

PRODUCT INFORMATION

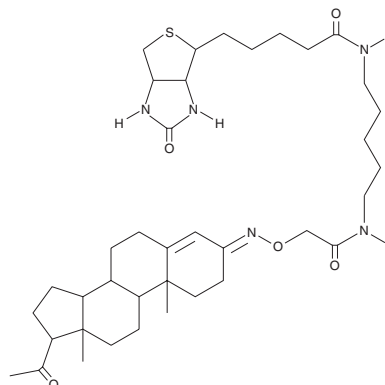


Progesterone 3-biotin

Item No. 9000644

Formal Name: N-(5-(2-(((E)-(17-acetyl-10,13-dimethyl-7,8,9,11,12,13,14,15,16,17-decahydro-1H-cyclopenta[a]phenanthren-3(2H,6H,10H)-ylidene)amino)oxy)acetamido)pentyl)-5-(2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamide

MF: C₃₈H₅₉N₅O₅S
FW: 698.0
Purity: ≥98%
Supplied as: A 5 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Progesterone, along with pregnenolone, is the biosynthetic precursor of all other steroid hormones. 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.¹ The main function of progesterone is to prepare the uterine lining for implantation of a fertilized ovum and to maintain pregnancy. Measurement of serum or plasma progesterone levels are used as an index to monitor ovulation and investigate luteal function. Progesterone 3-biotin is an affinity probe that allows progesterone to be detected or immobilized using the biotin ligand.

Reference

1. Erickson, G.F. The ovary: Basic principles and concepts. A. Physiology, Chapter 17, *in* Endocrinology and Metabolism. Felig, P., Baxter, J.D., and Frohman, L.A., editors, 3rd ed., McGraw-Hill, Inc. Health Professions Division, New York, 973-1015 (1995).

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, 11/24/2025

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