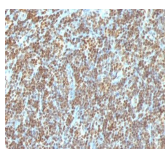


CD176 / T-F Ag Monoclonal Mouse Antibody (A68-B/A11)



Product Description

Recognizes a disaccharide epitope, Gal1-3GalNAc, of Thomsen-Friedenreich (TF) antigen. It is specific for both anomeric forms of the disaccharide (TF and TF, including related structures on the glycolipid) and shows no cross-reactivity with sialylated glycoprotein. The Thomsen-Friedenreich antigen acts as an oncofetal antigen, with low expression in normal adult tissues but increasing to fetal levels of expression in hyperplasia or malignancy. It is considered as a pan-carcinoma marker. This MAb is capable to agglutinate desialylated red blood cells. During metastasis, the ability of malignant cells to form multicellular aggregates via homotypic or heterotypic aggregation and their adhesion to the endothelium are critical. The tumor-associated carbohydrate Thomsen-Friedenreich antigen (Gal-GalNAc) is involved in tumor cell adhesion and tissue invasion. It also causes an immune response, and overexpression of the antigen causes cancer cells to be more sensitive to natural killer cell lysis. The Thomsen-Friedenreich antigen is suppressed in normal healthy cells and represents one of the few chemically well-defined antigens associated with tumor malignancy. The presence of the Thomsen-Friedenreich antigen on the surface of cancer cells may result from a divergence from the normal pathway for O-linked glycosylation in these cells, most likely caused by inappropriate localization of the enzymes involved in synthesis of the disaccharide.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#0937
Antibody reactivity (target)	CD176, T-F Ag
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	A68-B/A11
Isotype	IgM, kappa
Molecular weight of antigen	Multiple
Synonyms	T-F Antigen; TF Antigen; Asialoglycophorin; pan-carcinoma marker; CD176
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Neuraminidase-treated human red blood cells
Antibody target cellular localization	Plasma membrane
Species reactivity	Human, Mouse, Rat
Expected antibody applications	ELISA (published for clone), IHC (frozen) (published for clone)
Antibody application notes	Immunohistology (formalin), 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 tissues.
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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Tumor expression	Pan-carcinoma markers

This datasheet was generated on June 6, 2026 at 10:19:06 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd176-t-f-ag-thomsen-friedenreich-ag-a68-ba11/>