

**mFluor™ Green 620 Anti-human CD15
Antibody *AHC0040***Catalog number: 101510U0, 101510U1
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 | IgM |
| Immunogen | CD15 (3-FAL, X-hapten, Lewis X, SSEA-1) |
| Clone | AHC0040 |
| Conjugate | mFluor™ Green 620 |

Biological Properties

| | |
|-------------|--|
| Appearance | Purple liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------------|
| Conjugate | mFluor™ Green 620 |
| Excitation Wavelength | 525 nm |
| Emission Wavelength | 623 nm |

Applications

AHC0040 is an anti-human monoclonal antibody that targets the CD15 antigen. CD15 (sometimes called SSEA-1, fucosyltransferase or 3-FAL) is a glycoprotein that is found on the surface of cells such as granulocytes, stem cells and macrophages. CD15 is associated with a variety of biologically interesting macromolecules/ligands, for instance, selectins. CD15 is a fairly uncommon antibody target, with a little more than 5800

publications in the last decade. Even still, CD15 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of neuroscience cell markers, cell biology and immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 611/31 nm bandpass filter (for example, as in the Luminex Amnis CellStream).