

Purified Mouse Anti-Cyanine5 Antibody
MonoclonalCatalog number: V102055
Unit size: 1.0 mg**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	, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Cyanine5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	Flow Cytometry (FACS)

Applications

(2,5-Dioxopyrrolidin-1-yl)
8-[(2Z)-3,3-dimethyl-2-[(2E,4E)-5-(1,3,3-trimethylindol-1-ium-2-yl)penta-2,4-dienylidene]indol-1-yl]octanoate;hydrochloride has shown evidence of use in a variety of applications, particularly in the field of biomedical research and pharmacological research.