## PLAP Monoclonal Mouse Antibody (ALP/870)



## **Product Description**

Reacts with a 70 kDa membrane-bound isozyme (Regan and Nagao type) of Placental Alkaline Phosphatase (PLAP) occurring in the placenta during the 3rd trimester of gestation. It is highly specific for PLAP and shows no cross-reaction with other isozymes of alkaline phosphatase. Anti-PLAP reacts with germ cell tumors and can discriminate between these and other neoplasms. Somatic neoplasms e.g. breast, gastrointestinal, prostatic, and urinary cancers may also immunoreact with antibodies to PLAP. Anti-PLAP positivity in conjunction with anti-keratin negativity favors seminoma over carcinoma. Germ cell tumors are usually anti-keratin positive, but they regularly fail to stain with anti-EMA, whereas most carcinomas stain with anti-EMA. Anti-PLAP has been useful in the diagnosis of gestational trophoblastic disease.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0870, Anti-PLAP (ALPP/870)

## **Product attributes**

| Antibody number  | #0870  |  |  |
|--|--|--|--|
| Antibody reactivity (target)   | PLAP   |  |  |
| Antibody type  | Primary  |  |  |
| Host species   | Mouse  |  |  |
| Clonality  | Monoclonal   |  |  |
| Clone  | ALP/870  |  |  |
| Isotype  | IgG2b, kappa   |  |  |
| Molecular weight   | 70 kDa   |  |  |
| Synonyms   | Alkaline phosphatase placental type, Alkaline phosphatase<br>Regan Isozyme, ALP, Alp1, ALPP, Germ-cell alkaline<br>phosphatase, nagao Isozyme, PALP, Placental alkaline<br>phosphatase 1, placental heat-stable alkaline phosphatase,<br>PLAP-1, PLAP1   |  |  |
| Human gene symbol  | ALPP   |  |  |
| Entrez gene ID   | 250  |  |  |
| SwissProt  | P05187   |  |  |
| Unigene  | 284255   |  |  |
| Immunogen  | Recombinant full-length human ALPP protein   |  |  |
| Antibody target cellular<br>localization   | Plasma membrane  |  |  |
| Verified antibody applications   | IHC (FFPE) (verified), WB (verified)   |  |  |
| Species reactivity   | Human  |  |  |
| Antibody application notes   | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with   |  |  |
|  | secondary antibody, immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   |  |  |
| Positive control   | secondary antibody, İmmunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be  |  |  |
| Positive control<br>Shipping condition   | secondary antibody, İmmunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   |  |  |
|  | secondary antibody, Îmmunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  |  |  |
| Shipping condition   | secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,  |  |  |
| Shipping condition Storage Conditions  | secondary antibody, İmmunofluorescence: 0.5-1 ug/mL, Immunohistology fornalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when  |  |  |
| Shipping condition Storage Conditions Shelf life   | secondary antibody, İmmunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  |  |  |
| Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate                                    | secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA/0.05% azide, Purfied: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purfied: 1 mg/mL in PBS/0.05% BSA/0.05% BS |  |  |
| Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation                        | secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified; DS without azide  |  |  |
| Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Cell/tissue expression | secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.25-0.5 ug/mL, No special pretreatment is required for the immunohistochemical staining of formalin-fixed tissues, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user  HepG2 cells. Placenta or seminoma  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  Guaranteed for at least 24 months from date of receipt when stored as recommended  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide  Placenta  |  |  |

Ovarian cancer, Testicular cancer

Tumor expression

Call us: 800-304-5357 Email: btinfo@biotium.com

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel           | Dye Features     |
|-------------------|-------------|------------|------------|-----------------------------|------------------|
| BNC04             | CF®405S     | 404/431    | 405        | DAPI (microscopy),<br>AF405 | CF®405S Features |
| BNC88             | CF®488A     | 490/515    | 488        | GFP, FITC                   | CF®488A Features |
| BNC68             | CF®568      | 562/583    | 532, 561   | RFP, TRITC                  | CF®568 Features  |
| BNC94             | CF®594      | 593/614    | 561        | Texas Red®                  | CF®594 Features  |
| BNC40             | CF®640R     | 642/662    | 633-640    | Cy®5                        | CF®640R Features |
| BNC47             | CF®647      | 650/665    | 633-640    | Cy®5                        | CF®647 Features  |
| BNC74             | CF®740      | 742/767    | 633-685    | 775/50                      | CF®740 Features  |
| BNCB              | Biotin      | N/A        | N/A        | N/A                         |                  |
| BNUB              | Purified    | N/A        | N/A        | N/A                         |                  |
| BNUM              | Purified,   | N/A        | N/A        | N/A                         |                  |

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 1, 2025 at 11:57:58 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-plap-placental-alkaline-phosphatase-alp870/">https://biotium.com/product/monoclonal-anti-plap-placental-alkaline-phosphatase-alp870/</a>