

**XFD680 Anti-human CD120a Antibody \*H398\***

Catalog Number: 11200180, 11200181

Unit Size: 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      | Lot specific (please consult certificate of analysis for given lot)         |
| 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          | AF680                               |

**Biological Properties**

---

|                       |                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
|-----------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Appearance            | Blue liquid                                                                                                                                                                                                                                                                                                                                                                                                                                     |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF680 under optimal conditions                                                                                                                                                                                                                                                                                                                                            |
| Application           | Flow Cytometry (FACS), Fluorescence Imaging                                                                                                                                                                                                                                                                                                                                                                                                     |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.<br><br><i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i> |

**Spectral Properties**

---

Conjugate                    AF680

Excitation Wavelength   681 nm

Emission Wavelength    704 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 XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 680 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).