

## D(-)-Norgestrel

Norgestrel is a synthetic progesterone analog used as an oral contraceptive. It is a racemic mixture of levonorgestrel and dextronorgestrel. Levonorgestrel, the bioactive portion of Norgestrel, inhibits ovulation by causing cervical mucous to thicken. Levonorgestrel has 3 times higher affinity for the progesterone receptor than progesterone.

### Materials Provided

|                           |                           |
|---------------------------|---------------------------|
| <b>Catalog Number:</b>    | <b>92-1135</b>            |
| <b>Quantity Supplied:</b> | 1 vial(s), 10 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

|                          |   |
|--------------------------|---|
| <b>Molecular Weight:</b> | 312.45 g/mol                                    |
| <b>Source:</b>           | Synthetic                                       |
| <b>Purity:</b>           | ≥ 99%   |
| <b>Endotoxin Level:</b>  | N/A   |
| <b>Formulation:</b>      | N/A   |
| <b>Storage:</b>          | 2-8°C Please avoid multiple freeze/thaw cycles. |

### Reagent Preparation

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

|                                |        |
|--------------------------------|--------|
| <b>Stock Concentration:</b>    | 10 mM  |
| <b>Reconstitution Volume:</b>  | 3.2 mL |
| <b>Reconstitution Solvent:</b> | DMSO   |

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)