

9-cis-Retinoic acid

9-cis-Retinoic acid is a natural metabolite of vitamin A (retinol) and potent agonist of the retinoic acid and retinoid X receptors ($K_i=0.5-27\text{nM}$ and $3.8-12\text{nM}$, respectively). Mediates retinol functions for growth and development, and is active in embryonic development by intercellular signaling for the posterior portion of the embryo, and developmental patterning.

Materials Provided

Catalog Number:	92-1136
Quantity Supplied:	1 vial(s), 1 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:	300.44 g/mol
Source:	Synthetic
Purity:	≥ 98% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	1 mM
Reconstitution Volume:	3.328 mL
Reconstitution Solvent:	Ethanol

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