

Collagen II

Type II collagen makes up 50% of all protein in cartilage, particularly prevalent in articular and hyaline cartilage. Forms fibrils that allows cartilage to trap proteoglycan aggregate and provide tissue tensile strength.

Materials Provided

Catalog Number:	92-1090
Quantity Supplied:	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

Molecular Weight:	300 kg/mol
Source:	Bovine Articular Cartilage
Purity:	≥ 99% by SDS-PAGE, free of pepsin and proteoglycan
Endotoxin Level:	< 0.1 ng/μg (<1EU/μg).
Formulation:	N/A
Storage:	2-8°C in the dark Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	4 mg/mL
Reconstitution Volume:	2.5 mL
Reconstitution Solvent:	0.01M Acetic acid

Handling Instructions:

1. Upon arrival, the ligand should be stored at -20°C .
 2. Reconstitute 10 mg/vial of Collagen II with 2.5 mL of 0.01 M acetic acid to make a stock at 4 mg/mL. Mix gently to ensure Collagen II is in solution.
 3. This working stock solution can be stored at 4°C .
- Note: for optimal assay performance never freeze reconstituted ligand
4. The ligand is now ready for use.
 5. Serial dilution of the working stock should also be done in PBS+0.1% BSA.
 6. From the 4 mg/mL working stock, perform a 1:16 dilution in PBS + 0.1% BSA to obtain the highest dose reagent concentration (concentration of the ligand in the well) of 50 $\mu\text{g/mL}$ final (concentration of the dilution is at 5X = 250 $\mu\text{g/mL}$)
 7. Prepare eleven 2-fold serial dilutions of the 50 $\mu\text{g/mL}$ reagent in PBS + 0.1% BSA to obtain the lowest concentration of 24.4 ng/mL (final well concentration).
 8. Add a buffer control to complete the dose curve.
 9. Add 5 μL of each dilution into the assay plate.
 10. Aliquot unused ligand into vials and store at 4°C for 1 week or less. It is preferred to use a new aliquot of thawed ligand for best

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