

# PRODUCT INFORMATION

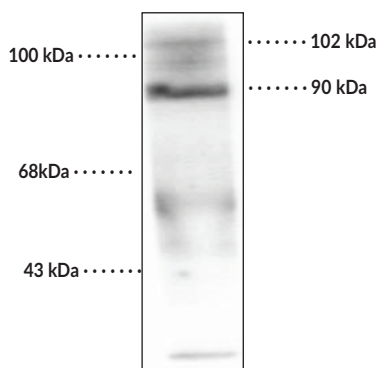
## Progesterone Receptor (Phospho-Ser<sup>190</sup>) Monoclonal Antibody (Clone 1154)

Item No. 10009762

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of protein G-affinity purified monoclonal antibody.
<b>Immunogen:</b>	Phosphopeptide from amino acid residues surrounding the phospho-Ser <sup>294</sup> of human progesterone receptor
<b>Molecular Weight:</b>	90/120 kDa
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
<b>Clone:</b>	1154
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG1
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:1,000 for IHC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



WB of T47D cell lysate prepared from cells that were incubated in the presence of the synthetic progestin agonist R5020 (500 nM) showing specific immunolabeling of the ~90 kDa PR-A isoform and the ~120 kDa PR-B isoform of the progesterone receptor phosphorylated at Ser<sup>190</sup>. The immunolabeling is specifically blocked by the phosphopeptide used as the antigen (not shown).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/30/2023

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

There is accumulating evidence to suggest that progesterone plays an essential role in the regulation of growth and differentiation of mammary glands and thus may play a key role in breast cancer.<sup>1</sup> The biological response to progesterone is mediated by two distinct forms of the human progesterone receptor (PR-A and PR-B forms). In most cell contexts, the B form functions as a transcriptional activator, whereas the A form functions as a transcriptional inhibitor of steroid hormones.<sup>2,3</sup> Recently it has been demonstrated that there is differential hormone dependent regulation of the phosphorylation of the A and B forms of the receptor.<sup>4</sup> Treatment of T47D breast cancer cells with progestin agonist increases the phosphorylation of Ser<sup>190</sup> and Ser<sup>294</sup> with different kinetics. These phosphorylation events may differentially affect the transcriptional activity of the receptor.

## References

---

1. Edwards, D.P. Regulation of signal transduction pathways by estrogen and progesterone. *Annu. Rev. Physiol.* **67**, 335-376 (2005).
2. Attia, G.R., Zeitoun, K., Edwards, D., *et al.* Progesterone receptor isoform A but not B is expressed in endometriosis. *J. Clin. Endocrinol. Metab.* **85(8)**, 2897-2902 (2000).
3. Lin, V.C.L., Woon, C.T., Aw, S.E., *et al.* Distinct molecular pathways mediate progesterone-induced growth inhibition and focal adhesion. *Endocrinology* **144(12)**, 5650-5657 (2003).
4. Clemm, D.L., Sherman, L., Boonyaratankornkit, V., *et al.* Differential hormone-dependent phosphorylation of progesterone receptor A and B forms revealed by a phosphoserine site-specific monoclonal antibody. *Mol. Endocrinol.* **14(1)**, 52-65 (2000).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM