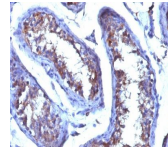


Progesterone Monoclonal Mouse Antibody (6-5E-3F)



Product Description

This MAb is specific for progesterone. It exhibits minimal cross reactivity with related compounds in ELISA. It reacts with Progesterone-11a-HMS-BSA: 100%; 5-beta-Pregnane-3,20-dione: 48%; 5-alpha-Pregnane-3,20-dione: 26.4%; 17-alpha-Hydroxyprogesterone: 2.5%; 20-alpha-Hydroxyprogesterone: 0.04%. Progesterone is a steroid hormone synthesized from the cholesterol derivative, pregnenolone, in the cortex of the adrenal gland. Progesterone is secreted by the corpus luteum and acts to prepare the endometrium for the implantation of a fertilized egg. During pregnancy, it is secreted by the placenta to prevent spontaneous abortion and to stimulate the development of mammary tissue to produce milk. Thus, progesterone plays a central role in the reproductive events associated with the establishment and maintenance of pregnancy. Luteinized theca cells of normal ovary secrete progesterone. The determination of progesterone concentrations in the body fluids is of great value for endocrinological investigations in women. This MAb may prove useful in identification of ovarian tumors.

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 0236, Anti-Progesterone (6-5E-3F)

Product attributes

Antibody number	#0236
Antibody reactivity (target)	Progesterone
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	6-5E-3F
Isotype	IgG1, kappa
Molecular weight	314.46Da
Synonyms	Pregn 4 ene 3 20 dione; Pregn 4 ene 3 20 dione
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Progesterone hemisuccinate conjugated to bovine serum albumin
Antibody target cellular localization	Cytoplasmic, Secreted (extracellular)
Species reactivity	All species
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	Ovary or 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
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	Endocrinology
Tumor expression	Ovarian cancer

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, or of LI-COR Bioscience.