

**PRODUCT: DiscoveryPak™ Stem Cell Fate Regulator Set VIII (VC6T)**

**CATALOG #:** K887-4

**AMOUNT:** 4 modulators

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

**DESCRIPTION:** A convenient set of four small molecule modulators (V6CT) useful for generation of iPSC's (induced pluripotent stem cells) from mouse fibroblasts using a single gene, Oct4. The four products are shown in the following table:

Product	Cat. No.	Size
Valproic Acid, sodium salt (VPA)	1647-200	200 mg
CHIR99021	1677-5	5 mg
RepSox (E-616452)	1894-5	5 mg
Tranylcypromine Hemisulfate (Parnate)	1816-25	25 mg

**REFERENCE:** Li, Y., *et al.* (2011). *Cell Research* **21**, 196-204.

**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. K867-10)  
DiscoveryPak™ Stem Cell Fate Regulator Set IX (VC6TFZ) (Cat. No. K888-6)  
(±) –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)  
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)

**USAGE:** FOR RESEARCH USE ONLY! Not to be used in humans.