

PRODUCT INFORMATION



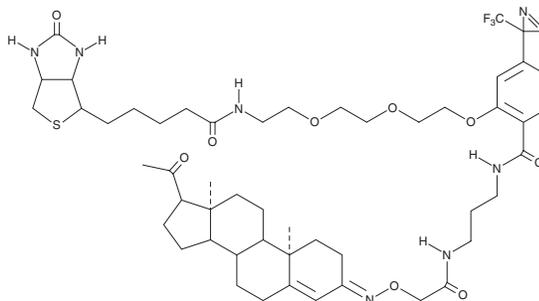
Progesterone-3-biotinyl Photoaffinity Probe

Item No. 9000891

Formal Name: N-(3-(2-(((10R,13S,E)-17-acetyl-10,13-dimethyl-1,2,6,7,8,9,10,11,12,13,14,15,16,17-tetradecahydro-3H-cyclopenta[a]phenanthren-3-ylidene)amino)oxyacetamido)propyl)-2-(2-(2-(5-(2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)ethoxy)ethoxy)-4-(3-(trifluoromethyl)-3H-diazirin-3-yl)benzamide

Synonyms:

MF: C₅₁H₇₁F₃N₈O₉S
FW: 1,029.2
Purity: ≥95% (mixture of isomers)
UV/Vis.: λ_{max}: 242 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Progesterone-3-biotinyl photoaffinity probe is supplied as a crystalline solid. A stock solution may be made by dissolving the progesterone-3-biotinyl photoaffinity probe in the solvent of choice. Progesterone-3-biotinyl photoaffinity probe is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of progesterone-3-biotinyl photoaffinity probe in these solvents is approximately 10, 1, and 15 mg/ml, respectively.

Progesterone-3-biotinyl photoaffinity probe is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Progesterone 3-biotinyl photoaffinity probe consists of active progesterone, a diazirine-UV (312 nm)-sensitive moiety, and a biotin tag. When activated by light, it binds covalently to its target and, thus, can be used to detect and observe the interaction of progesterone with target proteins. This probe has been used to help characterize the role of the human sperm progesterone receptor in male fertility.¹

Reference

1. Miller, M. R., Mannowetz, N., Lavarone, A. T., *et al.* Unconventional endocannabinoid signaling governs sperm activation *via* the sex hormone progesterone. *Science* **352(65285)**, 555-559 (2016).

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, 04/26/2024

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