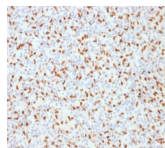


PU.1 Monoclonal Mouse Antibody (PU1/2146)



Product Description

PU.1 is a member of the ETS family of transcription factors and is important for normal B-cell development. It is expressed in the myeloid lineage and in immature as well as mature B-lymphocytes, with the exception of plasma cells. PU.1 is expressed in germinal center B-cells and mantle B-cells. Various lymphomas are also positive for this marker. It is essential during early B-cell differentiation. The absence of PU.1 results in total block of B-cell development at the pre-pro stage. PU.1 plays a key role in normal myeloid differentiation, and regulates the expression of immunoglobulin and other genes that are important for B-cell development. 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 to inquire about stock status and lead times before placing your order.

Product attributes

| | |
|---------------------------------------|--|
| Antibody number | #2146 |
| Antibody reactivity (target) | PU.1 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | PU1/2146 |
| Isotype | IgG2b, kappa |
| Molecular weight | 40 kDa |
| Synonyms | Transcription Factor spi1; 31kDa Transforming Protein; Hematopoietic Transcription Factor PU.1; SFP11; SPI1; SPIA; Spleen focus forming virus (SFFV) proviral integration oncogene spi1 |
| Human gene symbol | SPI1 |
| Entrez gene ID | 6688 |
| SwissProt | P17947 |
| Unigene | 502511 |
| Immunogen | Recombinant fragment (around aa 16-170) of human PU.1 protein (Exact sequence is proprietary) |
| Verified antibody applications | Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified) |
| Antibody target cellular localization | Nucleus |
| 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, Immunohistology (formalin): 1-2 ug/mL for 30 minutes at RT, Western Blot 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user |
| Positive control | K-562 Cells, THP-1 cells, Lymph Node, Spleen, Hodgkin's Lymphoma, Colon Carcinoma. |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 °C. Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Regulatory status | For research use only (RUO) |
| Antibody/conjugate formulation | 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 |
| Validated in protein | Monospecific |
| Antibody research areas | Immunology, Transcription factors |
| Cell/tissue expression | B-cells |

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel |
|-------------------|--------------------|------------|------------|--------------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 |
| 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 |

Dye Features

[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

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 LI-COR Bioscience.