

β-Estradiol

β-Estradiol is an estrogen and the primary female sex hormone and steroid, acting as a regulator of the estrous and menstrual female reproductive cycles. Directs the development of the female organs in embryogenesis and during puberty. It is an estrogen receptor agonist ($K_i=0.12$ and 0.13 nM for ER α and ER β , respectively).

Materials Provided

Catalog Number:	92-1128
Quantity Supplied:	1 vial(s), 100 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	272.38 g/mol
Source:	Synthetic
Purity:	≥ 99%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	23-25°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 mM
Reconstitution Volume:	3.671 mL
Reconstitution Solvent:	DMSO

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com