

DiscoveryPak™ Neuronal Cell Induction Small Molecules Set

CATALOG #:	K892-8
AMOUNT:	8 chemicals
FORM:	Solid
STORAGE:	-20 °C. Hygroscopic
DESCRIPTION:	A convenient set containing eight small molecules useful for induction of neuronal cells (hcnNs) from human fibroblasts. The eight products are shown in the following table:

Product	Cat. No.	Biological Activity	Size
Valproic Acid Sodium Salt (VPA)	1647	HDAC Inhibitor	200 mg
CHIR99021	1677	GSK-3 inhibitor	5 mg
RepSox	1894	TGFβ inhibitor	5 mg
Forskolin	1531	cAMP Activator	5 mg
SP600125	1669	JNK inhibitor	5 mg
Gö6983	9539	PKC inhibitor	1 mg
Y-27632 HCl	1596	ROCK inhibitor	5 mg
Dorsomorphin	1686	AMPK inhibitor	5 mg

USAGE: Dissolve all of the products except Valproic acid sodium salt (VPA) (Cat. No. 1647) in DMSO. VPA should be dissolved in D.I. water.

REFERENCES: Hu, W., *et al.* (2015). *Cell Stem Cell* 17, 204-212.

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 II (1000X), Sterile-Filtered (Cat. No. K870-1ML, 5ML, 1EA)
 StemBoost™ YPAC Cocktail Set (1000X), Sterile-Filtered (Cat. No. K871-1ML, 5ML, 1set)
 Thiazovivin (Cat. No. 1681-1,5)

FOR RESEARCH USE ONLY!



09/2015

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**)
Okadaic Acid, Sodium Salt (**Cat. No. 1764-025**)
Pepstatin A (**Cat. No. 1732-25, 100**)
Protease Inhibitor Cocktail (**Cat. No. K271-500**)
PMSF (**Cat. No. 1548-5**)

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