

## PRODUCT DESCRIPTION

AggreWell™800 plates (Catalog #27865/27965) have microwells 800 µm in diameter that can be used to aggregate human pluripotent stem cells (hPSCs) into embryoid bodies (EBs). EB formation is accomplished by adding a single cell suspension and then culturing for 24 hours. The resulting EBs are highly uniform in size. EBs generated using AggreWell™ plates are consistent in size and shape, and uniform within and between experiments.

## STABILITY AND STORAGE

This product has been aseptically manufactured using tightly controlled processes and sterilized by gamma irradiation. Please contact a STEMCELL Technologies representative if inner packaging is not intact.

Store at room temperature (15 - 25°C) away from direct light and heat. Product stable at 15 - 25°C for at least 5 years from date of receipt.

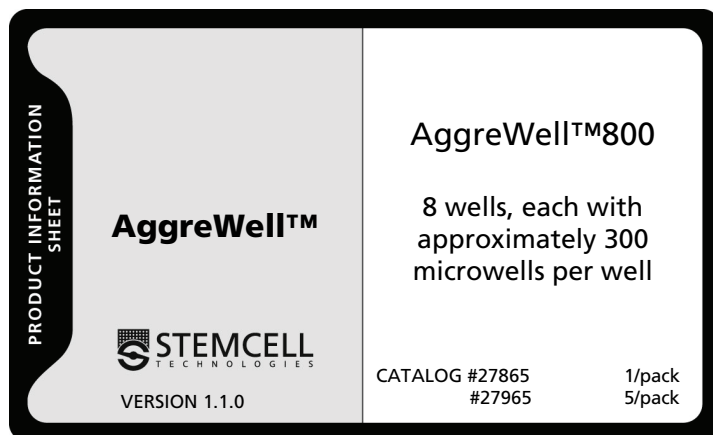
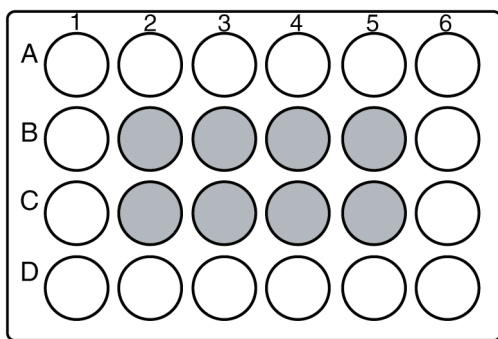
## DIRECTIONS FOR USE

For complete instructions on how to generate EBs from hPSCs using AggreWell™800 plates, see the Technical Manual "Reproducible and Uniform Embryoid Bodies Using AggreWell™ Plates" (Manual Catalog #29146) available on our website at [www.stemcell.com](http://www.stemcell.com).

AggreWell™800 wells are located in wells from B2 to C5 only, as shown below. Do not use empty wells for EB formation.

The non-reactive bonding agent may be visible underneath the microwell array, but it will not affect performance.

When generating EBs, it is essential to have a high quality starting population of undifferentiated hPSCs



## ADDITIONAL REQUIRED MATERIALS

PRODUCT	CATALOG #
PBS	37350
DMEM/F-12	36254
ACCUTASE™	07920
Y-27632 ROCK Inhibitor	07171 07172
Trypan Blue	07050
6-well ultra low adherence plate	27145
37 µm Reversible Strainer, small	27215
37 µm Reversible Strainer, large	27250
Conical tubes (50 mL)	e.g. BD Biosciences, Catalog #352070
Conical tubes (15 mL)	e.g. BD Biosciences, Catalog #352196
EB formation medium	Various <sup>a</sup>
EB suspension culture medium	Various <sup>b</sup>

ACCUTASE™ is a trademark of Innovative Cell Technologies, Inc. San Diego, CA

### Notes:

a, b: Please refer to the Technical Manual "Reproducible and Uniform Embryoid Bodies Using AggreWell™ Plates" (Manual Catalog #29146) available on our website at [www.stemcell.com](http://www.stemcell.com)



[WWW.STEMCELL.COM](http://WWW.STEMCELL.COM)

FOR RESEARCH USE ONLY. NOT FOR THERAPEUTIC OR DIAGNOSTIC USE.

MARCH 2011

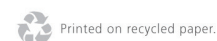
IN NORTH AMERICA TOLL-FREE T. 1 800 667 0322 TOLL-FREE F. 1 800 567 2899 T. 1 604 877 0713 F. 1 604 877 0704 E. [INFO@STEMCELL.COM](mailto:INFO@STEMCELL.COM)

IN EUROPE T. +33 (0)4 76 04 75 30 F. +33 (0)4 76 18 99 63 T. +33 (0) 4 76 04 75 30 F. +33 (0) 4 76 18 99 63 E. [INFO.EU@STEMCELL.COM](mailto:INFO.EU@STEMCELL.COM)

IN AUSTRALIA TOLL-FREE T. 1 800 060 350 F. +61 (03) 9338 4320 E. [INFO.AUS@STEMCELL.COM](mailto:INFO.AUS@STEMCELL.COM)

IN SINGAPORE T. (65) 6776 7754 F. (65) 6776 7114 E. [INFO.SG@STEMCELL.COM](mailto:INFO.SG@STEMCELL.COM)

#28508



Printed on recycled paper.

**RELATED PRODUCTS**

PRODUCT	CATALOG #
AggreWell™400	27845/27945
AggreWell™400Ex	27840/27940
AggreWell™ Medium	05893
mTeSR®1 Complete Kit for hES Maintenance Medium	05850/05870/05875/05857
TeSR™2 Complete Kit for hES Maintenance Medium	05860/05880
Y-27632 (Rock Inhibitor)	07171/07172
mFreSR™ Cryopreservation Medium	05854/05855
Collagenase IV (1 mg/mL)	07909
CryoStor™ CS10	07930
Dispase (1 mg/mL)	07923
Gelatin	07903
MEM Non-Essential Amino Acids, 10 mM (100X Concentrate)	07600
ACCUTASE™	07920
L-glutamine	07100
Trypsin-EDTA (0.05%)	07910
Trypsin-EDTA (0.25%)	07901
Anti-Oct 3/4 antibody	01550/01551
Anti-SSEA-1 antibody	01552
Anti-SSEA-3 antibody	01553
Anti-SSEA-4 antibody	01554
Anti-TRA-1-60 antibody	01555
Anti-TRA-1-81 antibody	01556

**REFERENCES**

1. Bauwens CL, Peerani R, Niebruegge S, Woodhouse KA, Kumacheva E, Husain M, Zandstra PW: Control of Human Embryonic Stem Cell Colony and Aggregate Size Heterogeneity Influences Differentiation Trajectories. *Stem Cells Express* doi: 10.1634/stemcells.2008-0183
2. Ungrin MD, Joshi C, Nica A, Bauwens C, Zandstra PW: Reproducible, Ultra-High-Throughput Formation of Multicellular Organization from Single Cell Suspension-Derived Human Embryonic Stem Cell Aggregates. *PLoS One* 3(2):e1565, 2008
3. Watanabe K, Ueno M, Kamiya D, Nishiyama A, Matsumura M, Wataya T, Takahashi JB, Nishikawa S, Nishikawa S, Muguruma K, Sasai Y: A ROCK inhibitor permits survival of dissociated human embryonic stem cells. *Nat Biotechnol.* 2007 Jun; 25(6):681-686

