

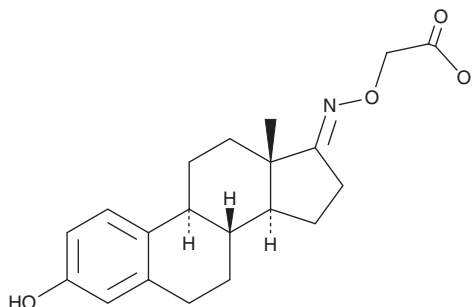
PRODUCT INFORMATION



Androgen Receptor Ligand 3

Item No. 45732

CAS Registry No.: 693227-00-8
Formal Name: (17E)-O-(carboxymethyl)oxime,
3-hydroxy-estra-1,3,5(10)-trien-
17-one
MF: C₂₀H₂₅NO₄
FW: 343.4
Purity: ≥98%
Supplied as: A 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

Androgen receptor ligand 3 is supplied as a solid. A stock solution may be made by dissolving the androgen receptor ligand 3 in the solvent of choice, which should be purged with an inert gas. Androgen receptor ligand 3 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Androgen receptor ligand 3 is an estrogen receptor (ER) ligand ($K_D = 0.39 \mu\text{M}$ in a radioligand binding assay) and an oxime derivative of estrone (Item No. 10006485).¹ Androgen receptor ligand 3 has been used in the synthesis of a specific and nongenetic IAP-dependent protein eraser (SNIPER) that decreases ER α levels in HT-1080 fibrosarcoma cells.² Biotin conjugates of androgen receptor ligand 3 have been used in yeast three-hybrid systems.³

References

1. Ratajczak, T., Sheppard, P.N., Capon, R.J., *et al.* The synthesis and study of some potential affinity labeling reagents for estrogen receptors. *Steroids* **38**(5), 537-555 (1981).
2. Itoh, Y., Kitaguchi, R., Ishikawa, M., *et al.* Design, synthesis and biological evaluation of nuclear receptor-degradation inducers. *Bioorg. Med. Chem.* **19**(22), 6768-6778 (2011).
3. Muddana, S.S. and Peterson, B.R. Facile synthesis of CIDS: biotinylated estrone oximes efficiently heterodimerize estrogen receptor and streptavidin proteins in yeast three hybrid systems. *Org. Lett.* **6**(9), 1409-1412 (2004).

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, 05/29/2026

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