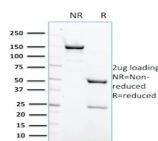


# Protein Tyrosine Phosphatase, non-receptor type 6 Monoclonal Mouse Antibody (CPTC-PTPN6-2)



## Product Description

PTPN6, which is also designated as SHP-1, is expressed primarily in hematopoietic cells and characterized by the presence of two SH2 domains N-terminal to the PTP domain. It maps to 12p13, a region commonly involved in leukemia-associated chromosomal abnormalities. SHP-1 plays a role in modulating intracellular signaling for various molecules. In T-cell lymphomas and other malignancies, SHP-1 phosphatase expression is lost, by a thus far undetermined mechanism. Preliminary studies suggest that it may be helpful in differentiating mantle/marginal zone lymphomas (SHP-1 ) from follicular lymphomas (SHP-1 negative).

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 [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 2259, Anti-Protein Tyrosine Phosphatase, non-receptor type 6 (CPTC-PTPN6-2)**

## Product attributes

|  |   |
|--|---|
| <b>Antibody number</b>                       | #2259   |
| <b>Antibody reactivity (target)</b>          | Protein Tyrosine Phosphatase, non-receptor type 6   |
| <b>Antibody type</b>                         | Primary   |
| <b>Host species</b>                          | Mouse   |
| <b>Clonality</b>                             | Monoclonal  |
| <b>Clone</b>                                 | CPTC-PTPN6-2  |
| <b>Isotype</b>                               | IgG2a, kappa  |
| <b>Molecular weight of antigen</b>           | 65 kDa  |
| <b>Synonyms</b>                              | PTPN6; SHP1; protein tyrosine phosphatase; non-receptor type 6; HCP; HCPH; HPTP1C; PTP-1C; SH-PTP1; SHP-1; SHP-1L; 70 kda SHP-1L protein; hematopoietic cell phosphatase; hematopoietic cell protein-tyrosine phosphatase; protein-tyrosine phosphatase 1C; PTPN6 / SHP   |
| <b>Human gene symbol</b>                     | PTPN6   |
| <b>Entrez gene ID</b>                        | 5777  |
| <b>SwissProt</b>                             | P29350  |
| <b>Unigene</b>                               | 63489   |
| <b>Immunogen</b>                             | Recombinant human full-length PTPN6 protein   |
| <b>Antibody target cellular localization</b> | Nucleus & cytoplasm   |
| <b>Species reactivity</b>                    | Human   |
| <b>Antibody application notes</b>            | For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry |
| <b>Positive control</b>                      | A431, Jurkat or K562 cell lysates; Lymph node or spleen.  |
| <b>Shipping condition</b>                    | Room temperature  |
| <b>Storage Conditions</b>                    | Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C   |
| <b>Regulatory status</b>                     | For research use only (RUO)   |
| <b>Antibody/conjugate formulation</b>        | 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   |
| <b>Shelf life</b>                            | Guaranteed for at least 24 months from date of receipt when stored as recommended   |
| <b>Product origin</b>                        | Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.   |

| Antibody # prefix | Conjugation        | Ex/Em (nm) | Laser line | Detection channel        | Dye Features                     |
|-------------------|--------------------|------------|------------|--------------------------|----------------------------------|
| BNC04             | CF®405S            | 404/431    | 405        | DAPI (microscopy), AF405 | <a href="#">CF®405S Features</a> |
| BNC88             | CF®488A            | 490/515    | 488        | GFP, FITC                | <a href="#">CF®488A Features</a> |
| BNC68             | CF®568             | 562/583    | 532, 561   | RFP, TRITC               | <a href="#">CF®568 Features</a>  |
| BNC94             | CF®594             | 593/614    | 561        | Texas Red®               | <a href="#">CF®594 Features</a>  |
| BNC40             | CF®640R            | 642/662    | 633-640    | Cy®5                     | <a href="#">CF®640R Features</a> |
| BNC47             | CF®647             | 650/665    | 633-640    | Cy®5                     | <a href="#">CF®647 Features</a>  |
| BNC74             | CF®740             | 742/767    | 633-685    | 775/50                   | <a href="#">CF®740 Features</a>  |
| BNCB              | Biotin             | N/A        | N/A        | N/A                      |                                  |
| BNUB              | Purified           | N/A        | N/A        | N/A                      |                                  |
| BNUM              | Purified, BSA-free | 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 June 3, 2026 at 05:18:14 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/protein-tyrosine-phosphatase-non-receptor-type-6-monoclonal-mouse-antibody-cptc-ptpn6-2/>