

CULTREX® Product Data

For Research Use Only. Not For Use in Diagnostic Procedures

Cultrex® R-Spondin1 (RSP01) Cells and Reagents

Catalog #: 3710-001-K

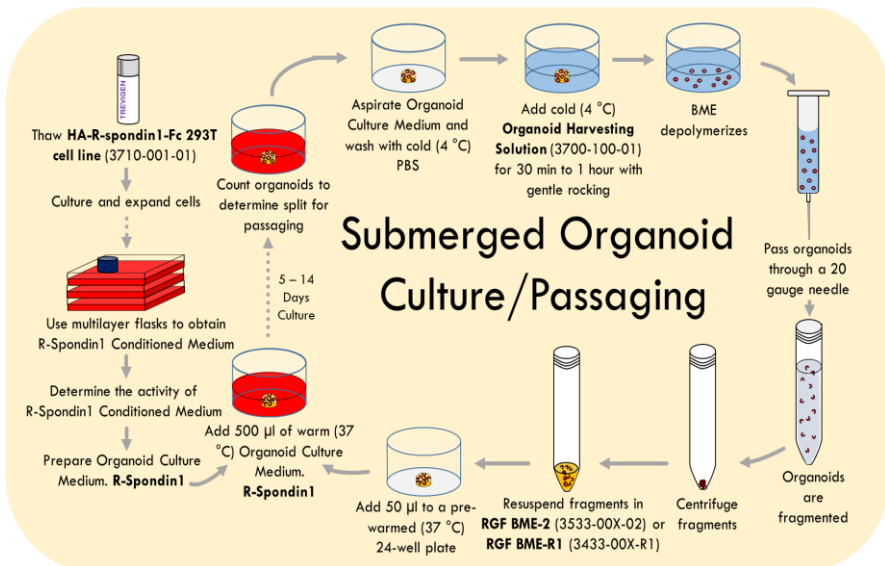
Contents:

<u>Cat#</u>	<u>Description</u>	<u>Size</u>	<u>Storage</u>
3710-001-01	HA-R-Spondin1-Fc 293T Cell Line	1 x 10 ⁶ cells	Liquid N ₂
3433-001-R1*	RGF BME-R1	1 ml	-80 °C
3533-001-02*	RGF BME-2	1 ml	-80 °C
3700-100-01*	Organoid Harvesting Solution	100 ml	4 °C

*Also included with Cultrex® R-spondin1 (RSP01) Cells at no additional charge

Description:

Organoid cultures represent the next generation of tissue culture models. These cultures are extracted directly from living tissues and, generally, they are never subjected to an artificial, tissue culture-treated plastic environment, providing more accurate disease modelling prospects. Our **Cultrex® R-spondin1 (RSP01) Cells** has been bundled in this presentation with 2 types of organoid culture qualified matrices, **RGF BME-2** and **RGF BME-R1**, and our **Organoid Harvesting Solution**, to provide your lab with essential tools to improve your organoid culture research. A summarized submerged organoid culture and passaging protocol is shown in the figure below. Visit our website for more information about each individual component.



© 2016 Trevigen, Inc. All Rights Reserved. Trevigen, Cultrex, and PathClear are registered trademarks, and 3-D Culture Matrix and DIVAA are trademarks of Trevigen, Inc.

TREVIGEN®

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 • www.trevigen.com

Related Products:

Catalog#	Description	Size
3432-005-01	Cultrex® Basement Membrane Extract, PathClear®	5 ml
3433-005-01	Cultrex® Reduced Growth Factor BME, PathClear®	5 ml
3533-005-02	Cultrex® Reduced Growth Factor BME, Type 2, PathClear®	5 ml
3632-005-02	Cultrex® Basement Membrane Extract, Type 3, PathClear®	5 ml
3445-005-01	Cultrex® 3-D Culture Matrix™ BME, PathClear®	5 ml
3446-005-01	Cultrex® 3-D Culture Matrix™ Laminin I	5 ml
3447-020-01	Cultrex® 3-D Culture Matrix™ Collagen I	100 mg
3434-005-02	Cultrex® Stem Cell Qualified RGF BME, PathClear®	5 ml
3415-001-03	Cultrex® Stem Cell Qualified Human BME, PathClear®	1 mg
3400-010-03	Cultrex® Stem Cell Qualified Laminin I, PathClear®	1 mg
3420-001-03	Cultrex® Stem Cell Qualified Human Fibronectin, PathClear®	1 mg
3420-001-03	Cultrex® Stem Cell Qualified Human Vitronectin, PathClear®	200 µg
3400-010-01	Cultrex® Mouse Laminin I	1 mg
3400-010-02	Cultrex® Mouse Laminin I, PathClear®	1 mg
3410-010-01	Cultrex® Mouse Collagen IV	1 mg
3440-100-01	Cultrex® Rat Collagen I	100 mg
3442-050-01	Cultrex® Bovine Collagen I	50 mg
3420-001-01	Cultrex® Human Fibronectin, PathClear®	1 mg
3416-001-01	Cultrex® Bovine Fibronectin, NZHD*	1 mg
3421-001-01	Cultrex® Human Vitronectin, PathClear®	50 µg
3417-001-01	Cultrex® Bovine Vitronectin, NZHD*	50 µg

Related Assays and Kits:

Catalog#	Description	Size
3500-096-K	Cultrex® 3D Spheroid Cell Invasion Assay	96 samples
3510-096-K	Cultrex® 3D Spheroid Fluorometric Proliferation/Viability Assay	96 samples
3511-096-K	Cultrex® 3D Spheroid Colorimetric Proliferation/Viability Assay	96 samples
3470-096-K	Cultrex® In Vitro Angiogenesis Assay, Tube Formation Kit	96 samples
3471-096-K	Cultrex® In Vitro Angiogenesis Assay, Endothelial Cell Invasion Kit	96 samples
3450-048-SK	Cultrex® Directed In Vivo Angiogenesis Assay (DIVAA™) Starter Kit	48 samples
3445-096-K	Cultrex® 3-D Culture BME Cell Proliferation Kit	96 samples
3448-020-K	Cultrex® 3-D Culture Cell Harvesting Kit	20 samples



**Cultrex® R-spondin1
(RSPO1) Cells**

Cat#: 3710-001-01
1-800-873-8443