

PRODUCT INFORMATION



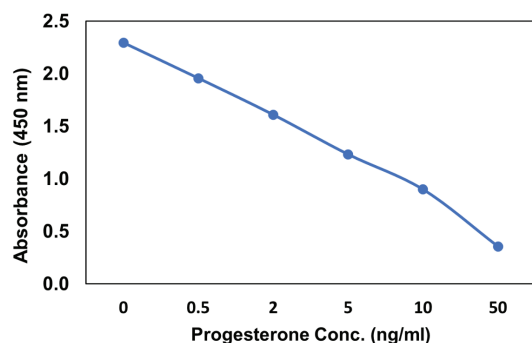
Progesterone Rabbit Monoclonal Antibody (Clone RM434)

Item No. 42801

Overview and Properties

Contents:	This vial contains 100 µg of protein-A affinity purified antibody.
Synonyms:	NSC 9704, NSC 64377
Immunogen:	Progesterone-KLH conjugate
Cross Reactivity:	(+) Progesterone
Species Reactivity:	(+) Species-independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM434
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA; the recommended starting dilution is 1-10 ng/well for ELISA. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



ELISA using Progesterone Rabbit Monoclonal Antibody (Clone RM434).

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, 01/14/2025

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

Progesterone is the biosynthetic precursor of all other steroid hormones.^{1,2} Progesterone is synthesized from cholesterol by the sequential action of desmolase in the mitochondria, which produces pregnenolone, followed by $\Delta^{4,5}$ -isomerase in the outer mitochondrial membrane and smooth endoplasmic reticulum of steroid-secreting cells.³ The main function of progesterone is to prepare the uterine lining for implantation of a fertilized ovum and to maintain pregnancy. Serum or plasma progesterone levels have been used as an index to monitor ovulation and investigate luteal function. Formulations containing progesterone have been used in support of early pregnancy in infertile women, prevention of endometrial hyperplasia, and the treatment of secondary amenorrhea and abnormal uterine bleeding. Cayman's Progesterone Rabbit Monoclonal Antibody (Clone RM434) can be used for ELISA.

References

1. Hormone action. *Biochemistry* 2nd ed., 1045-1085 (1988).
2. Disorders of steroidogenesis guide to steroid profiling and biochemical diagnosis. **1**, (2019).
3. Erickson, G.F. The ovary: Basic principles and concepts. A. Physiology. Endocrinology and *Metabolism* 973-1015 (1995).

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