

**Purified Mouse Anti-human LARGE1
Antibody *LARGE-02, monoclonal***Catalog number: V1031635
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	LARGE1
Clone	LARGE-02

Biological Properties

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

Applications

LARGE xylosyl- and glucuronyltransferase 1 is a 88 kDa transmembrane protein that can be expressed in the Golgi membrane, Golgi apparatus and integral component of Golgi membrane of cells. It is also known as glycosyltransferase-like protein and acetylglucosaminyltransferase-like 1A. In humans, glycosyltransferase-like protein binds to manganese ion. Sequencing of glycosyltransferase-like protein has demonstrated it contains 2 types of conserved structural units: cytoplasmic and lumenal domain. Glycosyltransferase-like protein has been closely linked to key functions such as acetylglucosaminyltransferase, xylosyltransferase and UDP-xylosyltransferase activity, and additionally, it takes part in organismal processes, for example, muscle cell cellular homeostasis, N-acetylglucosamine metabolic process and skeletal muscle tissue regeneration. Glycosyltransferase-like protein is clinically significant because abnormalities in its function have been closely linked to diseases like muscular dystrophy-dystroglycanopathy congenital with brain and eye anomalies type A6 (MDDGA6) and muscular dystrophy-dystroglycanopathy congenital with mental retardation type B6 (MDDGB6). Muscular dystrophy-dystroglycanopathy congenital with mental retardation B6, an autosomal recessive inheritance disorder characterized by lower limb hyperreflexia, babinski sign and elbow flexion contracture, has in specific been of interest to researchers.