

iFluor™ 532 Anti-human CD318 Antibody
CUB1Catalog number: 13180070, 13180071
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 IgG2b
Immunogen	CD318 (CDCP1, SIMA135)
Clone	CUB1
Conjugate	iFluor™ 532

Biological Properties

Appearance	Dark red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 532 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 532
Excitation Wavelength	537 nm
Emission Wavelength	560 nm

Applications

CUB1 is an anti-human monoclonal antibody that recognizes the CD318 antigen. CD318 (also known as CDCP1) is a 93 kD single-pass membrane protein that is found on the surface of cells like stem cells. CD318 is associated with a variety of biologically interesting macromolecules/ligands, in particular, N-Cadherin and syndecan-4. CD318 is a relatively rare antibody target, with fewer than 60 publications in the last decade. Even still,

CD318 is vital to neuroscience cell markers, neuroscience and stem cells research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 532 (ex/em = 537/560 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).