

Maximize Your Pluripotential

With the TeSR™ Family of hPSC Culture Media

Maintenance of high quality pluripotent stem cells (PSCs) is critical to success in all applications of PSC research. The TeSR™ family of feeder-free maintenance media can help you minimize variation in your research. Each TeSR™ medium is based on published formulations¹-³ from the laboratory of James Thomson and offers unique features to fit your research needs.

Which TeSR[™] Feeder-Free Medium is Right for You?

mTeSR™1: Most Published



- >1100 peer-reviewed publications (www.stemcell.com/mTeSR1publications).
- Published protocols for a wide variety of applications including bioreactor expansion and single-cell cloning.
- Published protocols for lineage-specific differentiation of PSCs that have been maintained in mTeSR™1.



- Used with thousands of human ES and iPS cell lines. PSCs maintained for more than 7 years and in >50
- Contains pre-screened BSA to stabilize medium, aid in lipid/nutrient transport, and protect cultures from cellular toxins and stresses.1

TeSR™2: More Defined



- Modified formulation is similar to mTeSR™1, but with all xeno-free components, to produce a more defined medium.²
- Compatible with published mTeSR™1 protocols for a wide variety of applications.
- Contains recombinant human albumin to aid in lipid/nutrient transport and protect cultures from cellular toxins and stresses.

TeSR™-E8™: Simplified



- Contains only the 8 most essential components required for PSC maintenance. 3,4
- Cutting edge, xeno-free formulation.
- 433X less protein than mTeSR™1.



Scientists Helping Scientists™ | WWW.STEMCELL.COM

DOCUMENT #28221 VERSION 2.1.1 DEC 2016

Maximize Your Pluripotential

With the TeSR™ Family of hPSC Culture Media

The TeSR™ family of PSC media also includes optimized formulations for cryopreservation, reprogramming or use in differentiation assays. These products have the same base formulation, allowing researchers to establish a continuous TeSR™ media-based workflow for their PSC culture system.

Which TeSR™ Medium is Right for You?

TeSR™-E7™ & ReproTeSR™: Reprogramming



- Feeder-free, xeno-free media for human iPS cell induction.
- Generates iPS cells with high quality colony morphology for easy identification and rapid subcloning.
- TeSRTM-E7TM is based on the E8 formulation,³ with TGFβ removed for improved reprogramming efficiency
 of fibroblasts.
- ReproTeSR™ was optimized for reprogramming blood-derived cell types and seamlessly integrates with STEMCELL products for isolation and expansion of hematopoietic cells.
- ReproTeSR[™] can also be used to reprogram other somatic cell types, and can be paired with ReproRNA[™] for reprogramming fibroblasts.

TeSR™-E6 & **TeSR™-E5**: Screening and Differentiation

Neutral base medium

- Based on the E8 formulation,³ but do not contain bFGF or TGFβ.
- Lineage neutral formulation makes these media ideal for differentiation, screening assays, and other applications.
- With insulin removed in TeSR™-E5, this formulation faciliates differentiation to lineages in which insulin is a known inhibitor, such as cardiomyocyte.

mFreSR™ & FreSR™-S: Compatible Cryopreservation



- Defined, serum-free medium optimized for cryopreservation of hPSCs cultured in TeSR™ maintenance media.
- hPSCs cryopreserved in mFreSR[™] have thawing efficiencies higher than reported with conventional serum-containing media.
- FreSR™-S is animal component-free and optimized for cryopreservation of cells as single cells.

For a complete list of related products, including specialized cell culture and storage media, matrices, antibodies, cytokines and small molecules, visit www.stemcell.com/hPSCworkflow or contact us at techsupport@stemcell.com.

References:

- 1. Ludwig TE et al. (2006) Feeder-independent culture of human embryonic stem cells. Nat Methods 3(8): 637-646.
- 2. Ludwig TE et al. (2006) Derivation of human embryonic stem cells in defined conditions. Nat Biotechnol 24(2) 185-187.
- 3. Chen G et al. (2011) Chemically defined conditions for human iPS cell derivation and culture. Nat Methods 8(5) 424-429.
- 4. Beers J et al. (2012) Passaging and colony expansion of human pluripotent stem cells by enzyme-free dissociation in chemically defined culture conditions. Nat Protocols 7 2029-2040.

Copyright © 2016 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, Scientists Helping Scientists, mFreSR and ReproRNA are trademarks of STEMCELL Technologies Canada Inc. mTeSR, TeSR, E8, and E7 are trademarks of WARF. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

STEMCELL TECHNOLOGIES INC.'S QUALITY MANAGEMENT SYSTEM IS CERTIFIED TO ISO 13485. PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES UNLESS OTHERWISE STATED.