

PE Mouse Anti-human FOLR2 Antibody
EM-35, monoclonal, Cross AdsorbedCatalog number: V103650
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	FOLR2
Clone	EM-35
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

Applications

Folate receptor β (alternatively called folate receptor, fetal/placental, placental folate-binding protein, FBP or folate receptor 2) is a protein with a molecular weight of 29 kDa, expressed in the anchored component of external side of plasma membrane, extracellular region and cell surface of cells. In Homo sapiens, folate receptor 2 has been associated with key functions such as folic acid receptor and signaling receptor activity. Folate receptor 2 aids in organismal processes, namely, folic acid transport, inflammatory response and folic acid metabolic process. Folate receptor 2 is thought to be essential to thymidylate synthase biosynthetic process.