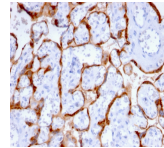


PAPP-A / Pappalysin-1 Monoclonal Mouse Antibody (PAPPA/2715)



Product Description

Pregnancy Associated Plasma Protein (PAPP-A) is found in maternal blood that increases as pregnancy progresses, although it is not specific to pregnancy. It is principally expressed in the syncytiotrophoblast of the placenta, which forms the main source of circulating maternal PAPP-A. It cleaves insulin-like growth factor binding proteins (IGFBPs), IGFBP-4 and IGFBP-5. IGFBP-4 cleavage is enhanced significantly in the presence of bound IGF, whereas IGFBP-5 cleavage is inhibited slightly by IGF presence. It is thought to be involved in local proliferative processes such as wound healing and bone remodeling. Low plasma level of this protein has been suggested as a biochemical marker for pregnancies with aneuploid fetuses. PAPPA has also been suggested as a potential biomarker of acute myocardial infarction and coronary artery disease (CAD).

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.

Product attributes

Antibody number	#2715
Antibody reactivity (target)	PAPP-A, Pappalysin-1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PAPPA/2715
Isotype	IgG2b, kappa
Molecular weight	181 kDa
Synonyms	ASBABP2; Aspecific BCL2 ARE binding protein 2; Differentially placenta 1 expressed protein; DIPLA1; IGF-dependent; IGFBP4ase; Insulin-like growth factor-dependent IGF-binding protein 4 protease (IGFBP-4 protease); PAPA; PAPP A; PAPPA1; Pappalysin-1; Pregnancy Associated Plasma Protein A (PAPP-A)
Human gene symbol	PAPPA
Entrez gene ID	5069
SwissProt	Q13219
Unigene	643599
Immunogen	Recombinant fragment (within aa 351-523) of human PAPP-A protein (exact sequence is proprietary)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Human
Positive control	Placenta
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
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
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

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, and of LI-COR Bioscience.