

Progesterone Receptor antibody (mAb)

Catalog Nos: 61023, 61024

RRID: AB_2793476

Clone: 6A

Isotype: IgG2a

Application(s): ChIP, ICC, IF, IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 0.73 µg/µl

Molecular Weight: 100 kDa

Background: Progesterone Receptor (PR, NR3C3) is a nuclear receptor that binds the steroid hormone progesterone, translocates into the nucleus and activates transcription.

Immunogen: This Progesterone Receptor antibody was raised against a peptide within the C-terminus of human Progesterone receptor.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Validated Applications:

ChIP: 0.5 - 2 µg/ml dilution

Published Applications:

ICC/IF

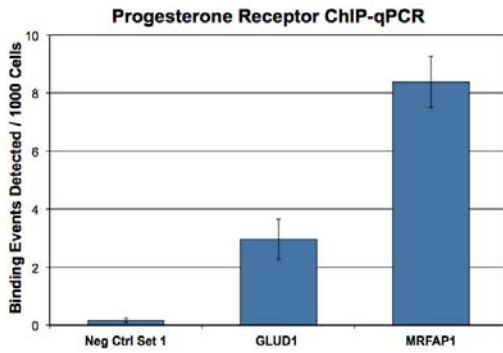
IHC

WB

See references for more information. Individual optimization may be required.

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Progesterone Receptor antibody (mAb) tested by ChIP.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 4.5 million treated MCF-7 breast cancer cells and 4 µl of Progesterone Receptor antibody. ChIP DNA was used in qPCR with the negative control primer pair or gene-specific primer pairs as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme, which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.