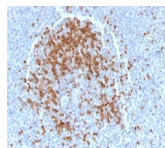


## TCL1 Monoclonal Mouse Antibody (TCL1/2078)



### Product Description

T-cell leukemia/Lymphoma Protein 1A (TCL-1A), also known as p14TCL1 is a product of the TCL1 gene that is involved in T-cell prolymphocytic leukemia (TPLL). T-PLL is a rare form of mature T-cell leukemia, which is consistently associated with chromosomal rearrangements characterized by the juxtaposition of the TCRA locus on chromosome 14q11 and the TCL1A gene on 14q32.13. TCL1 is overexpressed in Burkitt lymphoma, the majority of AIDS-related non-Hodgkin lymphoma-designated immunoblastic plasmacytoid lymphoma, lymphoblastic lymphoma, chronic lymphocytic leukemia, mantle cell lymphoma, follicular lymphoma, diffuse large B-cell lymphoma, and primary cutaneous B-cell lymphoma. 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 2078, Anti-TCL1 (TCL1/2078)**

### Product attributes

|                                       |                                                                                                                                                                                                                                                                                                                                                                                                                                                              |
|---------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Antibody number                       | #2078                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| Antibody reactivity (target)          | TCL1                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Antibody type                         | Primary                                                                                                                                                                                                                                                                                                                                                                                                                                                      |
| Host species                          | Mouse                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| Clonality                             | Monoclonal                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| Clone                                 | TCL1/2078                                                                                                                                                                                                                                                                                                                                                                                                                                                    |
| Isotype                               | IgG1, kappa                                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Molecular weight                      | 14 kDa                                                                                                                                                                                                                                                                                                                                                                                                                                                       |
| Synonyms                              | TCL1A; Lymphoma/leukemia, T-cell; Oncogene TCL1; P14 TCL1; T cell leukemia 1; T cell lymphoma 1; T cell lymphoma 1A; T-cell leukemia/lymphoma 1A; TCL1; TCL1A                                                                                                                                                                                                                                                                                                |
| Human gene symbol                     | TCL1A                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| Entrez gene ID                        | 10232                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| SwissProt                             | P56279                                                                                                                                                                                                                                                                                                                                                                                                                                                       |
| Unigene                               | 2484                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Immunogen                             | Recombinant human TCL1 protein fragment (around aa 2-109) (exact sequence is proprietary)                                                                                                                                                                                                                                                                                                                                                                    |
| Antibody target cellular localization | Nucleus & cytoplasm                                                                                                                                                                                                                                                                                                                                                                                                                                          |
| Verified antibody applications        | IHC (FFPE) (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. Immunohistochemistry (formalin): 1-2 ug/mL for 30 minutes at RT. 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                      | Ramos or Raji cells. Tonsil or T-cell Lymphoma.                                                                                                                                                                                                                                                                                                                                                                                                              |
| 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 array            | Monospecific                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Antibody research areas               | Cancer, Immunology                                                                                                                                                                                                                                                                                                                                                                                                                                           |

| 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>  |
| 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, LI-COR Bioscience.