

# CULTREX<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

---

## CellSpere<sup>™</sup>

**Catalog #:** 3450-048-05

**Size:** 15 ml

**Description:** CellSpere is a 36 kDa, bacillus-derived neutral metalloprotease that is recommended for recovering cells cultured on Cultrex<sup>®</sup> Basement Membrane Extracts. CellSpere will yield a single cell suspension far more gently and effectively than trypsin, collagenase, or other proteolytic enzymes, and it will not harm cells harvested for subcultivation or bioassays. In addition, CellSpere may be used for tissue dissociation. CellSpere cleaves fibronectin, collagen IV, and to a lesser extent Collagen I, but it does not cleave collagen V or Laminin.

**Source:** *Bacillus polymyxa*.

**Formulation:** 50 caseinolytic units/mL (5,000 caseinolytic units total) provided frozen in Phosphate Buffered Saline (PBS), pH 7.4. CellSpere is filtered through a 0.2 µm filter.

**Storage:** Store at -20°C in a manual defrost freezer in working aliquots (sterile plastic tubes); avoid freeze thaws.

### Quality Control:

- Solubilization of 1 mm thick layer of gelled Cultrex<sup>®</sup> Basement Membrane Extract (BME) at 37°C for two hours.
- No bacterial or fungal growth detected after incubation at 37°C for 14 days following USP XXIV Chapter 71 sterility test.
- No mycoplasma contamination detected by PCR.

**TREVIGEN<sup>®</sup>**

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**Instructions:**

1. Thaw CellSpers<sup>™</sup> in a 37°C waterbath and allow solution to equilibrate.
2. Using sterile technique, add two volumes of CellSpers<sup>™</sup> for every one volume of BME.
3. Incubate at 37°C for up to two hours for maximum dissolution.
4. Pipet mixture up and down to disperse cell suspension, and transfer suspension to a sterile centrifuge tube; record volume (one volume).
5. Add one volume of 10 mM EDTA to tube to stop CellSpers<sup>™</sup> reaction, and centrifuge at 250 g to pellet cells.
6. Wash cells with one volume of PBS, pH 7.4, and centrifuge at 250 g. Repeat this step twice.
7. Cells may be resuspended in fresh medium for continued culture or prepared for further analysis.

**Related Products:**

Catalog#	Description	Size
3455-024-K	Cultrex <sup>®</sup> 24 Well BME Cell Invasion Assay	24 inserts
3480-024-K	CultreCoat <sup>®</sup> 24 Well BME-Coated Cell Invasion Assay	24 inserts
3456-024-K	Cultrex <sup>®</sup> 24 Well Laminin I Cell Invasion Assay	24 inserts
3457-024-K	Cultrex <sup>®</sup> 24 Well Collagen I Cell Invasion Assay	24 inserts
3458-024-K	Cultrex <sup>®</sup> 24 Well Collagen IV Cell Invasion Assay	24 inserts
3455-096-K	Cultrex <sup>®</sup> 96 well BME Cell Invasion Assay	96 samples
3465-024-K	Cultrex <sup>®</sup> 24 well Migration Cell Assay	96 samples
3465-096-K	Cultrex <sup>®</sup> 96 well Migration Cell Assay	96 samples
3456-096-K	Cultrex <sup>®</sup> Laminin I Cell Invasion Assay	96 samples
3457-096-K	Cultrex <sup>®</sup> Collagen I Cell Invasion Assay	96 samples
3458-096-K	Cultrex <sup>®</sup> Collagen IV Cell Invasion Assay	96 samples
3490-096-K	CultreCoat <sup>®</sup> BME 96 Well Cell Adhesion Assay	96 samples
3496-096-K	CultreCoat <sup>®</sup> 96 Well Adhesion Protein Array	96 samples
3450-048-SK	Cultrex <sup>®</sup> Directed In Vivo Angiogenesis Assay (DIVAA <sup>™</sup> ) Starter	48 samples
3450-048-K	Cultrex <sup>®</sup> DIVAA <sup>™</sup> Kit	48 samples
3450-048-IK	Cultrex <sup>®</sup> DIVAA <sup>™</sup> Inhibition Kit	48 samples



**CellSpers<sup>™</sup>**  
 Cat#: 3450-048-05  
 Storage: ≤ -20 °C  
 (Manual Defrost)  
 1-800-873-8443