

FGF23 Monoclonal Mouse Antibody (FGF23/638)

Product Description

Fibroblast growth factor-1 (FGF-1, acidic FGF) and fibroblast growth factor-2 (FGF-2, basic FGF) are members of a family of growth factors that stimulate proliferation of cells of mesenchymal, epithelial and neuroectodermal origin. Additional members of the FGF family include the oncogenes FGF-3 (Int2) and FGF-4 (hst/Kaposi), FGF-5, FGF-6, FGF-7 (KGF), FGF-8 (AlGF), FGF-9 (GAF) and FGF-10 through FGF-23. Members of the FGF family share 30-55% amino acid sequence identity and similar gene structure. They are capable of transforming cultured cells when overexpressed by transfection. Cellular receptors for FGFs are members of a second multigene family, including four tyrosine kinases designated Flg (FGFR-1), Bek (FGFR-L), TKF and FGFR-3. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0638, Anti-FGF23 (FGF23/638)**

Product attributes

Antibody number	#0638
Antibody reactivity (target)	FGF23
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FGF23/638
Isotype	IgG1, kappa
Molecular weight	12-32 kDa
Synonyms	ADHR; FGF-23; FGFN; Fibroblast growth factor 23; HPDR2; HYPF; Phosphatonin; PHPTC; Tumor-derived hypophosphatemia-inducing factor
Human gene symbol	FGF23
Entrez gene ID	8074
SwissProt	Q9GZV9
Unigene	287370
Immunogen	Recombinant human FGF23 protein
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Human PBL cells or brain tumors
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. 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	Developmental biology, Signal transduction

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.

This datasheet was generated on May 4, 2024 at 04:46:59 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-fgf23-fgf23638/>