

PE Mouse Anti-human DDIT4L Antibody
DDIT-03, monoclonalCatalog number: V103570
Unit size: 0.1 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	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	DDIT4L
Clone	DDIT-03
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

Applications

DNA damage-inducible transcript 4-like protein (alternatively called Protein regulated in development and DNA damage response 2, HIF-1 responsive protein RTP801-like or REDD-2) is a protein with a molecular weight of 22 kDa, found in the cytoplasm of cells. In humans, DNA damage-inducible transcript 4-like protein negatively regulates signal transduction.