

iFluor™ 660 Anti-human CD120a Antibody
H398Catalog number: 112000G0, 112000G1
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	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	iFluor™ 660

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 660
Excitation Wavelength	663 nm
Emission Wavelength	678 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 iFluor™ 660 (ex/em = 663/678 nm). It is compatible with the 640 nm laser and 670/30 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).