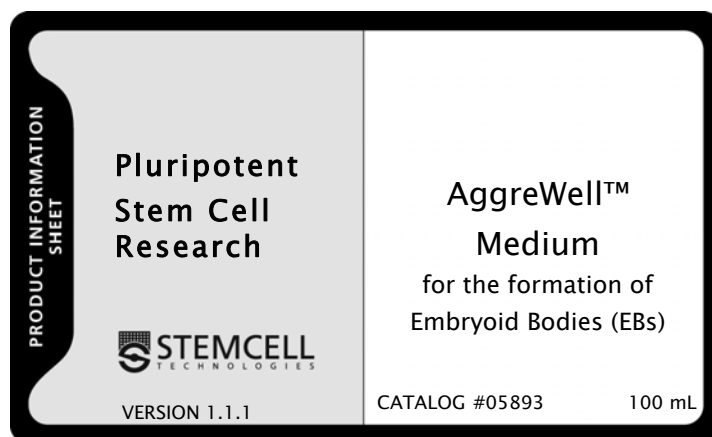


PRODUCT DESCRIPTION

AggreWell™ Medium (Catalog #05893) is designed for use with AggreWell™ plates for the generation and subsequent culture of embryoid bodies (EBs) formed from human pluripotent stem cells.

AggreWell™ Medium is suitable for the formation and culture of EBs generated from cells grown previously in mTeSR®1 or TeSR™2. AggreWell™ Medium does not contain pluripotency factors, but does contain basic fibroblast growth factor (bFGF, 40ng/ml).



STABILITY AND STORAGE

AggreWell™ Medium is stable at 2 - 8°C for up to 2 weeks.

AggreWell™ Medium can be stored at -20°C and is stable for up to 6 months.

DIRECTIONS FOR USE

1. Thaw AggreWell™ Medium (Catalog #05893) at room temperature (15 - 25 °C) or overnight at 2 - 8 °C.

If desired, AggreWell™ Medium can be aseptically dispensed into working aliquots and stored frozen at -20°C. Use aliquots within 3 months. Thawed aliquots should be used within 2 weeks. Do not refreeze aliquots after thawing.

2. For information on the formation of embryoid bodies using AggreWell™ plates, refer to the technical manual "Reproducible and Uniform Embryoid Bodies Using AggreWell™ Plates" (Manual Catalog #29146) available on our website at www.stemcell.com.
3. Depending on the differentiation protocol of choice, additional growth factors and/or serum may be added for subsequent directed differentiation.

RELATED PRODUCTS

PRODUCT	CATALOG #
AggreWell™ plates for the reproducible formation of uniformly sized embryoid bodies	27845/27945/27865/27965
Y-27632	07171/07172
mTeSR®1 Maintenance Medium for hESCs & hiPSCs	05850/05870
TeSR™2 Animal Protein-Free Maintenance Medium for hESCs & hiPSCs	05860/05880
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
FITC-conjugated goat anti-mouse IgG	10210
FITC-conjugated goat anti-mouse IgM	10211
APC-conjugated goat anti-rat IgM	10215
mFreSR® defined cryopreservation medium for hESCs and hiPSCs	05855/05854
ACCUTASE®	07920
Dispase (1 mg/mL)	07923



WWW.STEMCELL.COM

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

MAY 2010

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

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

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

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

#29939

