

Estradiol (E2)-6-CMO-BSA Antigen

Estradiol (E2) linked to bovine serum albumin (BSA) via a 6-carboxymethyloxime (6-CMO) linker

This product combines estradiol with bovine serum albumin (BSA) via a stable 6-carboxymethyloxime (6-CMO) linker, enhancing immunogenicity and enabling reliable use in various assays.

Estradiol, a potent form of estrogen, plays a crucial role in regulating reproductive health, bone density, and neural function. It exerts its effects by binding to estrogen receptors, influencing gene expression and cellular responses across multiple organ systems. Abnormal estradiol levels are associated with various health issues, including osteoporosis, cardiovascular diseases, and hormone-dependent cancers.

Ordering Information

[Order Online »](#)

ENZ-PRT333-1000	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	3,17 β -Dihydroxy-1,3,5(10)-estratriene 6-(O-carboxymethyl)oxime-Bovine Serum Albumin
Appearance	White to faint yellow powder.
CAS	35048-47-6 (w/o BSA)
Formulation	Lyophilized.
Reconstitution	Reconstitute with PBS
Solubility	Soluble in PBS at 0.90 – 1.10 mg/ml.

Last modified: November 11, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)