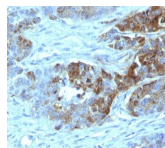


Blood Group H Type 2 Monoclonal Mouse Antibody (19-OLE)



Product Description

This antibody is applicable for staining ABO blood group O. The histo-blood group ABO involves three carbohydrate antigens: A, B, and H. This antibody recognizes the blood group H type 2 antigens, trisaccharide Fuc-a-1-2Gal1-4GlcNAc-a-1. A, B, and AB individuals express a glycosyltransferase activity that converts the H antigen to the A antigen (by addition of UDP-GalNAc) or to the B antigen (by addition of UDP-Gal), whereas O individuals lack such activity. It is expressed on endothelial cells, epithelial cells and granulocytes. Increased expression of this antigen has been observed on some tumor tissues such as gastric carcinomas, urothelial carcinomas, and colon carcinomas. This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Product attributes

Antibody number	#0947
Antibody reactivity (target)	Blood Group H Type-2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	19-OLE
Isotype	IgM, kappa
Molecular weight	Multiple
Synonyms	ABO Type O; Blood Group O; Blood Group H
Human gene symbol	ABO
Entrez gene ID	28
SwissProt	P16442
Unigene	654423
Immunogen	Mucinous colonic adenocarcinoma
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Immunohistology (formalin). Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes. Immunofluorescence 0.5-1 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	KG1 cells or human colorectal carcinoma.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Cancer, Hematology