

StemBoost™ Reprogramming Cocktail Set II (1000X), Sterile-Filtered

CATALOG #: K870-1ML 1 ml
K870-5ML 5 ml
K870-set 5 x 1 ml

STORAGE: Store at -20°C. Protect from light.

DESCRIPTION: A convenient, sterile-filtered cocktail solution (1000X, in DMSO) containing four small molecules useful for enhancing the reprogramming efficiency of human somatic cells to induced pluripotent stem cells (iPSCs). The four products are shown in the following table:

Product	Cat. No.	Concentration in the vial (1000X)
CHIR99021	1677	3.0 mM
PD0325901	1643	0.5 mM
SB-431542	1674	2.0 mM
Thiazovivin	1681	0.5 mM

REFERENCES: Yu, J., *et al.* (2011). *PLoS ONE* **6**, e17557; Zhu, S., *et al.* (2010). *Cell Stem Cell* **7**, 651-655; Kawamata, M., and Ochiya, T. (2010). *Proc. Natl. Acad. Sci. USA* **107**, 14223-14228; Lin, T., *et al.* (2009). *Nature Methods* **6**, 805-808; Li, W., *et al.* (2009). *Stem Cells* **27**, 2992-3000.

HANDLING: Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

RELATED PRODUCTS:

- DiscoveryPak™ Stem Cell Fate Regulator Set I (Cat. No. K852-3)
- DiscoveryPak™ Stem Cell Fate Regulator Set II (Cat. No. K853-8)
- DiscoveryPak™ Stem Cell Fate Regulator Set III (Cat. No. K854-3)
- DiscoveryPak™ Stem Cell Fate Regulator Set IV (Cat. No. K855-3)
- DiscoveryPak™ Stem Cell Fate Regulator Set V (Cat. No. K865-5)
- DiscoveryPak™ Stem Cell Fate Regulator Set VI (Cat. No. K866-5)
- DiscoveryPak™ Stem Cell Fate Regulator Set VII (Cat. No. K866-10)
- (±) –Bay K 8644 (Cat. No. 1682-5)
- BIX01294 (Cat. No. 1678-5, 25)
- CHIR99021 (Cat. No. 1677-5, 25)
- EZSolution™ CHIR99021 (Cat. No. 1748-5)
- PD0325901 (Cat. No. 1643-2)
- PS-48 (Cat. No. 1869-5)
- Pyrintegrin (Cat. No. 1729-1,5)
- EZSolution™ Pyrintegrin (Cat. No. 1737-1)
- RepSox (Cat. No. 1894-5, 25)
- Reversine (Cat. No. 1851-1, 5)
- RG 108 (Cat. No. 1679-10,30)
- SB-431542 (Cat. No. 1674-1)
- EZSolution™ SB-431542 (Cat. No. 1872-1)
- Sodium Butyrate (Cat. No. 1609-100, 1000)
- StemBoost™ Reprogramming Cocktail Set I (1000X), Sterile-Filtered (Cat. No. K869-1ML, 5ML, 1EA)
- StemBoost™ YPAC Cocktail Set (1000X), Sterile-Filtered (Cat. No. K871-1ML, 5ML, 1set)
- Thiazovivin (Cat. No. 1681-1,5)
- EZSolution™ Thiazovivin (Cat. No. 1736-1)
- Tranylcypromine Hemisulfate (Parnate) (Cat. No. 1816-25,100)
- Valproic Acid, Sodium Salt (Cat. No. 1647-200)
- Y-27632,2 HCl (Cat. No. 1596-1, 5, 50)
- EZSolution™ Y-27632, HCl (Cat. No. 1784-5)

FOR RESEARCH USE ONLY! Not to be used on humans.