent for membrane blocking, dilution of the Anti-HA-Biotin and dilution of the secondary detection antibody. • Reduce Anti-HA-Biotin antibody concentration |
| Staining of the protein of interest is too weak. | <ul style="list-style-type: none"> • Inadequate amounts of protein loaded onto the gel. • Inadequate conditions used for detection. | <ul style="list-style-type: none"> • Increase the amount total protein loading. • Increase the concentration of Anti-HA-Biotin. • Prolong exposure time used during detection. |
| Staining of the protein of interest is too strong. | <ul style="list-style-type: none"> • Inadequate amounts of protein loaded onto the gel. • Inadequate conditions used for detection | <ul style="list-style-type: none"> • Decrease the amount total protein loading. • Decrease the concentration of Anti-HA-Biotin. • Decrease the concentration of the secondary detection reagent. • Shorten exposure time used during detection. |

4.2 References

- Kolodziej, P.A. & Young, R.A. Epitope Tagging and Protein Surveillance. *Meth. Enzymol.* **194**, 508–511.
- Wilson, I.A., Niman, H.L., Houghten, R.A., Chersonson, A.R., Connolly, M.L. & Lerner, R.A. (1984): The structure of an antigenic determinant in a protein. *Cell* **37**, 767–778.
- Field, J., Nikawa, J.-I., Broek, D., MacDonald, B., Rodgers, L., Wilson, I.A., Lerner, R.A. & Wigler, M. (1988): Purification of a RAS-Responsive Adenylyl Cyclase Complex from *Saccharomyces cerevisiae* by Use of an Epitope Addition Method. *Mol. Cell. Biol.* **8**, 2159–2165.
- von Zastrow, M. & Kobilka, B.K. (1992): Ligand-regulated Internalization and Recycling of Human β2-Adrenergic Receptors between the Plasma Membrane and Endosomes Containing Transferrin Receptors. *J. Biol. Chem.* **267**, 3530–3538.
- Emrich, T., Förster, R. & Lipp, M. (1994): Topological Characterization of the Lymphoid-Specific Seven Transmembrane Receptor BLR1 by Epitope-Tagging and High Level Expression. *Biochem. Biophys. Res. Com.* **197**, 214–220.
- Qian, N.-X., Winitz, S. & Johnson, G.L. (1993): Epitope-tagged Gq α subunits: Expression of GTPase-deficient α subunits persistently stimulates phosphatidylinositol-specific phospholipase C but not mitogen-activated protein kinase activity regulated by the M1 muscarinic acetylcholin receptor. *Proc. Natl. Acad. Sci.* **90**, 4077–4081.
- Chen, Y.-T., Holcomb, C. & Moore, H.-P.H. (1993): Expression and localization of two low molecular weight GTP-binding proteins, Rab8 and Rab19, by epitope tag. *Proc. Natl. Acad. Sci.* **90**, 6508–6512.
- Harlow, E. & Lane, D. (1988): *Antibodies: A Laboratory Manual*, Cold Spring Harbour Laboratory Press, Cold Spring Harbour, N.Y.

* available from Roche Diagnostics

COMPLETE is a trademark of Roche.

Nonidet is a trademark of Shell International Petroleum Company Limited, U.K.

Tween is a trademark of ICI Americas, Inc., USA.

Changes to previous version

Editorial changes.

Trademarks

All third party product names and trademarks are the property of their respective owners.

Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

4.3 Related products

| Product | Pack size | Cat. No. |
|---|--|----------------|
| Anti-HA, High Affinity (3F10) lyophilized, unconjugated | 50 µg | 11 867 423 001 |
| | 500 µg | 11 867 431 001 |
| peroxidase | 25 U (25 µg) | 12 013 819 001 |
| HA peptide | 5 mg | 11 666 975 001 |
| Anti-HA (12CA5) lyophilized, unconjugated | 200 µg | 11 583 816 001 |
| in solution, unconjugated | 5 mg | 11 666 606 001 |
| biotin | 100 µg (500 µl) | 11 666 851 001 |
| fluorescein | 100 µg (500 µl) | 11 666 878 001 |
| rhodamine | 100 µg (500 µl) | 11 666 959 001 |
| peroxidase | 50 µg (500 µl) | 11 667 475 001 |
| Anti-Biotin-Peroxidase | 150 units | 11 426 303 001 |
| Streptavidin-Peroxidase | 500 units | 11 089 153 001 |
| Anti-c-myc lyophilized, unconjugated | 200 µg | 11 667 149 001 |
| in solution, unconjugated | 5 mg | 11 667 203 001 |
| peroxidase | 500 µg (500 µl) | 11 814 150 001 |
| Anti-VSV-G lyophilized, unconjugated | 200 µg | 11 667 351 001 |
| Anti-His₆ | 100 µg | 11 922 416 001 |
| Anti-His₆-Peroxidase | 50 units | 11 965 085 001 |
| Anti-GFP | 200 µg | 11 814 460 001 |
| rGFP | 50 µg | 11 814 524 001 |
| c-myc peptide | 5 mg | 11 667 246 001 |
| Multi-Tag Marker | 250 µl | 11 828 649 001 |
| X-tremeGENE siRNA Transfection Reagent | 1 ml, 400 Transfections | 04 476 093 001 |
| | 5 × 1 ml, 2000 transfections | 04 476 115 001 |
| FuGENE 6 Transfection Reagent | 1 ml | 11 814 443 001 |
| DOSPER Liposomal Transfection Reagent | 2 ml (5 × 0.4 ml) | 11 781 995 001 |
| | 0.4 ml | 11 811 169 001 |
| DOTAP Liposomal Transfection Reagent | 2 ml (5 × 0.4 ml) | 11 202 375 001 |
| | 0.4 ml | 11 811 177 001 |
| Lumi Light^{PLUS} POD Western Blotting Kit | 1000 cm ² | 12 015 218 001 |
| Lumi Light^{PLUS} POD Western Blotting Substrate | 100 ml | 12 015 196 001 |
| Lumi Light POD Western Blotting Substrate | 400 ml | 12 015 200 001 |
| BM Blue POD Substrate, precipitating | 100 ml | 11 442 066 001 |
| Lumi-Film Chemiluminescent Detection Film | 100 films (8 × 10 inches 20.3 × 25.4 cm) | 11 666 657 001 |
| Complete, EDTA-free Protease Inhibitor Cocktail Tablets | 20 tablets (each sufficient for 50 ml extract) | 11 873 580 001 |
| Complete, Mini, EDTA-free Protease Inhibitor Cocktail Tablets | 25 tablets (each sufficient for 10 ml extract) | 11 836 170 001 |

Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit sigma-aldrich.com, and select your home country. Country-specific contact information will be displayed.



Roche Diagnostics GmbH
Sandhofer Strasse 116
68305 Mannheim
Germany