

**PE/iFluor™ 594 Anti-human CD120a  
Antibody \*H398\***Catalog number: 112001Y0, 112001Y1, 112001Y2  
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

---

|                    |   |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving  |
| Concentration      | 0.1 mg/mL   |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

---

|                    |                                     |
|--------------------|-------------------------------------|
| Species Reactivity | Human                               |
| Class              | Primary                             |
| Clonality          | Monoclonal                          |
| Host               | Mouse                               |
| Isotype            | Mouse IgG2a                         |
| Immunogen          | CD120a (TNFR-I, TNFRSF1A, P60, P55) |
| Clone              | H398                                |
| Conjugate          | PE/iFluor™ 594                      |

**Biological Properties**

---

|             |   |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 594 under optimal conditions |
| Application | Flow Cytometry (FACS)   |

**Spectral Properties**

---

|                       |                |
|-----------------------|----------------|
| Conjugate             | PE/iFluor™ 594 |
| Excitation Wavelength | 566 nm         |
| Emission Wavelength   | 606 nm         |

**Applications**

---

H398 is an anti-human monoclonal antibody that is specific for the CD120a antigen. CD120a (alternatively called P55, TNFR-I, P60 or TNFRSF1A) is a 55 kD member of the TNFR superfamily that is expressed on the surface of cells such as B cells and T cells. CD120a is associated with a variety of biologically interesting macromolecules/ligands, for instance, TRAF, TRADD and Itα. CD120a is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD120a is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated

to PE/iFluor™ 594 (ex/em = 566/606 nm). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).